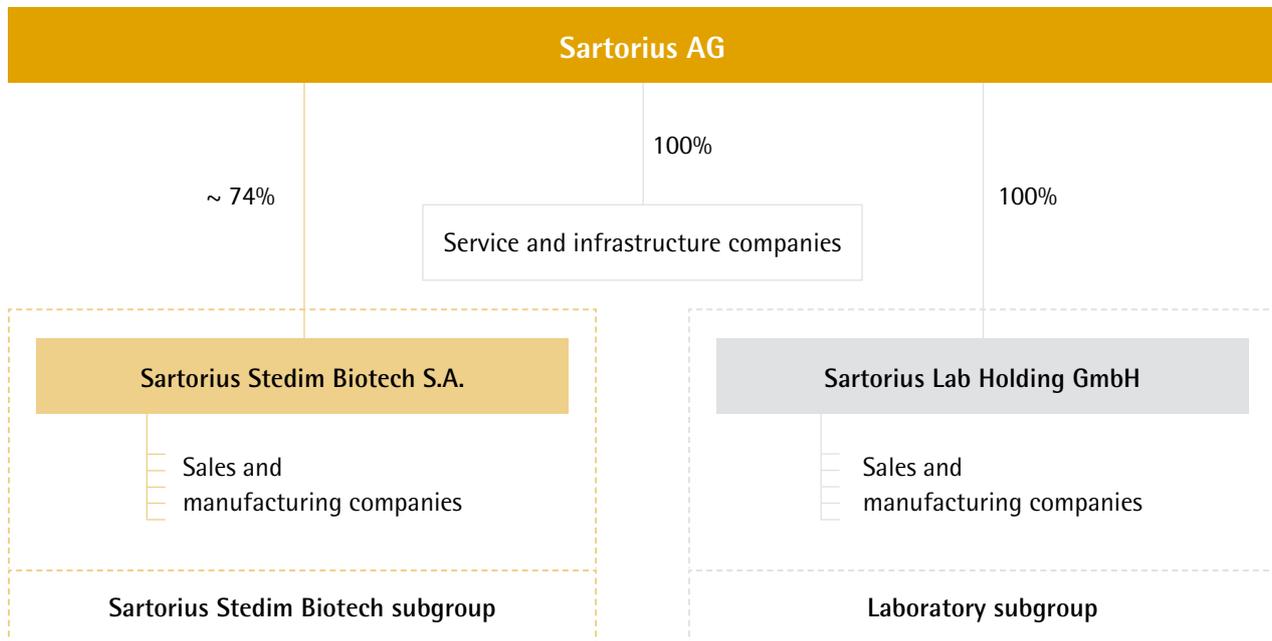


Combined Group Management Report

02

Structure and Management of the Group



Group Legal Structure

Sartorius is a globally operating company with subsidiaries in more than 30 countries. The holding company Sartorius AG is the parent corporation of the Sartorius Group. The corporation is headquartered in Göttingen, Germany, and is listed on the German Stock Exchange.

Sartorius manages its bioprocess business as a legally independent subgroup whose parent corporation is Sartorius Stedim Biotech S.A., which is listed on Euronext Paris. As of December 31, 2017, Sartorius AG held around 74% of the shares of Sartorius Stedim Biotech S.A. The Group's lab business is legally combined in a further subgroup whose parent company is Sartorius Lab Holding GmbH, in which Sartorius AG holds a 100% stake.

The consolidated financial statements include Sartorius AG and all major affiliates in which Sartorius AG has a controlling interest pursuant to IFRS 10.

Organization and Management of the Group

The Sartorius Group conducts its operating business in two divisions: Bioprocess Solutions and Lab Products & Services. The divisions each combine their respective businesses for the same fields of application and user groups, and share part of the infrastructure and central services.

The Group's central management entity is the Sartorius Group Executive Committee (GEC), which currently has six members. These are the Executive Board members of Sartorius AG and further senior managers with global responsibility.

The Sartorius Group is largely organized by function worldwide. Accordingly, the respective managerial responsibilities for the individual functions at the top management levels are performed across all sites and regions. Implementing the Group's various strategies and projects at the local level is the responsibility of the national affiliates. The management bodies of the local companies run their organizations in accordance with the applicable statutory provisions, articles of association and rules of procedure and in keeping with the principles of corporate governance that apply throughout the Sartorius Group worldwide.

Changes in the Group Portfolio

In the reporting year, Sartorius acquired three companies.

The portfolio of the Lab Products & Services Division was expanded in the area of bioanalytics in March 2017 by the acquisition of Essen BioScience Inc. Headquartered in Ann Arbor, Michigan, USA, this company develops and markets innovative real-time live-cell imaging and analysis systems for drug research.

Sartorius through its subgroup Sartorius Stedim Biotech additionally acquired MKS Instruments AB (Umetrics) based in Umeå, Sweden. This company, a leading specialist in data analytics software for modeling and optimizing biopharmaceutical development and manufacturing processes, had already been cooperating with Sartorius for around five years before the takeover. Umetrics was initially consolidated upon completion of the acquisition in April 2017.

Financial Controlling and Key Performance Indicators

The Sartorius Group is managed using a number of key performance indicators, which are also decisive for the determination of the variable remuneration component for the Executive Board and managers.

A key management parameter that Sartorius uses to measure the development of its size is currency-adjusted growth of sales revenue. The key profitability measure is EBITDA adjusted for extraordinary items, i.e. underlying EBITDA, and the corresponding margin.

Regarding the debt financing potential of the Sartorius Group, the key indicator is the ratio of net debt to underlying EBITDA for the last twelve months. Furthermore, the capex ratio, i.e., capital expenditures in proportion to sales revenue, represents a key control parameter.

In addition, the following financial and non-financial indicators are reported on a regular basis:

- Order intake
- Relevant net profit | Earnings per share
- Annual net profit | Earnings per share
- Equity ratio
- Net working capital
- Net cash flow from operating activities
- Number of employees

As a rule, the annual financial forecast that is published at the beginning of a fiscal year for the Group and the divisions refers to the development of sales revenue and of the underlying EBITDA margin. The expected capex ratio, as well as a directional forecast for the ratio of net debt to underlying EBITDA, is additionally indicated for the Group.

Strategy and Goals

The Sartorius Group is a leading international supplier of products and services covering a range from research and development to safe and efficient production of medications and vaccines using biotech methods.

Sartorius is increasingly focusing on the attractive biopharma market that is characterized by long-term and stable trends. For years, the global pharmaceutical market has been showing positive development primarily driven by the so-called pharmerging markets such as China, India and Russia in which access to medications was fueled by the expansion of national healthcare as well as higher private-sector spending. Additional significant trends are the demographic development in the Western industrialized countries and increasing life expectancy, which entail a greater need for medications.

The market for biotechnologically produced medications and vaccines has been expanding overproportionately for many years within the growing pharmaceutical market as a result of the introduction of new biopharmaceuticals and their further market penetration, among others. Biosimilars, or biological medications highly similar to already approved reference medicines, are considered factors for continued positive development in the future.

As part of its strategy as a total solutions provider, Sartorius has built up a broad and innovative portfolio of products and technologies across the entire value-added chain of its biopharmaceutical customers, both through the company's proprietary research and acquisitions. With this range, Sartorius aims to make complex, lengthy and expensive development of biopharmaceuticals and their production more efficient and less expensive for its customers. Moreover, Sartorius addresses the needs of laboratories of public research, defined as "academia."

The Sartorius Group conducts its operating business in two divisions: Bioprocess Solutions and Lab Products & Services. The divisions' market positioning and strategy are outlined as follows:

Bioprocess Solutions

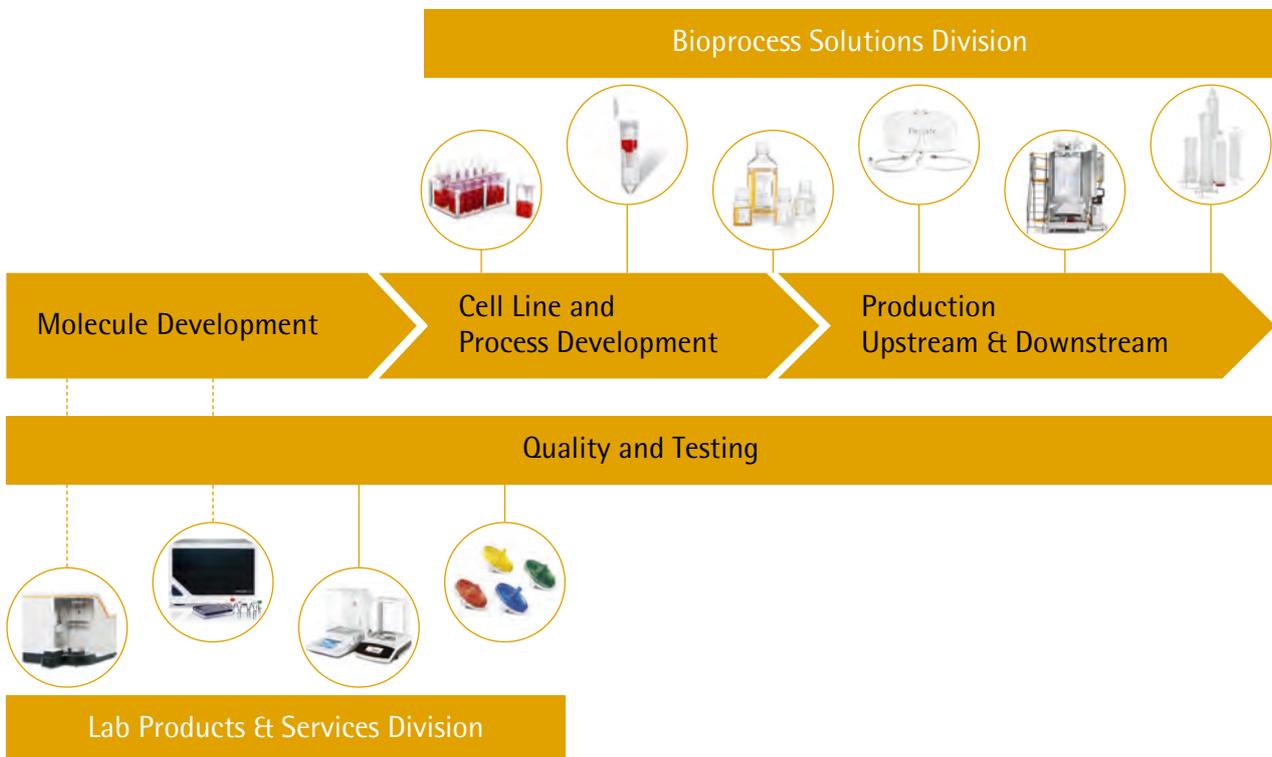
In the Bioprocess Solutions Division, Sartorius offers its customers innovative and efficient technologies and products extending from the development of cell lines and processes to manufacturing. These essentially cover cell culture media, cell line technologies, bioreactors, a wide range of products for separation, purification and concentration, as well as solutions for storage and transportation of intermediate and final biological products.

The division generates around three-quarters of its sales revenue with single-use products. Due to their cost advantages and higher flexibility compared with reusable technologies, such single-use products are increasingly being employed by the pharmaceutical industry. Through the acquisition of Umetrics, Sartorius is now a globally leading specialist in data analytics software for modeling and optimizing biopharmaceutical development and manufacturing processes.

With its global, specialized sales force, Bioprocess Solutions addresses an attractive market with above-average growth rates. As our customers' manufacturing processes are validated by the respective health authorities, product quality and assurance of supply are essential. We see the leading international market positions of this division as good stepping stones for sustained dynamic and profitable growth in the future. Beyond realizing our organic growth potential coupled with a continuous increase in profitability, we also aim to further expand the division's portfolio through complementary acquisitions and alliances.

Details on this division are provided in the chapter on Business Development of Bioprocess Solutions.

Strategic Focus on Biopharma Applications from Molecule Development to Production of Biopharmaceuticals



Lab Products & Services

In the laboratory business, Sartorius is positioned as a premium supplier of a wide array of instruments, consumables and services for sophisticated analyses, particularly in the laboratories of the pharmaceutical and. The division mainly addresses laboratories in the research and quality assurance sectors of the pharmaceutical and biopharmaceutical industries, as well as academic research institutes.

Primarily used in cell analysis and sample preparation, the product range of Lab Products & Services covers instruments such as laboratory balances, pipettes, bioreactors and lab water systems, as well as laboratory consumables, for example, filters and microbiological tests.

Based on its existing core expertise and considering its key customer requirements and the face-paced innovation in the industry, Sartorius decided in 2016 to expand its portfolio into the field of bioanalytics and to offer innovative technologies for the early phase of molecule development. More powerful procedures for analysis of cells are considered essential prerequisites for further medical progress, for example, in immunoncology and in antibody discovery and stem cell research. Apart from high growth in this segment, customers' digitalization efforts are expected to further propel progress in order to drive down R&D costs.

It was in this context that Sartorius acquired Essen BioScience in the reporting year. The innovative real-time live-cell imaging and data analysis systems designed by this U.S.-based company can contribute toward significantly accelerating discovery and development of new drugs as these automate steps, making high volumes of high-quality data available much faster for analysis. Essen BioScience offers high growth and margin potential.

The Lab Products & Services Division has a strong global market share in major product segments. Based on the company's history, market shares vary in part, depending on the region and product group. To realize its organic growth potential, the division is continuously expanding its direct sales organization.

With its comprehensive portfolio for sample preparation and innovative products for bioanalytics, the Lab Products & Services Division has a strong foundation for continuing strong organic growth. Based on the scalability of the division's business, this growth is projected to be accompanied by a steady increase in profitability. In addition, Sartorius plans to extend the lab division's portfolio by acquisitions.

Details on this division are provided in the chapter on Business Development of Lab Products & Services.



Sartorius 2020 Strategy

In 2011, Sartorius defined its strategy and long-term targets for 2020 to achieve profitable growth. At the beginning of 2016, close to the mid-point of this timeline, this 2020 plan was reviewed and updated.

The company's sales target of around €2 billion has been maintained despite the divestiture of the Industrial Technology Division in the interim. Growth is predominantly expected to be generated by Sartorius' existing portfolio, i.e., organically, and supplemented by acquisitions. In view of its profitability, Sartorius upgraded its margin target related to underlying EBITDA from around 23% to about 26% to 27% for 2020. This is assuming that the profitability of any future acquisitions would be at a level comparable to that of the existing business and that no significant changes in key exchange rates would occur.

Sartorius' 2020 targets are being implemented by various growth initiatives with the following areas of focus:

Regional Growth Initiatives

Regionally, North America and selected countries in Asia are at the focus of Sartorius' growth strategy.

North America is the world's largest market for the manufacture of biopharmaceuticals and laboratory products. Because North America is home to the main competitors for both company divisions, Sartorius has historically lower market shares in this region than in Europe and Asia. Accordingly, the company is striving to further gain market share, primarily by strengthening its sales and service capacities. Our second regional focus is on Asia, especially on China, South Korea and India. These markets have tremendous growth potential due to their expanding healthcare systems and increased spending by private households. Moreover, major production facilities for biosimilars are being set up in these countries, and contract manufacturers are strengthening their presence in this region. To participate in the momentum of the Asian markets in the best possible way, Sartorius already invested substantially in its sales infrastructure in this region.

Expansion of Product Portfolio

Regarding the further development of our portfolio, the 2020 strategy also provides for making acquisitions for both Group divisions. Such acquisitions will be primarily focused on adding complementary technologies and products that enhance the attractiveness of the company's portfolio even more from a customer perspective.

Infrastructure

The foundation for future profitable growth is constituted by efficient business processes, a powerful IT infrastructure and sufficient production capacities. Based on standardized business processes worldwide, Sartorius will be rolling out a new ERP system up to 2019 and is additionally investing continuously in the digitalization of its processes. Moreover, the company is considerably extending its production capacities at various locations, especially for filter and bag products. At the end of 2017, the two sites located in separate areas were combined for the most part into one location at the company's headquarters in Göttingen, Germany. This consolidation is expected to be completed by the end of 2018.

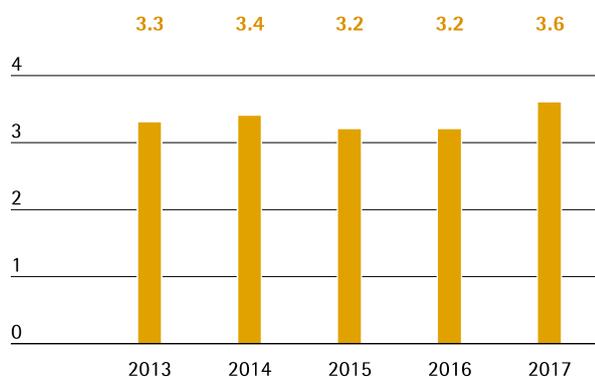
Macroeconomic Environment and Conditions in the Sectors

The sectors in which the Sartorius Group is active differ in their dependence on the economy. The Bio-process Solutions Division, for instance, operates in an environment that is largely independent of economic fluctuations. The Lab Products & Services Division, in contrast, is active in sectors whose development is more strongly affected by economic factors.

Macroeconomic Environment

According to the International Monetary Fund (IMF), growth of the global economy accelerated in 2017. Gross domestic product rose by 3.6% worldwide, an increase of approximately 0.4 percentage points over the prior year figure representing the most significant growth of the past five years. Whereas the industrialized countries reported a gain of 2.2%, growth in the emerging markets was at 4.6%.

Global Development GDP (2013 to 2017)
in %



Source: International Monetary Fund

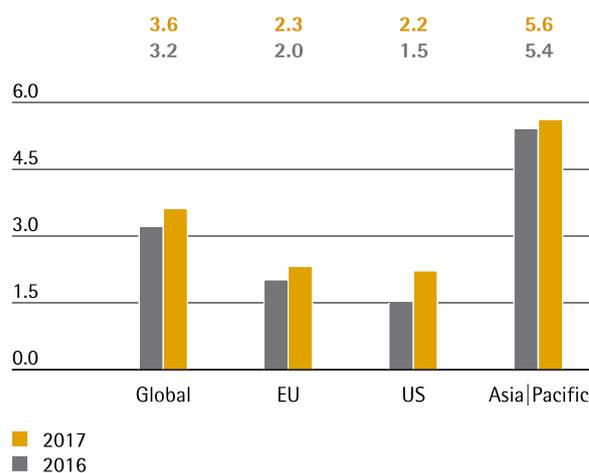
Nearly all regions participated in the global economic upturn. In the European Union, the increase amounted to 2.3%, a rise of around 0.3 percentage points on the previous year's figure. Growth in the reporting period was driven by the general revival in global trade and sustained strong domestic demand. France benefited in particular from the positive effects with growth of 1.6% in 2017 – significantly above the prior-year figure of 1.2%. In Germany, which is Sartorius' most important European market in terms of revenue, growth rose slightly to 2.0%. In the United Kingdom, consumer spending suffered from the depreciation of

the British pound. Thus, the increase of 1.7% fell just below the prior-year level of 1.8%.

The United States experienced an upturn in growth after the prior year's weak performance. Economic output rose approximately 2.2%, up from 1.5% in 2016. Growth was fueled by favorable financing conditions and rising business investments.

The Asia-Pacific economic area likewise registered an upward trend with an overall increase of 5.6% (previous year: 5.4%). Growth in the region received a boost from the revival in import demand in China. The world's biggest economy proved once again to be driving global growth with a rise of approximately 6.8% in total economic output. Sartorius also achieved significant revenue in South Korea, India and Japan. While economic output in South Korea grew by 3.0% (previous year: 2.8%), India posted a decline of 0.4 percentage points to 6.7%. In Japan, fiscal stimulus programs pushed up growth to 1.5% (previous year: 1.0%).

Gross Domestic Product by Region
in %



Source: International Monetary Fund

Exchange Rate Trends

In addition to the euro, the currencies relevant to the Sartorius Group include the U.S. dollar in particular as well as the South Korean won, the Chinese renminbi, the Swiss franc and the British pound.

The U.S. dollar lost considerable ground against the euro in the year under review. The exchange rate for the currency pair rose sharply in the second and third quarters of 2017 to reach a multi-year high of U.S. \$ 1.20 in September. The EUR/USD exchange rate fell back slightly after the U.S. Federal Reserve announced, as expected, plans to reduce its bond buying program, thus taking the first concrete steps towards initiating a reversal in monetary policy. Over the course of December the temporary drop was fully recovered. The exchange rate for the currency pair ended the year at U.S. \$ 1.20 as of December 31, 2017 compared with U.S. \$ 1.06 at the end of 2016.

The euro also made headway during the reporting year against the other currencies of relevance for Sartorius. For instance, it gained 1.2% on the South Korean won year-on-year. Compared with the Chinese renminbi and the Swiss franc, the euro rose 7.3% and 9.0%, respectively. The British pound fell 4.3% against the euro in 2017.

Interest Rate Trends

Interest rates remained at a very low level on average throughout the reporting year. The European Central Bank kept its key interest rate at 0.00%. The 3-month EURIBOR – i.e., the rate of interest on fixed-term deposits denominated in euros in interbank business – stood at -0.33% on December 31, 2017 (December 31, 2016: -0.32%).

Sources: International Monetary Fund: World Economic Outlook October 2017; Bloomberg.

Conditions in the Sectors

Sartorius' key customer groups include the biopharmaceutical and pharmaceutical industries as well as public research institutions. In addition, the company counts quality assurance laboratories in the chemicals and food industries among its customers. The progress of the Group's business accordingly depends on developments in these industries.

Continued Growth in the World's Pharmaceutical Markets

According to estimates from several market observers, the global pharmaceutical showed a positive development once again in 2017, with growth of approximately 4% to 5%. The market was again fueled by better access to health services in emerging markets and rising median ages in industrial countries, which led to an increasing demand for medications. By contrast, growth of the pharmaceutical market was dampened by government initiatives directed at reducing health spending as well as expiring patents for high-margin pharmaceutical products.

All the regions contributed to sector growth in the year under review. The United States continues to represent the biggest single market, followed by China, Japan and Germany. The strongest growth, though with a slower tempo, took place as in previous years in the "pharmerging markets", which include China, India, Brazil and Russia, among others. Performance of the European pharmaceutical market was impacted by increased efforts to control costs in the health sector. In the United States, expiring patents and tough competition among pharmaceutical manufacturers led to a moderate slowdown in market growth. Rising expenditures for newly launched medications had a positive impact.

Above-Average Growth in the Biopharmaceutical Market

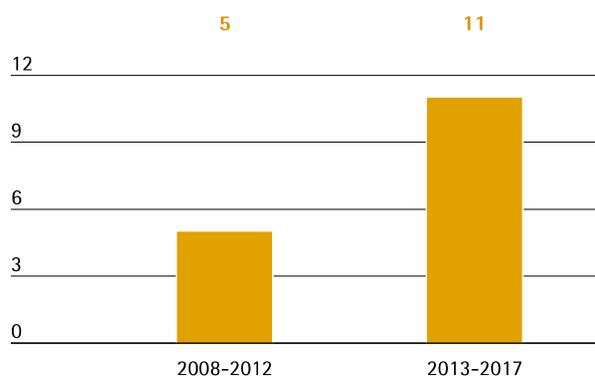
Given the company's highly specialized product portfolio, the trend in the biopharmaceutical market is especially relevant for Sartorius. The market for medications and vaccinations manufactured using biotech methods has grown faster than the rest of the pharmaceutical market for many years now. In 2017, the biopharmaceutical market was estimated by industry observers to have a volume of €214 billion, an increase of approximately 6% over the previous year. Overall, the proportion of sales revenue of the world's pharmaceutical market from medications manufactured using

biotech methods rose from around 25% to approximately 26%.

This growth is especially attributable to the launch of new biopharmaceutical drugs as well as additional market penetration of existing medicines, in part through expanded indications. A growing number of active substances manufactured using biotech production methods is being approved for the treatment of rare illnesses that have been untreatable thus far. Development activities in the pharmaceutical industry also reflect the steadily growing significance and rising acceptance of biopharmaceutical substances. For example, biopharmaceutical compounds account for more than 40% of the R&D pipeline.

In spite of market prospects that continue to be very positive, several large biopharmaceutical manufacturers reduced their inventories during the reporting year. The supply industry thus charted temporarily more subdued demand in North America and parts of Europe following several dynamic years.

Average Number of New Approvals of Biotech Medications in the USA per Year



Biosimilars, which are biological copycat medications, continued to play only a minor role in the growth of the biotechnology market in 2017. However, the market still advanced in the reporting year as new biosimilars were approved and launched on the market. At the time of this report, nine such drugs had already been approved in the United States and 36 in the European Union. China has currently approved 96 biosimilars and India 66. The biosimilars market is expected to see continual growth throughout the coming years due to expiration of the patents for a number of high-margin biopharmaceuticals. In addition, the regulatory, patent law-related and marketing hurdles that have hindered faster market penetration of biosimilars to date are likely to decrease step by step.

Single-Use Systems for Biopharmaceutical Production Continue to Gain Importance

Biotech production methods are much more complex and cost-intensive than traditional methods for producing medications. Consequently, manufacturers and suppliers are continuously looking to develop more efficient technologies. Single-use products play a decisive role in this effort: They require significantly less capital expenditure, reduce costs for cleaning and validation and minimize downtime. They also offer greater flexibility in production and help accelerate time to market.

Owing to these advantages, single-use technologies are already an established part of many process steps. Single-use systems are primarily employed in pre-commercial development activities and production phases and in small batch manufacturing. It can be expected that single-use technologies will become increasingly used for the production of high-volume commercial quantities. This is particularly relevant to the production of biotech substances whose clinical development takes place in single-use systems.

Moderate Growth in the Global Laboratory Market

Demand for laboratory instruments and supplies primarily stems from the biopharmaceutical and pharmaceutical industries as well as from public research initiatives. Laboratory products from Sartorius are also used by quality assurance laboratories in the chemicals and food sectors.

The global laboratory market grew by approximately 2.8% in the reporting year according to Frost & Sullivan. Europe reported higher growth of 1.9% thanks to the increased rate of economic growth. The United States, the largest market for laboratory products, registered growth of 3.2%. On the demand side, uncertainty regarding both planned legislation and the trend in government spending for academic and public research activities led to weak demand in the first quarter of 2017 in particular. These factors became increasingly less significant as the year progressed, however, resulting in a positive overall trend for the entire year – especially due to strong demand from the biopharmaceutical industry.

Significant growth was once again reported in Asian countries such as China and India, in which the laboratory market enjoyed an above-average expansion of 7.6% (China) and 8.6% (India).

Competitive Position

The competitive environment in each of our two divisions is as follows:

The Bioprocess Solutions Division operates as a total solution provider, covering the core process steps in biopharmaceutical production and upstream process development. It holds a leading global position in key technological fields, and offers the sector's most extensive portfolio of single-use technologies. The Bioprocess Solutions Division's principal competitors include Merck KGaA, Danaher, General Electric and Thermo Fisher.

The Lab Products & Services Division is positioned as a premium provider of laboratory instruments along with excellent services, and serves both R&D laboratories and quality assurance laboratories, with a focus on the biopharmaceutical industry. The product range includes laboratory balances, pipettes, and instruments for bioanalytics as well as a wide range of consumables. The division ranks among the leading providers worldwide in most of these areas. Principal competitors include Mettler Toledo Intl. Inc., Thermo Fisher Inc., Merck KGaA, Danaher Corp. and Eppendorf AG.

Sources: QuintilesIMS Institute: Outlook for Global Medicines through 2021, December 2016; Evaluate Pharma: World Preview 2017, Outlook to 2022, June 2017; BioPlan: 14th Annual Report and Survey of Biopharmaceutical Manufacturing Capacity and Production, April 2017; Deloitte: 2017 Global Life Science Outlook, October 2016; Frost & Sullivan: 2017 Mid-year Report: Forecast and Analysis of the Global Market for Laboratory Products, October 2017; www.fda.gov; FDA-Approves-Fifth-Biosimilar-in-US-First-for-Amgens-Blockbuster-Enbrel, www.raps.org

Group Business Development

Sales Revenue and Order Intake

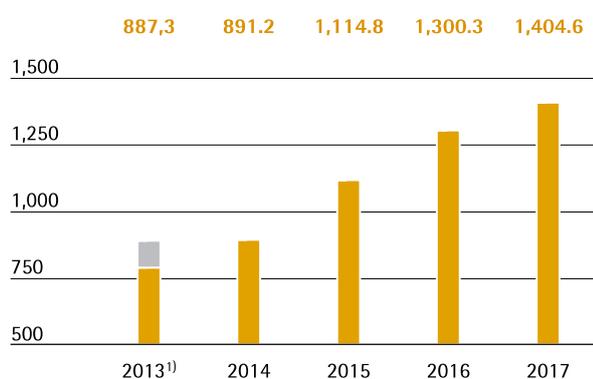
In the reporting year, the Sartorius Group continued to show positive development. Sales revenue thus rose 9.3% in constant currencies to €1,404.6 million, with the amounts contributed by organic growth and by acquisitions being approximately equal. On a reported basis, i.e., including currency effects, sales grew by 8.0%. Sartorius therefore reached its adjusted forecast for the third quarter of 2017.

Order intake also showed positive development in the reporting period, considerably exceeding sales. In 2017, orders increased 13.7% in constant currencies to €1,501.4 million. On a reported basis, this gain was 12.5%.

For a full comparison of the Group's business development with its forecast, see page 50.

Sales Revenue 2013 – 2017

€ in millions



¹⁾ Including sales revenue of discontinued operations

Sales Revenue and Order Intake

€ in millions	2017	2016	in % reported	in % wb ¹⁾
Sales revenue	1,404.6	1,300.3	8.0	9.3
Order intake	1,501.4	1,334.7	12.5	13.7

¹⁾ In constant currencies

Sales Driven by Both Divisions

After two years of exceptionally high growth in 2015 and 2016, sales revenue for the Bioprocess Solutions Division increased in the reporting year at a slightly more moderate rate of 4.9% to €1,010.3 million. The division's business performance was impacted by several simultaneous, temporary effects, primarily in the Americas region. Therefore, transient bottlenecks in the business for cell culture media, inventory destocking by a few relatively large customers and interruption of production at the plant in Puerto Rico in the wake of Hurricane Maria all had a dampening effect. As a result of these impacts, the division's growth was lower than originally expected. Acquisitions of the two companies kSep and Umetrics contributed around one percentage point to the division's growth.

The Lab Products & Services Division also saw exceptionally dynamic growth due to strong development of organic sales and the expansion of its portfolio into bioanalytics. Sales revenue rose year over year by 22.0% to €394.2 million, with around 14 percentage points attributable to the consolidation of IntelliCyt and ViroCyt acquired in 2016, as well as of Essen BioScience taken over in 2017. Growth was driven by all regions and all product areas.

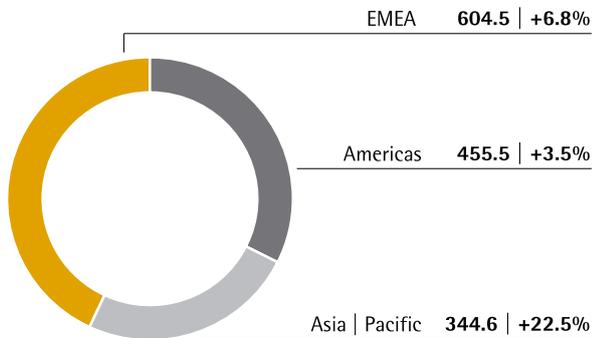
Sales by Division

€ in millions	2017	2016	in % reported	in % cc
Bioprocess Solutions	1,010.3	975.0	3.6	4.9
Lab Products & Services	394.2	325.3	21.2	22.0

Further information on the business development of the Group divisions is given on pages 42 et seq. for the Bioprocess Solutions Division and on pages 46 et seq. for the Lab Products & Services Division.

Significant Gains in All Regions

Sales Revenue and Growth¹⁾ by Region²⁾
 € in millions unless otherwise specified



¹⁾ In constant currencies

²⁾ Acc. to customers' location

In the reporting year, Sartorius recorded gains in all regions. In EMEA, the region contributing the highest share of revenue of around 43%, sales rose 6.8% to €604.5 million. While the Bioprocess Solutions Division achieved a gain of 2.5%, sales revenue for the Lab Products & Services Division was up 17.7% year over year.

The Americas region, in which sales grew at an especially strong rate over the past years, accounted for around 32% of consolidated revenue in 2017. Sales were up 3.5% to €455.5 million. Development of the Bioprocess Solutions Division was especially impacted by the temporary effects mentioned above; as a result, the division reported that sales decreased by 3.3% to €357.1 million. By contrast, the Lab Products & Services Division showed exceptionally dynamic development. Fueled by strong organic growth and by the acquisition of Essen BioScience, the division's sales were up 39.5%.

Business development for the Asia|Pacific region, which accounted for around 25% of Group revenue, showed the highest dynamics in the reporting year, with sales growing by 22.5% to €344.6 million. Sales revenue for the Bioprocess Solutions Division rose sharply by 25.3%, partly due to equipment shipments to fill relatively large orders. The Lab Products & Services Division also reported a double-digit gain of 16.6%. All growth rates are in constant currencies unless otherwise stated.

Sales by Region

€ in millions	2017	2016	in % reported	in % cc
EMEA	604.5	569.1	6.2	6.8
Americas	455.5	445.4	2.3	3.5
Asia Pacific	344.6	285.8	20.6	22.5

Development of Costs and Earnings

In the reporting year, the cost of sales stood at €697.7 million. In comparison with sales revenue growth of 8.0%, the cost of sales increased underproportionately by 4.4%, which was due to product mix effects and economies of scale. The cost of sales ratio was 49.7% relative to 51.4% a year ago.

Selling and distribution costs rose overproportionately relative to sales revenue, by 15.7% to €296.8 million. Accordingly, the ratio of selling and distribution costs to sales revenue increased from 19.7% in the previous year to 21.1%.

Expenses for research and development rose year over year by 15.8% to €68.8 million. The ratio of R&D expenses to sales revenue was 4.9%, slightly above the prior-year level of 4.6%.

Concerning general administrative expenses, Sartorius reported an 11.5% increase to €83.3 million, which can be primarily attributed to the expansion of the IT functional area. In relation to sales revenue, general administrative expenses rose from 5.7% in the previous year to 5.9% in the reporting year.

In fiscal 2017, the balance of other operating income and expenses was -€38.6 million relative to -€20.6 million a year earlier. This year-on-year change was driven, inter alia, by extraordinary items that totaled -€35.5 million in the reporting year (2016: -€30.7 million). These items were essentially incurred due to various cross-divisional projects, expenses related to the most recent acquisitions and to the consequences of Hurricane Maria.

In the year under review, the Group's EBIT edged down slightly by 0.5% to €219.4 million, especially due to higher PPA amortization in connection with the acquisition of Essen BioScience. The Group's EBIT margin was 15.6% because of higher depreciation as well as extraordinary items (2016: 17.0%).

The financial result was –€20.8 million in 2017 relative to –€16.4 million in 2016. This was essentially attributable to increased interest expenses.

In the reporting year, tax expenses of €39.3 million were substantially lower than the prior-year total of €59.1 million. This can essentially be attributed to the U.S. tax reform recently approved, which in the Sartorius Group resulted in positive extraordinary items due to reclassification of deferred tax liabilities of around €16 million in the USA. Because of this effect, the company's tax rate was just below 19.8%, down from 29.0% in the year before.

Net profit for the period rose 9.9% to €159.3 million (2016: €145.0 million).

Net profit attributable to shareholders of Sartorius AG amounted to €114.7 million in the reporting year relative to €102.9 million in 2016. Non-controlling interest stood at €44.6 million (2016: €42.1 million), which essentially reflected shares in Sartorius Stedim Biotech S.A. not held by the Sartorius Group.

Statement of Profit or Loss

€ in millions	2017	2016	in %
Sales revenue	1,404.6	1,300.3	8.0
Cost of sales	-697.7	-668.5	-4.4
Gross profit on sales	706.8	631.8	11.9
Selling and distribution costs	-296.8	-256.6	-15.7
Research and development costs	-68.8	-59.4	-15.8
General administrative expenses	-83.3	-74.7	-11.5
Other operating income and expenses	-38.6	-20.6	-87.3
Earnings before interest and taxes (EBIT)	219.4	220.5	-0.5
Financial income	11.0	5.7	93.5
Financial expenses	-31.7	-22.0	-44.1
Financial result	-20.8	-16.4	-27.0
Profit before tax	198.6	204.1	-2.7
Income taxes	-39.3	-59.1	33.6
Net profit for the period	159.3	145.0	9.9
Attributable to:			
Equity holders of Sartorius AG	114.7	102.9	11.5
Non-controlling interest	44.6	42.1	6.0

Underlying EBITDA

The Sartorius Group uses underlying EBITDA – earnings before interest, taxes, depreciation and amortization and adjusted for extraordinary items – as its key profitability indicator to provide a better picture of its operating development, also in an international comparison. More information on extraordinary items is provided on page 138.

Reconciliation from EBIT to Underlying EBITDA

€ in millions	2017	2016
EBIT	219.4	220.5
Extraordinary items	35.5	30.7
Amortization depreciation	98.3	74.2
Underlying EBITDA	353.2	325.4

In fiscal 2017, the Sartorius Group increased its earnings yet again. Its underlying EBITDA thus rose by 8.5% to €353.2 million. The Group's respective underlying EBITDA margin was 25.1% (2016: 25.0%), thus reaching our forecast adjusted for the third quarter.

	Underlying EBITDA € in millions	Underlying EBITDA margin in %
Group	353.2	25.1
Bioprocess Solutions	282.4	28.0
Lab Products & Services	70.8	18.0

Underlying EBITDA for the Bioprocess Solutions Division totaled €282.4 million, up 3.3% from €273.5 million a year ago. These earnings would have been even slightly higher without unfavorable currency effects. Moreover, the product mix effects in the reporting year also dampened underlying earnings. The division's margin reached the previous year's level of 28.0%. The Lab Products & Services Division substantially increased its earnings: its underlying EBITDA rose 36.3% from €51.9 million a year earlier to €70.8 million. The division's earnings margin of 18.0% was higher than the prior-year figure of 16.0% due to the positive effects resulting from the acquisition of Essen BioScience.

Relevant Net Profit

The relevant net profit attributable to the shareholders of Sartorius AG rose 8.6% from €132.6 million in 2016 to €144.0 million in 2017. This figure is the basis for determining the profit to be appropriated and is calculated by adjusting for extraordinary items and eliminating non-cash amortization, and is based on the normalized financial result as well as the corresponding tax effects for each of these items. The underlying earnings per ordinary share climbed by 8.6% to €2.10, up from €1.93 a year earlier, and by 8.6% per preference share to €2.11 euros, up from €1.94 euros a year ago.

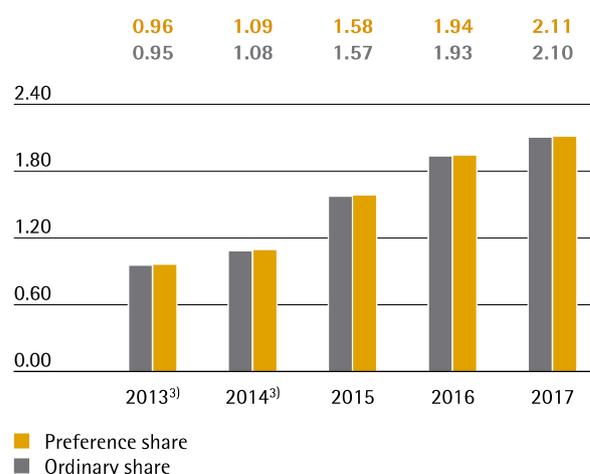
€ in millions	2017	2016
EBIT	219.4	220.5
Extraordinary items	35.5	30.7
Amortization	34.8	19.9
Normalized financial result ¹⁾	-17.7	-13.6
Normalized income tax (2016: 30%, 2015: 30%) ²⁾	-78.9	-77.3
Underlying earnings	193.1	180.3
Non-controlling interest	-49.1	-47.7
Underlying earnings after taxes and non-controlling interest	144.0	132.6
Underlying earnings per share		
per ordinary share (in €)	2.10	1.93
per preference share (in €)	2.11	1.94

¹⁾ Financial result excluding fair value adjustments of hedging instruments and non-periodic expenses and income

²⁾ Underlying income tax, based on the underlying profit before taxes and non-cash amortization

Underlying Earnings per Share¹⁾²⁾

in €



¹⁾ Adjusted for extraordinary items and non-cash amortization, based on the normalized financial result, including the corresponding tax effects

²⁾ 2013-15 adjusted for stock split; rounded values

³⁾ Including discontinued operation

Further information on earnings development and extraordinary items for the Group divisions is given on pages 42 et seq. and 46 et seq.

Research and Development

Sartorius continually invests in both new and further development of its products as well as the integration of new products through alliances. In 2017, the Sartorius Group spent €68.8 million on R&D, corresponding to an increase of 15.8% over the previous year's investment of €59.4 million. Owing to significantly higher sales, the ratio of R&D costs to sales revenue stood at 4.9%, slightly above last year's 4.6%.

The IFRS require that certain development costs be capitalized on the balance sheet and then amortized over subsequent years. During the reporting year, these development investments amounted to €27.4 million compared with €18.3 million the year before. This equates to a share of 28.5% (previous year: 23.5%) of the Group's total R&D expenses. Amortization related to capitalized development costs totaled €8.4 million during the reporting period (previous year: €9.4 million). These expenses are disclosed in the cost of sales.

We pursue a strategic intellectual and industrial property rights policy across our divisions to protect our expertise. Sartorius systematically monitors compliance with these rights on a cost|benefit basis to determine which specific individual rights are to be maintained.

In 2017, we filed a total of 161 applications for intellectual and industrial property rights (2016: 158). As a result of these applications, including those of prior years, we were issued 263 patents and trademarks during the reporting year (2016: 257). As of the balance sheet date, we had a total of 3,699 patents and trademarks in our portfolio (2016: 3,416).

Further information is provided in the sections covering the individual divisions on pages 42 et seq. and 46 et seq.

Capital Expenditures

Against a backdrop of strong organic growth in recent years, Sartorius is currently making above-average investments in the development of new capacity. In the reporting year, these development investments amounted to €209.4 million compared with €152.1 million the year before. At 14.9% (previous year: 11.7%), the ratio of investment to sales revenue was at the upper end of the forecast.

Our largest investment project is the expansion of our Group headquarters in Göttingen. As part of the consolidation of two formerly separate sites, our capacities for the production of filters and laboratory instruments will be extended there. New administration buildings were also opened during the reporting year.

And at our location in Yauco, Sartorius is investing in the expansion of its production capacities for filters and single-use bags. (Further information is provided in the sections covering the individual divisions on page 45.)

To extend the capacities in cell line development that we gained when Cellca was acquired in 2015, we began construction during the reporting year on a new building in Ulm.

Sartorius also made progress on the implementation of a new ERP system: in 2017 it was introduced at further sites in German, Switzerland and India and implementation is scheduled to be completed in 2019.

Employees

The numbers of employees reported include all staff members of the Sartorius Group, except for vocational trainees, interns, employees on extended leaves of absence and those participating in an early retirement plan. Numbers are reported as head counts.

On December 31, 2017, the Sartorius Group employed 7,501 people in 29 countries – 590 or 8.5% more than in the previous year. These figures include 214 employees who joined the Group in the reporting year through the acquisitions of the two companies Essen BioScience and Umetrics.

Employees

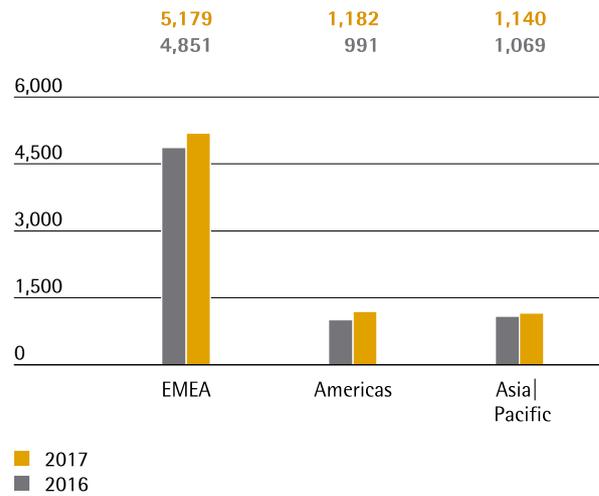
	2017	2016	Growth in %
Group	7,501	6,911	8.5
Bioprocess Solutions	4,980	4,584	8.6
Lab Products & Services	2,521	2,327	8.3

In the Bioprocess Solutions Division, 4,980 people were employed as of the end of the reporting year, 386 or 8.6% more than at year-end 2016. The acquisition of Umetrics resulted in a gain of 36 new staff.

The Lab Products & Services Division employed 2,521 people as of the year ended December 31, 2017; this was an increase of 194 or 8.3% compared with the year-earlier period. The acquisition of Essen BioScience contributed significantly to this growth of the workforce, with 178 new people joining Group staff.

Employees in central administrative functions were allocated to the divisions in proportion to cost of the services they performed during the year.

Employees by Region

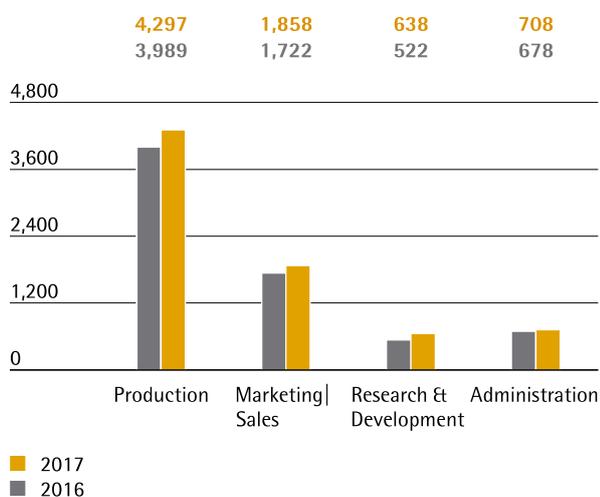


Regionally, EMEA with the highest number of employees added 6.8% or 328 new people. As of the end of the reporting year, Sartorius employed 2,807 people in Germany, or 37.4% of the total Group workforce, of which most staff members work at the Göttingen-based headquarters.

The Americas region charted an increase of 19.3% or 191 new people, again reporting the highest percentage of staff hired in the year under review.

The number of employees in the Asia|Pacific region grew by 6.6% or 71 people.

Employees by Function



Most employees at Sartorius work in production; at the end of 2017, they numbered 4,297, thus 7.7% more than at the end of 2016. In our production head count, we also include service staff and employees who perform services such as testing as part of our customers' core processes.

In marketing and sales, 1,858 people were employed at year-end, close to 8% more than a year ago.

In R&D, head count was 638 for the year ended December 31, 2017, thus around 22% higher than twelve months earlier. In this case, the company's acquisitions considerably increased the workforce, with 39 employees from Essen BioScience and 17 staff members from Umetrics joining the Sartorius Group.

As of the reporting date, 708 employees worked in administration, 4.4% more than in the prior year.

Expressed in percentages, 57.3% of Sartorius employees worked in production, close to 25% in marketing and sales, 8.5% in research and development and slightly more than 9% in administration (nearly 10% in 2016) at the end of 2017.

For more information on employees, see the Non-financial Group Declaration starting on page 86.

Net Worth and Financial Position

Cash Flow

In fiscal 2017, Sartorius reported a net cash flow from operating activities of €206.5 million relative to €170.4 million a year ago. This equates to growth of €21.2% and essentially reflects the increase in underlying EBITDA.

Cash outflows from investing activities were €199.1 million, 33.8% higher than a year ago. The investments in the reporting period were related, inter alia, to the expansion of production capacities at our site in Yauco, Puerto Rico, and to consolidation and extension of Group headquarters in Göttingen, Germany.

Cash outflows of €356.0 million were related to acquisitions of the companies Essen BioScience and Umetrics in the fiscal year ended December 31, 2017 (December 31, 2016: €119.5 million). In the prior year, cash outflows were due to the acquisitions of the three companies IntelliCyt, ViroCyt and kSep.

On the whole, net cash flow from investing activities and acquisitions|divestitures was -€555.1 million relative to -€268.2 million in the comparable period.

Net cash flow from financing activities, which amounted to €347.9 million (2016: €104.7 million), is predominantly attributable to the financing of the acquisitions mentioned above. Furthermore, this figure includes payment of dividends for fiscal 2016.

Cash Flow Statement Summary

€ in millions	2017	2016
Net cash flow from operating activities	206.5	170.4
Net cash flow from investing activities and acquisitions	-555.1	-268.2
Net cash flow from financing activities	347.9	104.7
Cash and cash equivalents	59.4	62.0
Gross debt	955.0	547.9
Net debt	895.5	485.9

Consolidated Statement of Financial Position

The balance sheet total of the Sartorius Group was €2,297.7 million as of the end of fiscal 2017 or €544.7 million above the previous year's figure of €1,753.0 million. This increase is predominantly attributable to the acquisitions of Essen BioScience and Umetrics.

Non-current assets rose €460.3 million to €1,625.7 million due to investments and acquisitions made in the reporting year.

Current assets at €672.0 million were €84.4 million higher than in the previous year, particularly because of the buildup in working capital.

Key Figures for Working Capital

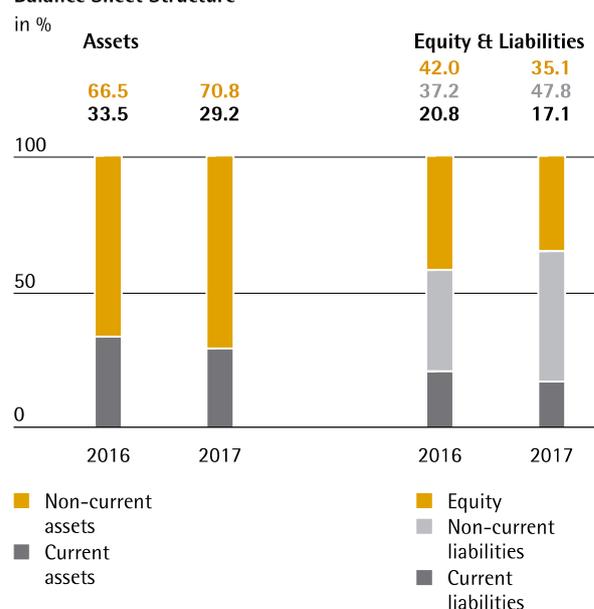
in days		2017	2016
Days inventory outstanding			
Inventories Sales revenue	x 360	63	61
Days sales outstanding			
Trade receivables Sales revenue	x 360	72	66
Days payables outstanding			
Trade payables Sales revenue	x 360	35	33
Net working capital days			
Net working capital ¹⁾ Sales revenue	x 360	99	94

¹⁾ Sum of inventories and trade receivables less the trade payables

Equity increased, driven by earnings, from €736.8 million to €806.6 million. The equity ratio for the Sartorius Group of 35.1% (2016: 42.0%) continued to remain at a comfortable level.

Current and non-current liabilities for the Sartorius Group rose from €1,016.2 million a year ago to €1,491.2 million in the reporting year. Above all, the acquisitions previously mentioned, as well as the mainly growth-induced increase in working capital, contributed to this figure.

Balance Sheet Structure

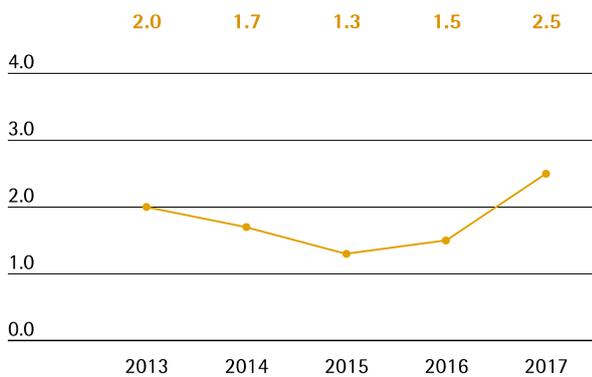


At the end of fiscal 2017, gross debt stood at €955.0 million relative to €547.9 million in fiscal 2016 and is comprised of liabilities to banks, including note loans ("Schuldscheindarlehen"), as well as liabilities from finance leases. The year-over-year increase is essentially attributable to the acquisitions of Umetrics and Essen BioScience, which were refinanced by placement of a note loan. Net debt, defined as gross debt less cash and cash equivalents, rose from €485.9 million a year ago to €895.5 million.

Ratio of Net Debt to Underlying EBITDA

Regarding the Sartorius Group's debt financing capacity, the ratio of net debt to underlying EBITDA is a key financial ratio that is calculated as the quotient of net debt and underlying EBITDA. As of December 31, 2018, it was 2.5 relative to 1.5 a year earlier and thus within the range of our forecast adjusted to allow for acquisitions.

Ratio of Net Debt to Underlying EBITDA^{1,2)}



¹⁾ Underlying

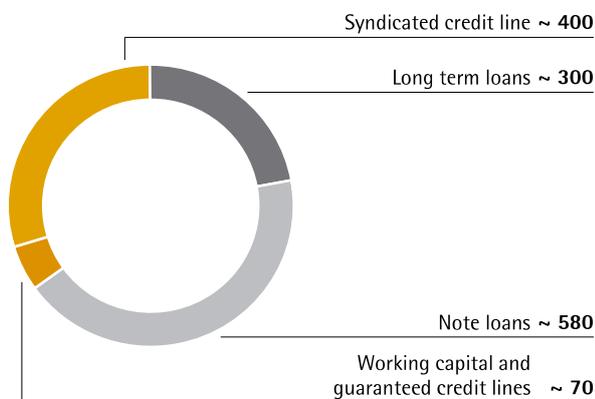
²⁾ Including discontinued operation 2013 and 2014

Financing | Treasury

The Sartorius Group is financed on a well-diversified basis, which covers both its short-term cash requirements and its long-term strategy.

Main Financing Instruments

€ in millions



A major pillar of this financing is the syndicated credit line of €400 million, whose maturity date will run until December 2021.

In addition to note loans ("Schuldscheindarlehen") that were placed in 2012 and 2016 and had a 2017 year-end volume of around €230 million, the company extended its financing in 2017 by a further €350 million note loan with maturities of 5, 7 and 10 years. In addition, several long-term loans totaling around €300 million are available to expand the company's production capacities, among other projects. Beyond these loans, Sartorius has diverse working capital and guaranteed credit lines totaling around €70 million.

The financing instruments mentioned above for the Sartorius Group comprise those with both fixed and variable interest rates. We pursue the strategy of using floating rate bank liabilities to hedge in part against an increase in the general interest rate level.

As a consequence of its global business activities, the Sartorius Group is exposed to fluctuations in foreign exchange rates. The Group's main currencies include the U.S. dollar as well as several others, such as the South Korean won, the Chinese renminbi, the Swiss franc and the British pound. Using its global manufacturing network with production facilities in North America, the U.K., China and India, among other places, Sartorius can compensate for the majority of currency fluctuations by natural hedging.

We generally hedge around two-thirds of our remaining net currency exposure over a time horizon of up to about 1.5 years through suitable currency transactions.

Business Development of Bioprocess Solutions

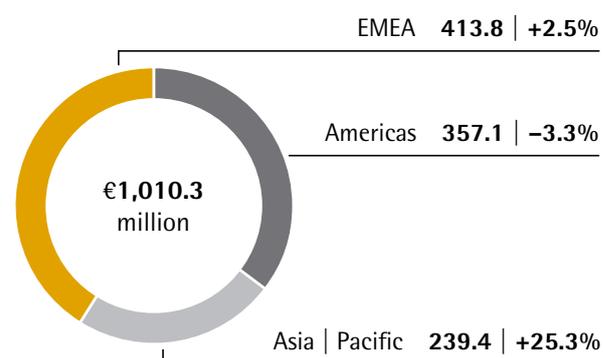
- > 4.9% sales growth relative to a very high revenue base
- > Development impacted by temporary effects in America
- > Portfolio expanded by the acquisition of Umetrics

Key Figures

€ in millions	2017	2016	Δ in %
Sales revenue	1,010.3	975.0	4.9 ¹⁾
Order intake	1,091.3	1,006.5	9.7 ¹⁾
Underlying EBITDA	282.4	273.5	3.3
as a % of sales	28.0	28.0	
Employees as of Dec.31	4,980	4,584	8.6

Sales Revenue and Growth¹⁾ by Region²⁾

€ in millions, unless otherwise specified



¹⁾ In constant currencies

²⁾ Acc. to customers' location

Products for Biopharmaceutical Manufacture



Filters for sterilization of biopharmaceutical media



Single-use bag for cell cultivation and storage of biopharmaceuticals



Single-use-based, multi-parallel bioreactor system for process development



Culture media and buffers for cell cultivation



Fully automated single-use centrifuge for cell harvesting



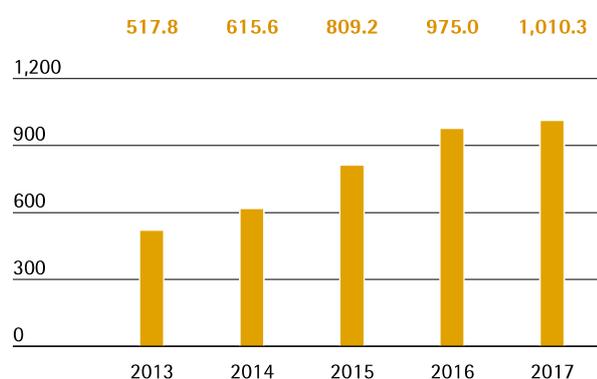
Single-use bioreactor systems for cell cultivation processes of up to 2,000 L

Sales Revenue and Order Intake

After two years of exceptionally strong growth in 2015 and 2016, the Bioprocess Solutions Division recorded sales growth in constant currencies of 4.9% to €1,010.3 million in the reporting year. Therefore, revenue was slightly lower than originally expected, but at the level of the sales forecast adjusted for the third quarter. In fiscal 2017, the division's development was impacted by a few simultaneous, temporary effects. For the year under review, the division thus reported relatively soft demand in North America and in parts of Europe, which was caused by inventory destocking by various customers. Moreover, no deliveries were meanwhile possible from the company's production facility in Puerto Rico due to Hurricane Maria, and ongoing bottlenecks in a partner's delivery capacity that lasted longer than expected impacted business performance in cell culture media. Consolidation of the acquisitions of kSep and Umetrics contributed around one percentage point in the reporting year to sales growth.

Sales Revenue 2013 to 2017

€ in millions



Sales Revenue and Order Intake

	2017	2016	in % reported	in % cc
€ in millions				
Sales revenue	1,010.3	975.0	3.6	4.9
Order intake	1,091.3	1,006.5	8.4	9.7

The division's sales development was quite different in the regions. EMEA, the region generating the highest revenue for the division and representing around 41% of its total sales, recorded a gain of 2.5% to €413.8 million. In the Americas region, which represented around 35% of revenue, sales decreased slightly by 3.3% to €357.1 million due to the temporary effects already mentioned above. The Asia | Pacific region, which accounted for around 24% of the division's sales, performed very dynamically. Partly driven by equipment shipments to fill relatively large orders, sales for this region rose significantly by 25.3% to €239.4 million.

Order intake showed positive development in the reporting year. Orders rose more strongly than division sales, and were up 9.7% in constant currencies to €1,091.3 million. On a reported basis, this increase was 8.4%.

Sales by region

	2017	2016	in % reported	in % cc
€ in millions				
EMEA	413.8	406.9	1.7	2.5
Americas	357.1	374.6	-4.7	-3.3
Asia Pacific	239.4	193.5	23.7	25.3

Overproportionate Increase in Earnings

In the reporting year, the Bioprocess Solutions Division yet again slightly increased its underlying EBITDA, which rose by 3.3% from €273.5 million to €282.4 million. The division's respective margin was at the previous year's level of 28.0%.

Underlying EBITDA and EBITDA Margin

	2017	2016
Underlying EBITDA in € millions	282.4	273.5
Underlying EBITDA margin in %	28.0	28.0

The division recorded extraordinary items of -€21.6 million in the reporting year, relative to -€20.1 million a year ago.

Product and Sales

The product portfolio of the Bioprocess Solutions Division covers virtually all steps in biopharmaceutical production process, and increasingly its upstream process development as well. The portfolio includes cell culture media, cell lines, bioreactors, a wide range of products for separation, purification and concentration, and systems for the storage and transport of intermediate and finished biological products. We also offer an extensive range of services to support our customers in complying with regulatory requirements.

Well-Developed Portfolio Supports Customers

The Bioprocess Solutions Division expanded its portfolio in the reporting year via product upgrades and new generations of existing product lines as well as through the acquisition of Umetrics.

Furthermore, Umetrics provides software solutions that enable biopharmaceutical customers to increase efficiency as well as assisting them in digitalizing their processes, especially during scale-up.

In the product area, Sartorius presented the second generation of its BIOSTAT STR single-use bioreactor series in the year under review. The second generation is designed for the new single-use bags and can handle a working volume of 12.5 to 2,000 liters. The scalable series saves customers a significant amount of time in developing production processes.

In addition, a newly developed single-use system offers users an innovative, comprehensive solution for the filtration of biopharmaceutical media on a large scale. The new system makes filtration processes even easier and less time consuming.

The area of membrane chromatography was expanded as well. Now also available as cassettes, membrane adsorbers can be combined on a large scale to form a production system for the purification of biopharmaceutical media, thus eliminating the previous limitations inherent in the smaller sizes.

Sales Activities Expanded

The Bioprocess Solutions Division markets its product portfolio directly through its own field sales representatives. Sales activities for key accounts are coordinated and supported by global key account management. As part of our regional initiatives for implementing our mid-term strategy, we stepped up our sales activities in North America during the year under review.

Research and Development

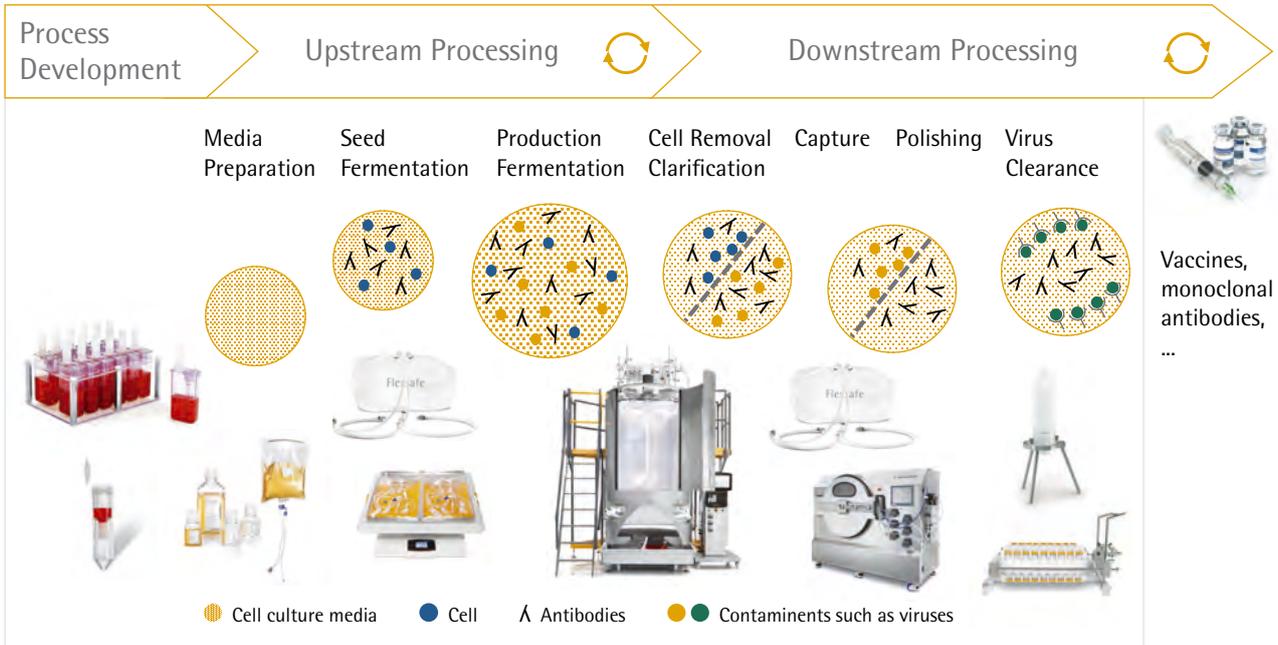
The Division's research and development (R&D) activities encompass both new and advanced in-house product developments in our own core technologies as well as the integration of new products through alliances and acquisitions. Sartorius aims to help customers continually optimize their processes and steadily increase their efficiency.

Development activities at Sartorius focus in particular on the following technology areas: membranes, which are the core component of all types of filter products; various base technologies such as single-use containers and sensors; and control technologies, for instance for fermentation.

During the reporting year, one focus of the R&D activities was the development of a new software platform for the control of individual process steps. It enables the generation of consistent data in real time for optimized monitoring and control of both upstream and downstream processes from laboratory scale all the way to commercial production.

Another key task is ongoing integration of the software portfolio acquired in the takeover of Umetrics. Even prior to the acquisition, Sartorius had already spent several years working with Umetrics data analysis specialists for the purpose of modeling and optimizing biopharmaceutical development and production processes. Additional data-based software analysis tools are planned for the future to enable customers to work more efficiently and cost-effectively.

From a regional perspective, the largest R&D site is located at our headquarters in Göttingen, with other key R&D activities taking place in Aubagne (France) Guxhagen (Germany) and Bangalore (India) as well as Bohemia (USA), Royston (UK) and Umeå (Sweden).



Simplified diagram

Production and Supply Chain Management

The Bioprocess Solutions Division operates a well developed global production network with plants in Europe, North America and Asia. The largest production sites are located in Germany, France and Puerto Rico. Moreover, Sartorius has manufacturing operations in the UK, Switzerland, Tunisia, India and the USA.

Expansion of Production Capacity

Against the backdrop of the dynamic growth of the bioprocess business in recent years, Sartorius increased and accelerated the expansion of its production capacities compared to its initial plans. Membrane production in Göttingen was further expanded.

In September, Hurricane Maria hit Puerto Rico. Thanks to the Yauco site's high construction standards, both existing buildings and those under construction suffered only minor damage. Heavy infrastructure damage made deliveries from the plant impossible for around four weeks, however. Even so, Sartorius considers the Yauco location to be highly suitable. In the future, however, production will focus more strongly on standard products that can be stored on the U.S. mainland to reduce possible future disruptions in the supply chain.

Business Development of Lab Products & Services

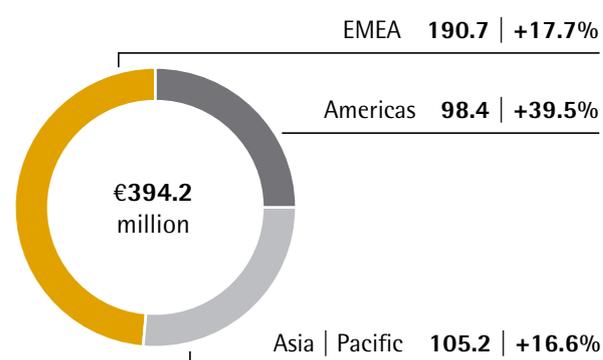
- > Exceptionally dynamic business development
- > Strong growth in all product areas and regions
- > Bioanalytics portfolio extended by the acquisition of Essen BioScience

Key Figures

€ in millions	2017	2016	Δ in %
Sales revenue	394.2	325.3	22.0 ¹⁾
Order intake	410.1	328.1	25.8 ¹⁾
Underlying EBITDA	70.8	51.9	36.3
as a % of sales	18.0	16.0	
Employees as of Dec.31	2,521	2,327	8.3

Sales Revenue and Growth¹⁾ by Region²⁾

€ in millions, unless otherwise specified



¹⁾ In constant currencies

²⁾ Acc. to customers' location

Products for Laboratories in Quality Assurance and Research



Laboratory balances for reliable and convenient weighing processes



Syringe filters for analytical sample preparation and sterile filtration



The lightest and smallest electronic pipette on the market



Real-time live-cell imaging system for cell analysis



Filter membranes for microbiological testing in the lab



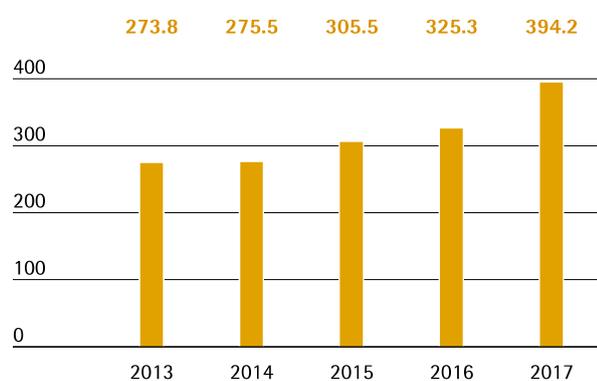
Cell screening platform that has already won several awards

Sales Revenue and Order Intake

In the reporting year, the Lab Products & Services Division recorded a substantial increase in sales revenue by 22.0% to €394.2 million. The division's portfolio expansion into bioanalytics by the acquisition of Essen BioScience at the end of March 2017, as well as of IntelliCyt and ViroCyt in 2016 contributed close to 14 percentage points to this increase. Driven by strong demand in all regions and in all product areas, organic growth for the division was around 8%. Therefore, sales development was within the range of the company's forecast that was adjusted in April 2017 due to its acquisitions.

Sales Revenue 2013 to 2017

€ in millions



Sales Revenue and Order Intake

€ in millions	2017	2016	in % reported	in % cc
Sales revenue	394.2	325.3	21.2	22.0
Order intake	410.1	328.1	25.0	25.8

In EMEA, the region contributing the highest share of revenue of around 48%, the division achieved a 17.8% increase in sales to €190.7 million. The Americas region, whose share of revenue rose to around 25%, reported the highest growth dynamics due to the acquisition of Essen BioScience, recording an uptick of 39.5% in sales to €98.4 million. Besides the division's excellent organic performance, the acquisitions previously mentioned contributed an above-average proportion to the region's growth. The Asia | Pacific region, which accounts for a good 27% of business for the Lab Products & Services Division, reported a double-digit gain of 16.6% to €105.2 million.

During the reporting period, order intake showed especially encouraging development and grew to €410.1 million at a considerably stronger rate, 25.8%, than did the division's sales revenue. On a reported basis, this growth rate was 25.0%.

Sales by region

€ in millions	2017	2016	in % reported	in % cc
EMEA	190.7	162.2	17.6	17.7
Americas	98.4	70.8	39.0	39.5
Asia Pacific	105.2	92.3	14.0	16.6

Earnings

The Lab Products & Services Division increased its underlying EBITDA in the reporting period by 36.3%, from €51.9 million a year ago to €70.8 million, which was due to economies of scale in connection with the division's strong organic growth, as well as to growth contributed by its acquisitions. Despite unfavorable currency effects, the division's corresponding margin improved from 16.0% in the previous year to 18.0%. This increase was likewise attributable to economies of scale related to strong organic growth as well as growth contributed by acquisitions.

Underlying EBITDA and EBITDA Margin

	2017	2016
Underlying EBITDA in € millions	70.8	51.9
Underlying EBITDA margin in %	18.0	16.0

The division recorded extraordinary items of -€13.9 million in the reporting year relative to -€10.6 million a year ago.

Products and Sales

The Lab Products & Services Division focuses on both research laboratories and quality assurance laboratories in the pharmaceutical and biopharmaceutical sectors, as well as on academic research institutes.

Significant additions were made to the division's portfolio in the area of bioanalytics in the year under review. The remaining product portfolio is used especially for sample preparation and other standard applications, and includes instruments such as laboratory balances, pipettes and bioreactors, in addition to consumables like laboratory filters and microbiological tests. Extensive services round out the portfolio.

The division sells its products through three channels: laboratory dealers, direct sales and e-business. The dealer sales channel is already well established; direct sales and the digital channels are being expanded continually.

Expanded Bioanalytics Portfolio

Sartorius significantly extended its expertise in the field of bioanalytics through the acquisition of Essen BioScience, which further improves the company's positioning among biopharmaceutical customers and life science research institutes in particular. The acquisition will lead to growing synergies between the Lab Products & Services and Bioprocess Solutions divisions.

Essen BioScience offers a platform consisting of instruments, software and reagents for fully automated, real-time analysis and visual depiction of biological processes using high-resolution imaging. The data obtained provides new insight and understanding into the mechanisms of healthy and diseased cells, which helps to significantly accelerate the often time-consuming development process for new substances.

Range of Services and Test Laboratories in All Regions

The Lab Products & Services Division offers a wide range of services covering the entire life cycle of laboratory products, from device installation and commissioning to verification, calibration, regular maintenance and repair. Our services are not limited to Sartorius instruments; they are offered for devices from other manufacturers as well. This extensive range enables our customers to minimize the number of service providers they use and to reduce complexity and costs.

Our application laboratories in all regions offer customers the opportunity to test Sartorius products, even using their own samples, and to take training courses.

Growing Importance of Bioanalytics

Sartorius enjoys a strong presence in the lab market in essential product areas, but for historical reasons, its market share varies by product group and region. This situation offers significant potential for growth. In particular, the company's products in the area of bioanalytics offer strong unique selling points. We aim to use our global sales platform to accelerate the penetration of these technologies across the laboratory market.

However, ongoing improvements in sales efficiency are just as important as expanded sales structures. The Lab Products & Service Division is also benefiting from the customer access established for biopharmaceutical customers by the Bioprocess Solutions Division.

Research and Development

Within the scope of its laboratory business, Sartorius has extensive technological expertise in the area of laboratory instruments and laboratory consumables. A further technology focus is bioanalytics.

Most of the division's R&D takes place centrally at the Göttingen site, but additional R&D activities are pursued at our facilities in Helsinki, Stonehouse, Ann Arbor and Albuquerque.

Products and services from Sartorius help users make laboratory processes more productive, efficient as well as safer. Compliance with regulatory requirements is critical. Product development priorities for Sartorius therefore include data management, connectivity and process automation.

As part of its increasing focus on the biopharmaceutical sector, Sartorius develops weighing technology and equipment optimized for pharmaceutical applications. In further development stages, its instruments are specially adapted to these applications, for example, by combining a comparatively high weighing capacity and exceptionally high accuracy.

For example, special software routines already guide laboratory employees through work processes and transfer measurement and process data to laboratory management systems. Intuitive user interfaces in laboratory instruments enable safer and faster work.

Production and Supply Chain Management

The Lab Products & Services Division operates plants in Germany, China, Finland, the U.K. and the U.S. The plants serve as centers of competence and tend to focus on one product group or a small set of product groups. Laboratory balances are manufactured in Göttingen and Beijing, for example, while pipettes are made at our Helsinki and Suzhou sites. Microbiological tests kits are produced in Stonehouse, and most membrane-based products are manufactured in Göttingen.

Assessment of the Economic Position

The Sartorius Group's Business Performance

Sartorius made further progress in fiscal 2017 and grew profitably yet again. In line with the ongoing implementation of Sartorius' strategic plan, the company expanded its product portfolio during the year under review through two acquisitions. The business of the Bioprocess Solutions Division was extended in the area of data analysis through the acquisition of Umetrics, and the Lab Products & Services Division strengthened its position in the high-growth segment of bioanalytics by the takeover of Essen BioScience. Due to consolidation of these acquisitions, Sartorius raised its forecast in April 2017 for the current fiscal year.

The Lab Products & Services Division saw exceptionally dynamic growth due to its strong organic development and to the expansion of its portfolio into bioanalytics. All regions and product areas contributed to this strong growth.

Business development of the Bioprocess Solutions Division was slightly lower than we had originally expected, especially in the second half, due to several simultaneous and temporary effects. Primarily impacting the Americas region, these were temporary bottlenecks in supply for the division's business with cell culture media, inventory destocking by a few large customers and an interruption in production at the plant in Puerto Rico following Hurricane Maria.

Because of the temporary effects previously mentioned, Sartorius adjusted its annual forecast for the Bioprocess Solutions Division and the Group in October 2017. The company's forecast for the Lab Products & Services Division remained unchanged.

Consolidated sales revenue rose during 2017 by 9.3% in constant currencies to €1,404.6 million. The Group's earnings margin, based on underlying EBITDA, increased to 25.1%, also meeting our expectations adjusted in the third quarter.

The ratio of net debt to underlying EBITDA for the last twelve months stood at 2.5 as of December 31, 2017 (2016: 1.5), which was in the range of our adjusted forecast. As a result, the Sartorius Group continues to have significant financing flexibility to further implement its strategy.

In the context of its 2020 strategy, Sartorius is making above-average investments in the expansion of its manufacturing capacities, its IT systems, and in the consolidation and expansion of its Group headquarters in Göttingen, Germany. Capital expenditures accordingly rose in the reporting year from €152.1 million to €209.4 million.

Projected | Actual Comparison for the Year 2017

	Actual 2016	Guidance January 2017	Guidance July 2017	Prognose October 2017	Actual 2017
Sartorius Group					
Sales growth ¹⁾	18.2%	~8% - 12%	~12% - 16%	~9%	9.3%
Underlying EBITDA margin in %	25.0%	~+0.5% ¹⁾	~+0.5% ¹⁾	~25% ³⁾	25.1%
Gearing (underlying)	1.5	< 1.5 ²⁾	~2.4	~2.4	2.5
CAPEX ratio	11.7%	~12% - 15%	~12% - 15%	~12% - 15% (upper end)	14.9%
Sartorius Sparten					
Bioprocess Solutions					
Sales growth ¹⁾	22.1%	~9% - 13%	~9% - 13%	~4%	4.9%
Underlying EBITDA margin in %	28.0%	~+0.5% ¹⁾	~+0.5% ¹⁾	~28% ³⁾	28.0%
Lab Products & Services					
Sales growth ¹⁾	7.9%	~6% - 10%	~20% - 24%	~22%	22.0%
Underlying EBITDA margin in %	16.0%	~+1% ¹⁾	Close to +2pp	~18% ³⁾	18.0%

¹⁾ In constant currencies

²⁾ Possible acquisitions are not considered

³⁾ In actual currencies

Annual Financial Statements of Sartorius AG

The retained profit of Sartorius AG is the key reference value for the payment of dividends to our shareholders. Whereas the Sartorius Group financial statements are drawn up according to the International Financial Reporting Standards (IFRS), the annual financial statements for Sartorius AG are prepared by applying the rules and regulations of the German Commercial Code (HGB).

The Management Report of Sartorius AG and the Group Management Report for fiscal 2017 are combined. The HGB annual financial statements of Sartorius AG and the combined management report are published simultaneously in the German Federal Gazette (Bundesanzeiger).

Business Operations, Corporate Strategy, Corporate Management and Oversight, Overview of Business Development

Sartorius AG has exercised only the functions of the strategic, group-leading management holding entity for the Sartorius Group since the beginning of fiscal 2011, and we refer in this connection to the explanatory reports concerning business operations, corporate strategy, corporate management and oversight, as well as the overview of business development, presented on pages 21 et seq. of the combined management report of Sartorius AG and the Group.

Earnings

Sales revenue of Sartorius AG essentially consists of cost transfers to affiliated companies within the Group for management services rendered as well as of the rental of buildings on Sartorius Campus.

Income from investments of €28.7 million relative to €22.8 million a year ago concern dividends paid out for the French subsidiary Sartorius Stedim Biotech S.A.

The total loss accepted under profit and loss transfer agreements was €40.0 million. Of this amount, €0.1 million was transferred to Sartorius Corporate Administration GmbH (2016: profit of €1.4 million) and €39.9 million to Sartorius Lab Holding GmbH, compared with a profit of €4.1 million that the latter had accepted in 2016.

The loss incurred by Sartorius Lab Holding GmbH essentially resulted from the non-cash valuation of a U.S. dollar-denominated loan needed for acquisitions executed in the USA.

Net Worth and Financial Position

The balance sheet total of Sartorius AG rose in the reporting year by €384.1 million to €1,225.3 million. This increase was primarily due to the acquisitions executed in the USA and investments made at the Göttingen site.

The balance sheet structure of Sartorius AG reflects its function as the management holding entity for the Sartorius Group. Fixed assets consist essentially of financial assets and, in the reporting year, stood at €507.4 million (2016: €474.8 million). Accordingly, fixed assets accounted for 41.4% of the balance sheet total (2016: 56.4%). The equity ratio for fiscal 2017 was 22.5% relative to 38.4% in 2016.

Statement of Profit and Loss of Sartorius AGBased on the total cost accounting method according to Section 275, Subsection 2, of HGB¹⁾

€ in K	2017	2016
1. Sales revenue	6,930	6,023
2. Other operating income	9,833	3,226
3. Employee benefits expense	-5,611	-5,108
4. Depreciation and amortization	-1,116	-879
5. Other operating expenses	-14,129	-7,228
6. Income from investments	28,749	22,817
7. Profit received under a profit and loss transfer agreement	0	5,569
8. Loss accepted under a profit and loss transfer agreement	-39,979	0
9. Interest and similar income	11,252	7,449
10. Interest and similar expenses	-12,964	-7,198
11. Income tax expense	-1,000	-2,135
12. Profit after tax	-18,035	22,536
13. Other taxes	-23	-23
14. Net profit for the period	-18,058	22,513
15. Profit brought forward	159,923	168,527
16. Retained profits incl. net profit for the period	141,865	191,040

¹⁾ HGB = German Commercial Code**Balance Sheet of Sartorius AG**According to HGB¹⁾, € in K

	Dec. 31, 2017	Dec. 31, 2016
Assets		
A. Fixed Assets		
I. Property, plant and equipment	63,731	33,190
II. Financial assets	443,622	441,622
	507,353	474,812
B. Current Assets		
I. Trade and other receivables	713,536	363,981
II. Cash on hand, deposits in banks	2,099	278
	715,635	364,259
C. Prepaid Expenses	2,298	2,109
	1,225,286	841,180

	Dec. 31, 2017	Dec. 31, 2016
Equity and Liabilities		
A. Equity		
I. Subscribed capital	74,880	74,880
Nominal value of treasury shares	-6,492	-6,492
Issued capital	68,388	68,388
II. Capital reserves	54,065	52,823
III. Earnings reserves	10,867	10,867
IV. Retained profits incl. net profit for the period	141,865	191,040
	275,185	323,118
B. Provisions	27,559	30,937
C. Liabilities	922,542	487,125
	1,225,286	841,180

¹⁾ HGB = German Commercial Code

Proposal for Appropriation of Profits

The Executive Board will submit a proposal to the Annual Shareholders' Meeting to appropriate the retained profit of €141,864,793.40 reported by Sartorius AG for the year ended December 31, 2017, as follows:

	in €
Payment of a dividend of €0.50 per ordinary share	17,106,112.00
Payment of a dividend of €0.51 per preference share	17,429,794.68
Unappropriated profit carried forward	107,328,886.72
	141,864,793.40

Research and Development

Detailed information about the research and development activities of the Sartorius Group and of its divisions is given on pages 42 et seq. and 46 et seq.

Employees

Sartorius AG as a holding company does not employ any staff to be disclosed pursuant to Section 285, No. 7, of HGB.

Risks and Opportunities

The opportunities and risks affecting the business development of Sartorius AG as the management holding entity are essentially equivalent to those of the Sartorius Group. Sartorius AG shares in the risks to which its investments and subsidiaries are exposed in proportion to the extent of its investment. Where expedient and feasible, we adopted countermeasures and/or arranged for balance sheet measures during the reporting year to cover all discernible risks within Sartorius AG that had the potential to negatively impact our net worth, financial position and profitability.

A detailed Opportunity and Risk Report for the Sartorius Group is provided on pages 54 to 62; a description of the internal control and risk management system, on pages 67 to 68.

Report on Material Events

Please refer to page 64 to read the Report on Material Events for Sartorius AG and the Sartorius Group.

Forecast Report

Earnings trends for Sartorius AG depend substantially on the progress of its subsidiaries and, hence, on the Sartorius Group.

The development of the Sartorius Group's business is discussed in the Forecast Report on pages 63 to 65.

Opportunity and Risk Report

Principles

Every business activity entails opportunities and risks, which have to be managed. The skill with which this is done goes a long way in determining the future development of a company's shareholder value. The central element of risk management is systematic identification and realization of opportunities, as well as avoidance of risks that could jeopardize the success of the company.

In managing risks and opportunities, Sartorius aims to identify and use business opportunities systematically, as well as to recognize and evaluate risks at an early stage and take measures to counter them where possible. It is not the task of risk management to eliminate all risks: rather, our approach is to intentionally take a certain measure of risk in our business activities in order to be successful in unlocking opportunities. However, in this endeavor, it is important to keep risks contained within acceptable limits and to control them carefully. Through appropriate guidelines, we ensure that risk assessments are taken into account in the decision-making processes from the very beginning.

Sartorius has no single organizational unit tasked with identifying and managing opportunities and risks. Instead, it prefers to make this a cross-functional component of Group management. To this extent, our risk management organization reflects a global functional matrix organization in which individuals heading a functional area, so-called "risk owners," are each responsible for their own management of opportunities and risks. The Finance & Controlling department is responsible for the organization of the respective reporting process, including the further development of the Group's risk management system.

Managing Opportunities

Our opportunity management centers on the analysis of target markets and sector environments, as well as the assessment of trends, both of which give strong indications as to future business opportunities. The identification of the potential for development in this context is one of the key roles of the relevant managers and initially takes place at the local rather than the central level. The market-facing functions, such as strategic marketing and product management in the individual divisions, play a leading role in this respect. The central Business Development unit additionally supports these areas with market monitoring, data analysis and the implementation of strategic projects.

As part of strategy reviews, the members of the Group Executive Committee regularly meet with the managers having operational responsibility and with the Business Development unit to discuss short-, medium- and long-term opportunity potential for the various business areas. The subsequent steps of prioritizing the opportunities and evaluating them from a business management perspective, deriving strategic measures and allocating resources proceed in accordance with a standardized decision-making process that applies throughout the Group. If the opportunities are short-term in nature, they are considered in annual budget planning. Medium- and longer-term opportunities are tracked systematically as part of strategic planning. The status of opportunity management as a permanent fixture of the corporate management system means that it also features in the discussions and decision-making processes of top-level management, such as the Executive Board and the Supervisory Board.

Key areas of opportunity are presented below. Where appropriate, reference is made to the relevant section of the Group Management Report in order to avoid repetition. Most of the risks we describe in the section on specific risks represent opportunities, should events develop in the opposite, positive direction. For this reason, we discuss these opportunities in the section on specific risks and opportunities at the end of this chapter.

Areas of Opportunity

As a supplier for the pharmaceutical and laboratory industries, Sartorius operates in future-oriented and high-growth sectors. The significant opportunities generated by the various market and technology trends are described in detail in the sections entitled "Sector Conditions" and "Outlook for the Sectors" on pages 29 and 63, respectively.

Our assessments rank the company as one of the global market leaders in many subsegments and product areas. We believe the high quality of our products, our strong brand recognition and our established customer relationships give us strong opportunities to continue extending our market leadership. The corresponding division strategies and the growth opportunities and initiatives based on them are discussed in the section on the strategy of the Bioprocess Solutions Division, which begins on page 25, and in the section on the strategy of the Lab Products & Services Division, which starts on page 26.

Strict management of processes and costs provides opportunities to further increase our profitability. Key target areas in this respect include continued enhancements of our procurement chain and ongoing efforts to optimize production, which we present on pages 45 and 49.

Other opportunities are discussed in the context of the presentation of specific risks and opportunities beginning on page 56.

Risk Management

Just as for opportunity management, overall responsibility for the maintenance of an effective risk management system ensuring comprehensive and consistent management of all material risks rests with the Executive Board. Coordinating and developing this system and combined risk reporting are the responsibilities of the Finance & Controlling department, while the particular functional areas are responsible for identifying and reporting risks, as well as for assessing their potential impact and for taking the appropriate countermeasures.

The Supervisory Board of Sartorius AG monitors the effectiveness of the risk management system, with the preparatory work necessary for it to do so being performed by the Audit Committee of this board. Furthermore, while carrying out their statutory audit mandate for the annual financial statements and consolidated financial statements, the independent auditors examine whether the early warning system in place is capable of prompt identification of risks that could jeopardize the future of the company.

Risk Management System and Risk Reporting

At the heart of the risk management system is the Sartorius Group Risk Management Handbook, which applies throughout the entire Group organization. The Handbook, which includes definitions of the framework, the structural organization, processes, risk reporting and monitoring and control of the effectiveness of the risk management system, is based on the internationally recognized COSO standard. There are also a number of other sources that contain stipulations for handling risks, including the articles of association and rules of procedure of the Group companies and other internal guidelines.

The Group-wide risk reporting system forms the cornerstone of internal risk communication. The object is to make it possible to address risks in a structured, continuous manner and to document them in accordance

with the relevant statutory and regulatory requirements. The strong growth of the Group over the past years and the rising demands of customers and regulators meanwhile require that we continue to adapt our guidelines and rules.

A key element of our internal communication of risks is Group-wide risk reporting. The objective of this is to enable structured, continuous tracking of risks and to document them in compliance with legal and regulatory requirements.

The prescribed reporting process in the risk categories subsequently described establishes the rules for the ongoing review of and information on risk situations. Those responsible for functional areas at the Group subsidiaries continually review and assess their respective risk situations. If any specific risks are discernible, these are documented with respect to their assessment, probability of occurrence and measures to be taken to eliminate such risks or to mitigate their impact. In addition, as soon as these risks reach defined size criteria, they are reported to the central risk management system. Appropriate insurance policies are taken out to reduce any remaining risk situations, where feasible. New organizational units joining the consolidated Group companies are successively integrated into our risk reporting process.

We have an urgent reporting procedure in place to ensure that when a new or emerging significant risk to our net worth, financial position and profitability is identified and estimated as involving €2.5 million or higher, the Executive Board of Sartorius AG receives all of the necessary details without undue delay.

Risk Classification

The first level of risk management relates to the four main risk categories defined by Sartorius: external risks, operating risks, financial risks and corporate governance risks.

The second level consists of additional subcategories that we classify within these main categories, such as supply chain risks, sales and distribution risks, and quality risks.

We categorize risks according to the scale of their implications too, and also perform a specific evaluation in which all risks are assigned the value of their maximum impact at the time of risk analysis. In other words, we record the maximum risks without considering the probability of occurrence or the effects of risk mitigation measures.

For the purposes of this report, we have assessed the probability of the risks as shown below and, in the adjacent columns, classify their particular significance for the entire Group.

Probability of Occurrence	
Low	< 5%
Medium	5% - 20%
High	> 20%

Significance

in thousands of €	Impact on Earnings*	Impact on Assets*
Of limited significance	< 1,000	< 5,000
Significant	> 1,000	> 5,000

Explanation of Principal Risks and Opportunities

General and Macroeconomic Risks and Opportunities

General Risks

In principle, our ability to foresee and mitigate the direct and indirect effects of risks entailed by life in general is limited, but we proactively take measures, whenever feasible, to ensure that we can respond appropriately and at short notice or are insured against any damage entailed by such risks that include, for instance, force majeure, natural catastrophes and their associated damage to commercially significant and critical infrastructure. After severe hurricanes hit Puerto Rico in the past fiscal year, we deployed a task force that defined the immediate measures to take, thus limiting the economic damage to the Group. At the same time, a project for updating our production network strategy was launched.

Political developments, such as the referendum on the United Kingdom's leaving the European Union or the change in politics in the USA, can basically have an impact on the Group's business. To cite such examples, such developments may involve changes to the tax system or customs duties, as well as impacts on the exchange rate of the euro to the British pound or the U.S. dollar (for more on the subject of exchange rates, see the section below on Exchange Rate Risks and Opportunities). These further developments are being closely observed so that measures can be taken to reduce such risks, as necessary.

Business Cycle Risks

The nature of our various business areas means that Sartorius as a whole is insulated to a certain extent from the full force of wider cyclical effects. While macroeconomic effects have very little bearing on the progress of the Bioprocess Solutions Division, they are particularly relevant to the Lab Products & Services Division and can represent a risk to the latter's growth in this context. If economic developments prove more positive than expected, this, in turn, can additionally stimulate stronger growth for Lab Products & Services.

Supply Chain Risks and Opportunities

Our supply chain extends from procurement to production to sales and distribution. Problems within this workflow can have consequential effects, including delays in deliveries. The global supply chain management system we have instituted throughout our production processes to prevent such problems largely minimizes the associated risks by analyzing and controlling all of the operations involved. The strongly international alignment of our organization opens up a whole series of opportunities too. The various risks and opportunities encountered within our supply chain are explained in detail below.

Procurement Risks and Opportunities

We purchase a wide range of raw materials, components, parts and services from suppliers and are consequently exposed to the risks of unexpected delivery bottlenecks and | or price increases.

Over the past years, we have implemented powerful tools and robust processes in our Materials Management unit to manage risks arising from critical materials. These means enable us to meet the needs of our customers with respect to assurance of supply and transparency. This represents a significant competitive advantage. Important measures in this respect are to maintain security stock and to define alternative suppliers as part of our second supplier policy.

In addition, we conduct regular supplier reviews and also use early warning systems.

Risks arising from raw material prices play a rather subordinate role in our business. On the one hand, the proportion of raw materials in our production costs is comparatively low. On the other hand, we purchase a wide range of materials so that price increases for certain materials do not represent any significant factor.

Opportunities can arise in the area of procurement when our growth enables us to increase order quantities and thereby strengthen our position with our suppliers, such as by receiving price discounts or preferential treatment as a "preferred customer." In addition, we maintain a list of preferred suppliers in parallel, which permits us to enter into long-term business relationships with key suppliers to our mutual benefit.

Increased globalization of our supplier pool holds the prospect of purchasing on more favorable terms, moreover, and there is also a possibility of our expanded purchasing activities in the international markets leading us to identify suppliers with special product and technical expertise that could eventually enhance our own competitive edge.

Production Risks and Opportunities

Based on our core technology expertise, we ourselves manufacture a large proportion of the products that involve a high level of vertical integration. Examples include filters and laboratory balances. Other products, such as reusable fermenters and bioreactors, are manufactured in collaboration with suppliers so that some of the production risks are transferred to external third parties. When we manufacture products ourselves, we also bear the associated risks of capacity bottlenecks or overcapacity, production downtimes, excessive reject rates and high levels of tied-up working capital, as well as dependency on individual manufacturing sites. We contain and reduce these risks by planning production capacities carefully, using versatile machines, semi-automated individual workstations and flextime work schedules, and by continuously monitoring production processes. Moreover, our global manufacturing network enables us to compensate for any capacity bottlenecks by shifting production to other regional plants and to minimize our dependency on individual local manufacturing sites.

Furthermore, we have taken out policies for business interruption insurance to compensate for any possible losses due to production downtimes.

Beyond this, we work closely together with our customers to gain a better understanding of their needs and to schedule our production capacities optimally.

We consider it an opportunity that our investments in infrastructure and production resources, among other things, have given us high flexibility in our manufacturing operations and that we are capable of meeting our customers' requirements and regulatory standards with respect to business continuity concepts. In addi-

tion, this approach ensures that our individual production sites can concentrate on specific manufacturing technologies, gaining added efficiency as a result. Our international manufacturing network also makes it possible to capitalize on the cost advantages offered by individual sites. Furthermore, continuous improvements in production, such as simplifying processes and increasing levels of automation, help drive manufacturing efficiency even higher.

Over the past years, we have substantially enhanced our production workflows at our largest production facility at Göttingen corporate headquarters by consolidating individual manufacturing sites at Sartorius Campus and by optimizing process sequences when operations were started up at the new laboratory production building, thus minimizing the associated risks.

Sales and Distribution Risks and Opportunities

We use a variety of channels to sell and distribute our products around the world. The potential risks entailed are unexpected changes in the demand structure, growing price pressure and non-compliance with supply agreements concluded with customers. We employ targeted market analyses to identify emerging demand trends in individual segments early on so that we have time to respond appropriately. Our technical innovations and our focus on less price-sensitive sales markets, such as products for validated production processes in the biopharmaceutical industry, reduce our exposure to the risk of growing price pressure. We have minimized our risk exposure in the area of logistics in recent years by setting up and using central warehouses to optimize distribution logistics.

Opportunities arise in the area of sales and distribution when the increasing breadth of our product range – in both bioprocess and lab segments – puts us in a position to sell new products to existing customers. Our business relationships, most of which are established for the long term, and our global presence provide opportunities, moreover, and our ongoing project to strengthen direct sales, especially in the lab segment, also promises to enhance our sales prospects.

With our acquisitions in cell analytics, we have strengthened the synergies between both divisions and offer our biopharmaceutical customers comprehensive solutions for research laboratories all the way to production processes.

Quality Risks and Opportunities

Our customers use Sartorius products in a wide range of critical production processes, including the manufacture of pharmaceuticals, foods and chemicals, and in research and development laboratories. The main risk encountered in these areas is non-compliance with agreed quality criteria, which can lead to losses for our customers, or their customers, for which we may be made liable through compensation claims. We employ rigorous quality checks and advanced production methods and processes, such as cleanroom technology, to ensure that our products satisfy the most stringent quality standards and high regulatory requirements. These manufacturing methods and processes are subject to constant review under our continuous improvement processes, moreover, and are optimized as requirements evolve. Our successful completion of a host of annual audits by customers and our certification under ISO 9001 and ISO 13485 together document the high level of quality achieved in Sartorius products and processes. Irrespective of these measures, we also maintain significant insurance coverage against product liability risks. Sartorius has established a traceability system that enables us to recall an entire production batch immediately, if necessary, and minimize any adverse consequences in the event of defects being discovered in a product.

We have installed a complaints management system to deal with customer requests and to ensure full documentation.

Quality requirements are growing more and more stringent all the time, not least as a result of increasing requirements on protection of medical patients and on product safety by regulatory authorities, so we regard this first and foremost as an opportunity that opens up new market prospects. Also, challenging quality demands represent a considerable barrier to entry for potential new competitors and provide stimulus for further technical innovation to which we actively respond.

R&D Risks and Opportunities

We devote a considerable share of our resources to research and development. Potential risks in this area may arise from development results that diverge from market needs and application requirements and from exceeding planned development deadlines. Our advanced project management, intensive R&D controlling and early involvement of our customers in the development process substantially limit these R&D risks. Patents and continuous tracking of the technologies

and competitors relevant to us secure our technology and marketing position.

On the other hand, the R&D sphere also offers a number of opportunities. Our intensive collaboration with partners that rank among the global market leaders in their own fields opens up the opportunity for us to jointly develop products with an especially high level of innovation. In areas such as membrane technology and plastics technology, sensorics and biopharmaceutical process engineering, as well as weighing technology, in turn, the expertise of our own specialists puts us at the very forefront of global research and development, presenting us with an opportunity to turn this technical knowledge into potential sales and an even stronger position on the market.

Customer Risks and Opportunities

Sartorius sources its key customers from the pharmaceutical, chemical and food industries and from research and educational institutions of the public sector. These customers are usually relatively large organizations that have been in existence for some time and have strong credit ratings. Most of our business areas have a highly diversified customer base, so the Group as a whole is not dependent on individual key accounts to any significant degree.

Competitive Risks and Opportunities

Sartorius has a leading competitive position in most of its markets and competes with mainly larger rivals sharing our status as a globally operating company. As we serve a large number of customers from highly regulated sectors like the pharmaceutical and food industries, and the technology barriers to market entry are substantially high, we regard the probability of new competitors emerging within the short term as low. Overall, the competitive environment is therefore very stable.

The fact that many of our products are used in validated processes, especially those in the biopharmaceutical industry, ensures a robust revenue stream, reducing the risk of losing significant market share within a short timeframe. At the same time, it is not possible for us to quickly force out the competition that serves customers in this area.

Changes in the competitive environment, for example, consolidation in the markets, can pose opportunities. Our sectors find themselves in an ongoing process of change in which Sartorius has been actively participating. We have been continuously making acquisitions in recent years to reinforce our market position and open up new potential synergies.

Acquisition Risks and Opportunities

By nature, acquisitions provide many opportunities, such as sales growth, extension of our product portfolio and development of new markets. By contrast, the purchase and sale of companies or parts of companies entail a number of typical risks, such as incorrect valuation assumptions or insufficient usage of anticipated synergy effects. To prevent these risks, we take various measures, such as performing a standard due diligence review of important areas and carrying out comprehensive analysis of the market concerned. In addition, we involve external consultants and experts in the purchase or sales process as required. We especially focus on drafting transaction contracts so that they adequately counter such risks, especially by clauses assuring specific characteristics or by contractual warranty or guarantee provisions, as well as agreements on mechanisms for adjustment of the purchase price and on liability clauses or by taking out the appropriate insurance policies. Immediately after an acquisition has taken place, an integration phase is initiated in which any potential risks can likewise be detected as early as possible and prevented or minimized by taking the appropriate counteractions.

Personnel Risks and Opportunities

As an innovative technology group, Sartorius employs a large percentage of highly qualified people. We counter the risks of a possible scarcity of required specialists, especially those in key positions, and of demographic change by offering performance-related remuneration models, targeted continuing professional development options, further attractive social benefits, continuous education and training for junior staff members within our organization and interesting people development opportunities.

The success of these measures is apparent in the low attrition rates of recent years and the many years of seniority our people accumulate on average. Employment contracts in certain cases contain a clause prohibiting any move to a direct competitor.

Opportunities for Sartorius primarily arise in that it can further qualify its staff by offering its own training courses and retain such staff over the long term, thus covering company needs for qualified personnel particularly well.

Financial Risks and Opportunities

The global nature of the Sartorius Group's operations means that its business activities are inevitably exposed to financial risks. The most significant of these, aside from risks associated with Group accounting, are exchange rate risks, interest rate risks and liquidity risks, all of which are described below and addressed in detail in the Notes to the Consolidated Financial Statements. Vice versa, financial risks, most notably exchange rate risks and interest rate risks, are balanced by opportunities of approximately equal magnitude.

Risks Associated with Group Accounting

Except for the general, typical risks inherent in any accounting process, no specific risks concerning Group accounting are discernible. Typical accounting errors in this connection are, for example, incorrect discretionary decisions in the measurement of assets and liabilities. The use of various standardized, partially automated control mechanisms largely ensures that such errors are recognized and corrected at an early stage. A highly detailed monthly reporting and analysis process, as well as regular legal audits, essentially contributes to these targets.

Exchange Rate Risks and Opportunities

As we generate around half of consolidated sales revenue in foreign currencies and two-thirds of this total revenue in U.S. dollars or in currencies pegged to the U.S. dollar, we are positively or negatively impacted by currency effects, especially when converting the currencies of balance sheet items and profit or loss items, respectively. To largely compensate for the general risk resulting from the impact of individual foreign currencies, we have taken a number of measures besides hedging currencies. Our global production network thus enables us to offset the lion's share of sales revenues received in foreign currency within the Group against costs likewise incurred in foreign currency. For example, we manufacture many of our products for the North American market locally, and are not disadvantaged in competition with our U.S. rivals, insofar as this general currency risk is concerned. We continuously monitor both exchange rates and our net cur-

currency exposure – i.e. that proportion of our foreign currency sales revenue that remains after we have settled our costs, likewise in a foreign currency – and use derivative financial instruments for hedging. These instruments are primarily spot, forward and swap transactions, on the basis of current and anticipated net currency exposure and foreign currency levels. We make it a policy to hedge up to 70% of our exposure in advance for the following 18 months. Hedging transactions are set up by one group of staff and monitored by another, separate group.

Interest Rate Risks and Opportunities

We have concluded fixed interest agreements for approximately two-thirds of our outstanding loans to eliminate the risk posed by variable interest payments. The remaining portion of the financial instruments outstanding on the reporting date is subject to variable interest based on the market rate. Almost one-third of these is currently covered by interest rate swaps. We monitor interest rate trends and our interest rate exposure constantly and have the facility to arrange for additional hedging transactions where we consider it necessary and economically advisable to do so for individual loans.

Liquidity Risks and Opportunities

The Sartorius Group actively manages liquidity centrally in order to check and minimize liquidity risks and optimize liquidity management within the organization. For this purpose, we use various long- and short-term financial instruments. Regarding the maturities of our loans, we make it a policy to take a risk-averse approach.

For short-term liquidity procurement, we also employ various instruments. In addition to the syndicated credit line that can be accessed and repaid at short notice, we have a number of bilateral credit lines in place on a smaller scale for individual Group companies. Furthermore, we use cash pooling agreements between selected Group companies as the primary tool to manage liquidity within the Group.

Under the present loan agreements, the Group is required to comply with standard financial key ratios, or covenants. In this context, the ratio of net debt to underlying EBITDA may not be greater than 3.25 and 4.00, respectively. There is a basic risk in this connection that the Group might not comply with these covenants, which could lead to a termination of the loan agreements. Based on the current information available, this is considered unlikely.

Regulatory Risks

Our role as a supplier to the biopharmaceutical industry and health care providers means that Sartorius can also be affected by underlying developments in these areas. The possibility of the regulatory authorities (FDA, EMA) adopting a more restrictive approach to the approval of new medications remains the principal source of risk in this context. Such a move would reduce the number of new pharmaceutical products to be marketed and would consequently downgrade future prospects for Sartorius over the medium term.

Environmental Risks

Sartorius has established an environmental management system that encompasses, and is integrated into, all divisions and covers a whole series of environmentally relevant regulations to minimize environmental risks. This management system has been certified for compliance with ISO 14001 at a number of the company's relatively large manufacturing sites. The respective company organizational units ensure at the particular sites that the laws and regulations relating to environmental protection are observed and that further technical possibilities for limiting environmental risks are identified on an ongoing basis.

The increasing importance of sustainability considerations in many industries represents an opportunity. That is why this aspect is a key element in our supplier selection process for assessing the suitability of a particular company as a business partner. For further information on this group of topics, please refer to our Non-Financial Group Declaration.

IT Risks and Opportunities

Besides the risks already described, the Sartorius Group is exposed to potential risks in the area of IT as a result of its pronounced dependence on these systems, since their error-free operation is essential for the smooth functioning of the company's business processes. We reduce IT security risks by continuously enhancing and implementing IT security guidelines and policies. These rules and measures are based on the requirements of ISO 27001 and the standards of the German Federal Office for Information Security (BSI Standards). Furthermore, our company's existing IT applications and IT systems are checked for potential risks in regular external and internal IT audits, and appropriate measures are taken to minimize any risks identified. Continuous alignment of our IT strategy and business strategy, tracking of new technical developments and the use of advanced hardware and

software minimize the risk inherent in the operation of our IT system environment. A new ERP system commissioned by Sartorius at its Göttingen Group headquarters in 2012 has been successively rolled out to the Group sites around the world. Up to the end of 2017, approximately 75% of the Group's global sales transactions have been processed using the new system. In conducting this IT project, we have continued to focus on controlling the risks involved, such as by maintaining a precautionary backup system. The implementation of the new system brings with it a whole series of opportunities, especially in relation to efficiency gains and the standardization and harmonization of business processes worldwide.

Process Risks

Process risks for Sartorius can arise from pending or forthcoming legal disputes or from administrative proceedings. All judicial or extrajudicial disputes are attended to by the company's own attorneys and legal experts, who engage external lawyers as needed.

At present, there are no pending or discernible legal disputes or proceedings that lack any cost coverage allowances in the statement of financial position or that could have a substantial negative impact on Group.

Insurance

We have taken out insurance policies to cover a wide range of risks where possible and economically advisable. These insurance policies include coverage against liability, property damage, business interruption, transport, material and pecuniary damages and other risks, and provide comprehensive coverage for legal costs. An independent department working in conjunction with an external insurance broker regularly reviews the nature and extent of our insurance protection and makes any adjustments necessary.

When choosing our insurers, we particularly consider the credit rating of these entities as potential contractual partners, as well as aim to achieve a high degree of diversity in order to mitigate the related risks.

Assessment of the Overall Risk Situation and Risk Outlook

Where feasible, we adopted countermeasures and/or arranged for balance sheet measures during the reporting year to cover all discernible risks within the Sartorius Group, and those of a defined probability of occurrence, that had the potential to damage our net worth, financial situation and profitability.

For the purposes of this report, we have assessed the probability of occurrence of the risks as shown below and, in the adjacent columns, classify their particular significance for the entire Group.

Risk Category	Probability of Occurrence	Significance
General and macroeconomic risks		
Business cycle risks	Medium	Significant
General risks	Low	Significant
Supply chain risks		
Procurement risks	Low	Of limited significance
Produktion risks	Low	Significant
Sales and distribution risks	Medium	Significant
Quality risks	Low	Significant
R&D risks	Low	Significant
Customer risks	Low	Of limited significance
Competitive risks	Low	Of limited significance
Acquisition risks	Low	Significant
Personnel risks	Low	Of limited significance
Financial risks		
Risks associated with Group accounting	Low	Of limited significance
Exchange rate risks	Medium	Significant
Interest rate risks	Medium	Significant
Liquidity risks	Low	Significant
Regulatory risks	Low	Of limited significance
Environmental risks	Low	Of limited significance
IT risks	Low	Significant
Process risks	Low	Of limited significance

After thorough analysis of the entire risk situation and according to our current review, there are no discernible risks at present that could jeopardize the continued existence of the Group.

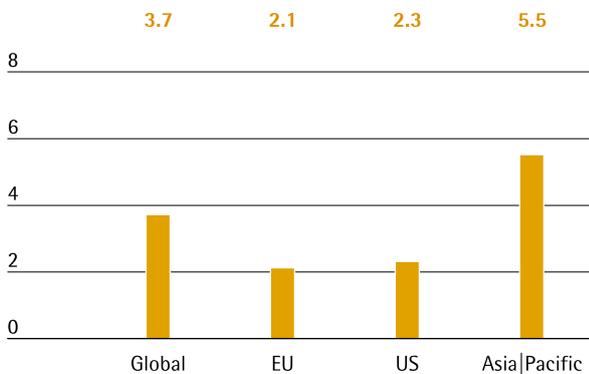
Similarly, based on our current review, there are no discernible risks that could jeopardize the future existence of the Group.

Forecast Report

Future Macroeconomic Environment

According to the October 2017 outlook by the International Monetary Fund (IMF), the global economy is expected to continue expanding at a high level in 2018. Economic growth is projected to reach 3.7%, up from 3.6% a year earlier, with the upswing supported by pickups in investment, global trade and industrial production.

Forecasted GDP Growth Rates for 2018
in %



Source: International Monetary Fund

The growth forecast for the European Union is moderate this year. Estimates call for growth of 2.1%, down from 2.3% in 2017. Medium-term growth could be negatively impacted by the overall weak productivity trend as well as excess levels of both public and private debt in some countries. The German economy in particular is expected to see a slight decline in 2018 compared with the historically high growth rates recently experienced. The IMF currently predicts growth of 1.8%, a decrease of around 0.2 percentage points on the prior year. Economists are projecting a similar decline for the United Kingdom, with growth forecast to drop to 1.5%. The depreciation of the British pound continues to negatively impact consumer spending.

For the U.S. economy, the IMF forecasts economic expansion of 2.3% in 2018, compared with 2.2% in the previous year. That projection does not account for possible effects resulting from planned corporate tax cuts by the U.S. administration. Economic researchers view the high confidence levels evidenced by both consumers and companies as fundamentally positive.

The Asia-Pacific economic area is likely to grow by around 5.5% in 2018, or approximately at the same level as the prior year. Growth will be fueled by a

revival of China's import demand. The world's largest economy is expected to expand by roughly 6.5% on the back of fiscal support programs (previous year: 6.8%). The Japanese economy is forecast to see a sharp decline in growth of around 0.8 percentage points to 0.7%, due to phase-outs of economic stimulus packages and decreasing consumer spending.

Future Exchange and Interest Rate Trends

Experts predict that base interest rates will remain at a very low level in 2018, as they did last year.

Market forecasts of the euro-U.S. dollar exchange rate for the course of 2018 range between 1.08 euro | U.S. dollar and 1.30 euro | U.S. dollar.

Sources: International Monetary Fund, World Economic Outlook October 2017; Reuters Forex Poll, December 2017

Outlook for the Sectors

Pharmaceutical Industry Continues to Grow

Strong, long-term trends are driving growth in the global pharmaceutical industry. Major growth factors include the growing and aging global population, combined with increasing access to healthcare, especially in the emerging and developing countries. However, the expiration of patents and measures aimed at limiting national healthcare expenditures will have a dampening impact. Market observers estimate growth of the world's pharmaceutical market to come in at between 4% and 7% per year during the period up to 2021.

The U.S. pharmaceutical market – the world's largest – is expected to grow at a rate of 6% to 9% on average during the period between 2017 and 2021. Expansion will be driven principally by new, innovative medications, while moderating effects will result from expiring patents for lucrative medications and increased use of biosimilars.

Growth in the European pharmaceutical market is likely to remain moderate over the next few years as continuing austerity measures affect its national healthcare systems. Thus, average growth of between 1% and 4% is projected for the region until 2021.

It is to be expected that the “pharmerging markets” and especially China will continue to expand at an above-average rate. Market observers estimate growth at around 6% to 9% annually from 2017 to 2021 owing to demographic trends, rising investment in government-led healthcare systems and increased private spending.

Biotech Sector Enjoys Above-Average Growth

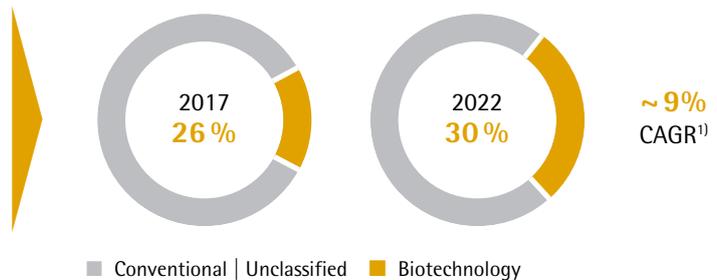
Market observers forecast that the biopharma segment of the pharmaceutical market, which has been enjoying particularly strong growth for years, will continue to outperform the market. Forecasts call for the global biotechnology market to grow by an average of around 9% annually during the period from 2017 to 2022, which would increase the market volume from the current €214 billion to €326 billion. The share of biological medications and vaccines in the total revenue generated by the pharmaceutical market is expected to rise from roughly 26% at present to around 30% in 2022.

This dynamic trend will be driven largely by the increasing market penetration of already approved biopharmaceuticals and an expanded range of indications. Thanks to technical and medical advances, the field of applications for substances manufactured using biotech production methods is growing steadily. For example, biopharmaceuticals are increasingly being used in yet-to-be fully explored therapeutic areas and in the treatment of rare diseases that have thus far been untreatable. This relatively young market segment is driving sector growth with its high innovate power, as reflected in the strong research and development pipelines. Overall, around 40% of the medications in R&D pipelines are based on biological manufacturing processes. These include more than 800 biosimilars and 500 “biobetters,” i.e. generic versions of biologic agents with better efficacy or fewer side effects than the original compounds.

The high level of innovation and R&D productivity in the biotechnology sector are mirrored by the rising number of new product approvals. In the USA for example, the number of newly approved biological medications during the last five years was approximately 120% higher than the number of approvals from the period 2008 to 2012.

Strong Growth in the Biopharma Market

- Growing & aging population
- Increasing access to healthcare
- Strong R&D pipelines
- Emerging biosimilars market



¹⁾ Source: Evaluate Pharma®: World Preview 2017, Outlook to 2022, June 2017; CAGR 2017 to 2022.

Biosimilars which are biological copycat medications, will continue to increase in importance during the years to come. They are less expensive than the reference compound, while having similar properties in terms of their application, dosage and efficacy. Increased use of biosimilars thus brings the prospect of lower individual treatment costs and financial relief for national healthcare systems. This is a goal likely to be pursued by the industrialized countries, where aging populations are leading to rising demand for medications and notably higher expenditures. In many developing countries, biosimilars offer the initial access to biopharmaceutical substances. The development of national production capacities is receiving political support. Regulatory requirements are generally much less strict in developing countries than in the industrialized countries, thus the biosimilars market could experience substantial growth in the emerging economies over the next few years.

Eight of the strongest selling biopharmaceuticals are due to lose their patent protection in the United States and Europe by the year 2020. Various pharmaceutical manufacturers have therefore announced the market launch of those biosimilars that have already been approved. At present, around 30 companies are actively working on the development of biopharmaceutical copycat products for 16 different molecules. At the same time, the biosimilars market will also see a gradual decline in regulatory and patent-related barriers and challenges to marketing the products – factors that have thus far slowed development of the market. Against this backdrop, current estimates call for accelerated growth over the next few years. By 2020, sales of biosimilars could multiply to between U.S. \$25 billion and U.S. \$35 billion.

Furthermore, some spectacular achievements have been made in the field of cell therapy lately. The goal of cell therapy is to prevent, treat, heal or mitigate human disease or injury through the use of autologous, allogenic or xenogenic cells that have been manipulated ex vivo or otherwise altered. Regardless of the type of cell therapy, the production technologies used for the manufacturing of the respective therapeutic products are similar to those used for the manufacturing of biopharmaceuticals. The growing number of cell-based therapies currently in clinical development and the rising number of approved cell therapies already in use could lead to increasing demand for the corresponding production capacities in the future.

Stable Growth Expected in the World's Laboratory Market

According to Frost & Sullivan, global demand for laboratory products is likely to remain stable, with growth of 2.8% in 2018. The key U.S. market is expected to generate growth of 3.0%, in part due to the sustained strength of the biopharmaceutical industry and the business-friendly climate. In Europe, experts anticipate an increase of just 1.7% owing to the comparatively moderate economic growth and as a result of uncertainties in the wake of the Brexit referendum. However, the European Union's increased budget for academic research will likely spur mid-term growth. As before, market observers expect the highest growth rates in Asian countries such as China and India, in which the individual labor markets are likely to grow around 7.5% to 8.8% in 2018.

Sources: QuintilesIMS Institute: Outlook for Global Medicines through 2021, December 2016; Evaluate Pharma: World Preview 2017, Outlook to 2022, June 2017; BioPlan: 14th Annual Report and Survey of Biopharmaceutical Manufacturing Capacity and Production, April 2017; Deloitte: 2017 Global Life Science Outlook, October 2016; Frost & Sullivan: 2017 Mid-year Report: Forecast and Analysis of the Global Market for Laboratory Products, May 2017

Outlook for 2018

The outlook for 2018 considers the sector environment, economic trends and the opportunities and risks outlined in this Annual Report. The following forecast is provided under the assumption that we have accurately anticipated the relevant trends for Sartorius:

Sartorius Group

Sartorius expects to grow profitably again in 2018. Specifically, management projects that Group sales revenue for the full year will grow by about 9% to 12% in constant currencies. Regarding profitability, management forecasts that the company's underlying EBITDA margin will increase by around half a percentage point over the prior-year figure of 25.1%.

In view of continued strong organic growth, Sartorius had already started in 2016 to expand its production capacities earlier and to a greater extent than initially planned. Accordingly, the Group projects that it will invest overproportionately in its worldwide infrastructure in 2018 as well, at a ratio of capital expenditures to sales revenue of around 15%. The focal areas of its investments will be to extend the Yauco site for manufacturing single-use filters and bags, as well as to combine and expand Group headquarters in Göttingen, Germany.

With regard to our financial position, we forecast that our ratio of net debt to underlying EBITDA will be slightly below the previous year's level (Dec. 31, 2017: 2.5) at the end of 2018. This ratio does not take any potential acquisitions into account.

Division Forecasts

For the Bioprocess Solutions Division, management expects significant increases for the full year of 2018 as well. The division's sales revenue is projected to grow by about 8% to 11%. The division's underlying EBITDA margin is anticipated to rise by around 0.5 of a percentage point over the prior-year figure of 28.0%.

The Lab Products & Services Division in some parts of its portfolio depends on general economic cycles. For this division, management projects that, assuming an overall stable economic environment, sales will increase by about 12% to 15%. The division's underlying EBITDA margin is forecasted to rise by about one percentage point from 18.0% in the previous year.

All forecasted figures are given in constant currencies. Due to the latest currency developments, especially in the exchange rate between the U.S. dollar and the euro, our reported figures in actual currencies might differ from our constant currency guidance in which we will provide additional details during the full year of 2018.

Sustained Reduction in Tax Rate

Sartorius' corporate tax rate is expected to be reduced by about 2 percentage points to around 27% from 2018 onwards as a future effect of the U.S. tax reform approved in December 2017, which essentially lowers the U.S. corporate tax rate from 35% to 21%. Since the changes in American tax legislation are very extensive, and further explanations and instructions for application by the U.S. tax authorities are still pending, this estimate is to be considered tentative.

Report on Material Events

No material events occurred after the close of fiscal 2017.

Description of the Key Features of the Internal Control Management System

In relation to the Group Accounting Process (Section 289, Subsection 4, and Section 315, Subsection 4, of the German Commercial Code [HGB])

Definitions and Elements of the Internal Control Management System at the Sartorius Group

The internal control system of Sartorius AG and the Sartorius Group encompasses all of the principles, procedures and measures adopted to ensure the organizational implementation of management decisions. The main priority of the system as it relates to Sartorius AG's and the Group's accounting process is to make sure that accounting is effective, cost-efficient and formally correct and that it complies with the pertinent legal provisions.

The internal control system of the Sartorius AG and the Sartorius Group consists of a combination of process-integrated and non-process-integrated monitoring measures. The process-integrated safeguarding measures are subdivided, in turn, into organizational measures and other control measures. The Supervisory Board, specifically in this case the Audit Committee of Sartorius AG, and the Legal Affairs & Compliance unit with its Group Auditing department are involved in the Sartorius Group's internal control system through their non-process-integrated audit activities. The Audit Committee regularly reviews quarterly reports besides the annual financial statements of the parent corporation and of the consolidated annual financial statements.

Moreover, to ensure systematic, early identification of risks across the entire Group, a "monitoring system for early group-wide detection of risks with the potential to jeopardize the company's continued existence" as defined in Section 91, Subsection 2, of the German Stock Corporation Law (AktG) is in place at the Sartorius Group. The efficacy of the early risk detection system, which the Sartorius Group adapts promptly in response to any relevant changes in circumstances, is assessed by the independent auditors of Sartorius AG in accordance with Section 317, Subsection 4, of the German Commercial Code (HGB). An integral component of this system is also operational risk management, which involves activities such as the transfer of risk to insurance companies through coverage for damage and liability risks and the arrangement of suitable hedges to limit currency and interest rate risks.

Organizational Measures

Accounting processes are strictly organized according to the principle of segregation of functions, taking into account the "four-eyes" principle – i.e. review by two individuals. Duties and responsibilities are clearly assigned to different specialized units, companies and regional units. The separation of administrative, executive, settlement and approval functions reduces the possibility of fraud. It also continues to play a significant role in ensuring that any possible errors are discovered early and any potential misconduct is prevented.

The IT applications used in the company's accounting processes have access restrictions, which allow only authorized persons to have controlled access to the accounting system and data. Each access right is assigned specifically according to the tasks to be performed and will be subject to annual review in the future. Furthermore, the four-eyes principle is also applied in IT process designing and assignment of access rights.

In addition, defined written local and global operating procedures exist, particularly the Group accounting guidelines, which are regularly updated and communicated throughout the Group. The scope of regulation at Group level also extends to the central definition of measurement rules and parameters, among other factors. Additional data for the presentation of external information in the notes to the financial statements and in the Group management report is also prepared and aggregated at Group level.

Continuous coordination of internal accounting during the year for planning and control with external accounting significantly contributes to the quality of Group financial reporting. Reporting itself is done through a standardized management reporting system implemented throughout the Group. This system visualizes all consolidation processes. Internal controls, on the one hand, and the Group auditors of Sartorius AG, on the other hand, ensure that Group financial reports are accurately generated from the consolidated Group companies' financial statements.

The employees involved in the accounting process meet qualitative standards and receive regular training. The Group Financial Reporting department assists the local units in resolving complex accounting issues, such as measuring fair value, to ensure consistent and

accurate reporting in the consolidated financial statements. Complex evaluations, such as actuarial reports and company valuations or purchase price allocations, are assigned to specialized service providers who involve the respectively qualified in-house staff.

Control Measures

Comprehensive control activities are performed by managers and staff to ensure effective and reliable accounting. As a result, this ensures compliance with legal requirements and internal guidelines, as well as properly conducted business transactions. Examples of such control activities include the analysis of situations and developments with reference to specific key indicators. Moreover, individual reporting units monthly comment on and explain special characteristics or variances using Group-wide standardized analytical tools as the basis. Further specific control activities performed to ensure effective and reliable Group accounting encompass the analysis and, where applicable, correction of the individual financial statements submitted by the Sartorius Group companies. A large number of control mechanisms already incorporated into the consolidated reporting system enable erroneous information to be identified and corrected at Group level. Impairment tests are conducted centrally for the specific cash-generating units, known as CGUs, from the Group's perspective to ensure that consistent, standardized evaluation criteria are applied.

The Legal Affairs & Compliance unit annually draws up a risk-based audit plan and reviews in spot checks whether basic legal requirements and internal group guidelines are complied with for the entire control and risk management system of the Group. This monitoring function covers, in particular, auditing the functional efficiency and effectiveness of defined control measures. The results of these audits are reported directly to the Executive Board and enable any deficiencies determined to be corrected and the company's internal control system (ICS) to be further enhanced.

A manual on the company's internal control system that focuses on the business processes of our company further contributes toward strengthening ICS. As a result, this manual combines all ICS-relevant requirements that we consider of material importance into one standardized document and will be supplemented by further appropriate rules as necessary.

Qualifying Statements

The internal control and risk management system enables the complete recording, processing and evaluation of company-related matters, on the basis of the organizational, control and monitoring structures defined in the Sartorius Group, as well as their accurate presentation in Group accounting.

In particular, decisions based personal judgment, erroneous controls, criminal acts and other circumstances that can impair the efficacy and reliability of the internal control and risk management system in place cannot be ruled out, however, so even the application throughout the Group of the systems adopted cannot provide an absolute assurance as to the accurate, complete and timely recording of matters in Group accounting.

The statements made relate solely to the subsidiaries included in the consolidated financial statements of Sartorius AG, whose financial and business policies Sartorius AG can directly or indirectly determine in order to obtain benefits from their activities.

Explanatory Report of the Executive Board

On the Disclosures Pursuant to Section 289a, Subsection 1, and Section 315a, Subsection 1, of the German Commercial Code (HGB)

Composition of the Issued Capital | Limitations to Voting Rights

Sartorius AG's capital stock totals €74,800,000. It comprises 74,800,000 no par value individual bearer shares, 37,440,000 of which are ordinary shares and 37,440,000 of which are non-voting preference shares. Each share certificate represents a calculated proportion of €1 of the issued capital.

The rights and obligations associated with these shares are governed by the provisions of the German Stock Corporation Law (Aktengesetz, abbreviated "AktG"). According to the company's Articles of Association, preference shares are entitled to a dividend payment that is one euro cent higher per share than that for ordinary shares. However, this entitlement to receive dividends shall be at least two euro cents per preference share. Apart from the cases provided for in Sections 140 and 141 of AktG, preference shares are non-voting. Beyond this, preference shares grant all other rights to which every shareholder is entitled.

The company holds 3,227,776 ordinary shares and 3,263,932 preference shares; these do not entitle the company to any membership rights.

Direct or Indirect Equity Ownership Exceeding 10% of Voting Rights

The community of heirs consisting of Mrs. U. Baro, resident of Munich, Germany; Mrs. C. Franken, resident of Bovenden, Germany; Mr. A. Franken, resident of Riemerling, Germany; Mr. K.-C. Franken, resident of Göttingen, Germany; and Mrs. K. Sartorius-Herbst, resident of Northeim, Germany, holds a voting percentage of approximately 50.1% in Sartorius AG (18,754,160 votes; source: according to the list of attendees at the Annual Shareholders' Meeting on April 6, 2017). The decedent Horst Sartorius ordered that his will be administered by an executor. Dr. Lothar Kappich, resident of Hamburg, Germany, was appointed executor to succeed Prof. Dr. Dres. h.c. Arnold Picot who unexpectedly passed away on July 9, 2017. The executor exercises the specified voting rights at his own discretion as defined by Section 22, Subsection 1, sentence 1, item no. 6, of the German Securities Trading Act (Wertpapierhandelsgesetz, abbreviated "WpHG").

According to a mandatory announcement dated April 1, 2011, Bio-Rad Laboratories Inc., 1000 Alfred Nobel Drive, Hercules, California 94547, USA, to which the voting rights of Bio-Rad Laboratories GmbH, Heidemannstr. 164,

80939 Munich, Germany, are ascribed according to Section 22, Subsection 1, sentence 1, item no. 1, of WpHG, holds 30.01% of the voting rights in Sartorius AG (11,237,196 votes, taking into account the share capital increase from retained earnings that went into effect as of June 1, 2016).

Appointment and Dismissal of Executive Board Members | Amendment to the Articles of Association

Executive Board members of Sartorius AG are nominated and/or appointed as well as removed from office in accordance with Sections 84 et seq. of the German Stock Corporation Law (AktG) and Sections 31 and 33 of the German Codetermination Law (Mitbestimmungsgesetz, abbreviated "MitBestG"). Amendments to Sartorius AG's Articles of Association are regulated by Sections 133 and 179 of the German Stock Corporation Law (AktG).

Powers of the Executive Board to Issue Shares

Subject to approval by the Supervisory Board, the Executive Board is authorized to sell treasury shares held by the corporation, including selling them through channels other than the stock exchange or by tendering an offer to all shareholders in proportion to their participation in the company, provided that these shares are transferred to third parties as contribution in kind, particularly in the (indirect) acquisition of companies, in return. Under these circumstances, the preemptive rights of the shareholders are excluded.

Material Agreements with Clauses Regulating the Event of a Change of Control

Sartorius AG has material loan agreements containing customary market clauses regulating the possible event of a change of control and giving the participating lenders the option of extraordinary termination of a particular agreement or to request complete repayment of the respective loan:

These are a syndicated loan agreement concluded in 2014 for a current volume of €400 million and a term until December 2021; two note loans ("Schuldscheindarlehen"), for the amount of €200 million placed in 2016 and 2017, respectively; and a €350 million note loan for a maturity term of five to ten years; two promotional loans, each for €90 million, concluded in 2015 and 2016 for maturity terms of 8 years; and a note loan originally for €100 million and placed in 2012 for a maturity term of five to ten years. The repayment sum outstanding for these material loan agreements stands at approximately €760 million as of December 31, 2017.

Report and Declaration on Corporate Governance

Corporate Governance Report

Corporate government aligned with the interests of stakeholders, lawful and responsible conduct, and constructive cooperation between the managerial bodies and within the company in a spirit of mutual trust constitute the essential cornerstones of Sartorius' corporate culture.

The Executive Board and the Supervisory Board report in the following declaration on the key aspects of corporate management and governance pursuant to § 289f of the German Commercial Code ("HGB") and to Article 3.10 of the German Corporate Governance Code.

Declaration of Compliance with Corporate Governance

Declaration of the Executive Board and of the Supervisory Board of Sartorius AG Concerning the Recommendations of the Government Commission on the German Corporate Governance Code Pursuant to § 161 of the German Stock Corporation Law ("Aktiengesetz"):

The Executive Board and the Supervisory Board declare that full compliance will be achieved with the recommendations made by the Government Commission on the German Corporate Governance Code and published by the German Federal Ministry of Justice in the official section of the German Federal Gazette ("Bundesanzeiger"), as amended on February 7, 2017.

Since last year's Declaration of Compliance was issued, Sartorius AG has complied with the recommendations in the valid version of the Government Commission on the German Corporate Governance Code to the full extent.

Göttingen, December 7, 2017

For the Supervisory Board For the Executive Board

Dr. Lothar Kappich

Dr. Joachim Kreuzburg

Further Remarks Concerning Corporate Governance

Sartorius AG is a joint stock corporation founded under German law and headquartered in Göttingen, Germany. With the Annual General Shareholders' Meeting, Supervisory Board and Executive Board, it has three corporate managerial bodies whose tasks and powers are essentially derived from the German Stock Corporation Law ("Aktiengesetz") and the company's Articles of Association.

As owners of the company, the shareholders exercise their rights at its Annual General Shareholders' Meeting, where they decide, in particular, on the appropriation of profits, measures concerning share capital, amendments to the Articles of Association, discharge of the Supervisory Board and the Executive Board and on the appointment of statutory auditors, as well as elect shareholder representatives to the Supervisory Board. The Annual General Shareholders' meeting is held at least once a year within the first eight months of the respective fiscal year. Any shareholder who registers in due time may attend the Annual General Shareholders' Meeting in person or can be represented by a proxy according to the various options provided by law.

In managing the company, the Supervisory Board and the Executive Board perform their tasks in a two-tier board structure, each with separate duties and powers.

The Supervisory Board appoints members to the Executive Board, determines their remuneration and monitors and advises the Executive Board in its management of the company. The Supervisory Board is not authorized to take any operational management measures for the business.

The Executive Board is responsible for independently managing the company. In particular, it defines corporate strategy, coordinates and agrees on this approach with the Supervisory Board and implements such corporate strategy. In line with established reporting obligations, the Executive Board regularly informs the Supervisory Board promptly and comprehensively, and requests the latter's approval for certain key business transactions.

Composition and Operating Mode of the Supervisory Board and Its Committees

The Supervisory Board has an equal number of shareholder representatives and employee representatives:

six shareholder representatives elected by the Annual General Shareholders' Meeting and six employee representatives elected according to the German Code-termination Law ("Mitbestimmungsgesetz"). Details on the members of the Supervisory Board and its committees are provided on pages 180 to 181.

The Supervisory Board Chairman coordinates the work within this board, as well as convokes and conducts Supervisory Board meetings. Furthermore, he is the first individual for the Executive Board to contact and externally represents the matters of the Supervisory Board.

The Supervisory Board holds at least two meetings every six months; as a rule, four or more conferences take place, as required. This board has established four committees: the Executive Task Committee, the Audit Committee, the Conciliation Committee and the Nomination Committee. The Executive Task Committee, Audit Committee and Conciliation Committee each have four members and have an equal number of shareholder representatives and employee representatives. The Executive Task Committee and Audit Committee hold regular meetings; the Conciliation Committee and the Nomination Committee meet only as necessary.

The Executive Task Committee carries out preparatory work for resolutions and issues to be addressed in the meetings of the Supervisory Board. It also oversees the preparations for appointments, including the remuneration and employment contract conditions of members of the Executive Board. The Audit Committee supports the Supervisory Board in performing its supervisory function. The chairperson of the Audit Committee is an independent member of the Supervisory Board and has detailed knowledge and extensive experience in the application of accounting standards and internal control systems from his or her own professional practice. The Conciliation Committee meets if the majority required in connection with the appointment of members to the bodies authorized to represent the company for legal purposes is not reached. The Nomination Committee comprises representatives of the shareholders only. Its function is to propose suitable candidates to the Supervisory Board for the latter's election proposals submitted to the Annual Shareholders' Meeting. More information on the individual meetings held in the reporting year by the Supervisory Board and its committees is given in the Report of the Supervisory Board on pages 12 to 15.

The Supervisory Board regularly reviews the efficiency of its work based on a questionnaire sent in advance and by a discussion of the results of this questionnaire in a plenary session.

Appointment Objectives for the Supervisory Board; Diversity Policy and Competence Profile

Members of the Supervisory Board of Sartorius AG are to be appointed such that they, on the whole, have the knowledge, skills and experience that are necessary to perform the board's duties properly.

For this purpose and based on the recommendations of the German Corporate Governance Code, the Supervisory Board decided on the following appointment objectives:

- Independent members shall account for no less than 25% of the Supervisory Board.
- The upper age limit of 70 at the time of election should be taken into consideration, but may be waived in exceptional cases, provided there are no reservations about the suitability of the persons proposed and their election is expedient to the interests of the company in spite of the age limit being exceeded.
- As a matter of principle, no member may serve on the Supervisory Board for longer than five elected terms. This limit can be exceeded in individual cases if the Supervisory Board does not have any reservations about the suitability of an individual for serving longer and election of this person is apparently in the interest of the company.
- Care must be taken to ensure that each Supervisory Board member has sufficient time available to perform his or her mandate.

In addition, the Supervisory Board also defined a competence profile for itself. This also covers aspects such as diversity, for example, with regard to professional background and international experience. In view of achieving an appropriate gender balance, the legal quotas of at least 30% women and at least 30% men apply to the Sartorius Supervisory Board. The shareholder representatives and the employee representatives decided to fulfill these legal targets separately.

- In its election proposals for membership, the Supervisory Board is required to consider whether candidates have international experience or an international background within the scope of its current members.
- The Supervisory Board must have members with expertise in one or several of the international markets relevant for Sartorius.

- Members of the Supervisory Board need to bring in knowledge of technologies and products relevant to the Group, as well as of digitalization, and experience in research and development, particularly in the biopharmaceutical industry.
- The Supervisory Board must have members who have expertise in setting up and developing innovative business models, as well as knowledge of corporate strategies.
- Members of the Supervisory Board need to have in-depth knowledge of financial business processes and competences in financial controlling and risk management; at least one independent member of the Supervisory Board must have expert knowledge of accounting or auditing (§ 100, Subsection 5, of AktG).
- Members of the Supervisory Board must have expertise in law, corporate governance and compliance.
- The Supervisory Board needs to have in-depth knowledge and a thorough background in issues concerning human resources.

According to the Supervisory's self-assessment, the members on its board meet the diversity and competency requirements. The appointment objectives previously mentioned regarding independence and the age and membership limits are also met:

Concerning the criterion of independence, this is met by all members of the Supervisory Board with one exception in the past and the present: the executor of Horst Sartorius' estate. This individual was Dr. Arnold Picot, Ph.D., until July 9, 2017, who was also the Chairman of the Supervisory Board, and is now Dr. Lothar Kappich, likewise its current Chairman. In the opinion of the Supervisory Board, an appointment as an employee representative to this board or an existing employment contract with the company alone does not preclude the independence of this member because an employee is protected by the laws in force in performing his or her duties on the Supervisory Board.

In one instance, the company diverged from the established statutory retirement age and the statutory limit of five elected terms when Dr. Arnold Picot was reelected. There were no doubts about his suitability, and his reelection was also in interest of the company in the context of his capacity as executor of Horst Sartorius' estate. Currently, there are no Supervisory Board members over the age of 70 or who have served for more than five elected terms.

In view of achieving an equal balance of gender, the Supervisory Board meets the quota of 30% set for the underrepresented gender. Since the election of members to the Supervisory Board in April 2017, seven men total (around 58%) are on the board, of whom four are shareholder representatives and three employee representatives. In addition, five women (around 42%) are members of this board, among them two representatives of the share owners and three representatives of the employees. As a result, the gender quota requirements are met on both sides of Supervisory Board representation and on the full Supervisory Board itself.

The proposal of the Supervisory Board submitted to the Annual General Shareholders' Meeting for election of shareholder representatives to this board was based on the target quotas stated above.

To facilitate comparison of the appointment objectives, brief resumés of the Supervisory Board members are available on the Sartorius website.

Composition and Operating Mode of the Executive Board

The Executive Board of Sartorius AG manages the company under its own responsibility, with the goal of increasing the company's sustainable value. It develops the company's strategy, coordinates it with the Supervisory Board and ensures implementation of this strategy. Beyond that, the rules of procedure for the Executive Board define the legal transactions requiring approval by the Supervisory Board in order for such transactions to be effected. The Executive Board is responsible for compliance with all provisions of the law and the company's internal policies, as well as for appropriate risk management.

Decision-making by the Executive Board is done at its regular meetings, which are convoked and conducted by the Chairman. As a rule, the members of the Group Executive Committee attend these meetings, and, if required, further specialists and managers are invited to provide advice.

The Executive Board members are jointly responsible as a collegiate body for matters of special significance. As for the board's remaining responsibilities, each member independently manages the area assigned to him or her according to the plan for allocation of areas of responsibilities, and is required to notify the Chairman of all material transactions and events.

Composition of the Executive Board, Diversity and Competency Requirements

In the opinion of the Supervisory Board, the basic qualification criteria for appointments to positions on the Executive Board are professional qualifications for heading each particular area of responsibility, a proven track record in the individual's career path and convincing managerial skills. In addition, the Supervisory Board also considers the aspect of diversity in its appointment decisions. Therefore, the Supervisory Board strives to appoint people with complementary profiles, professional and personal life experiences and in different age brackets to the Executive Board. Moreover, the latter board is required to have broad international experience.

Pursuant to the German Act on Equal Participation of Women and Men in Executive Positions in the Private and the Public Sectors (abbreviated as "German Act on Equal Participation of Women and Men" below), the Supervisory Board defined a target quota for the Executive Board that is defined in the following section.

Disclosures Concerning the German Act on Equal Participation of Women and Men

The Sartorius Executive Board is a committee that presently consists of three members and is therefore relatively small so establishment of a rigid gender quota can be problematic. The current members of this board of Sartorius AG are male. Given that the Executive Board contracts are still in force, the Supervisory Board had defined a quota of 0% up to June 30, 2022. For the current target achievement deadline by June 30, 2022, the Supervisory Board decided to appoint at least one woman to the Executive Board.

Also regarding the appointment of women to the Executive Board of Sartorius AG, the Supervisory Board supports the activities of the Executive Board to further increase the percentage of female executives at the first two management levels in the company. The Executive Task Committee responsible and the full Supervisory Board regularly receive reports on the development of the proportions of women in senior-level management positions.

First and Second Management Levels Below the Executive Board

Disclosures Concerning the German Act on Equal Participation of Women and Men

Over the past years, the percentage of women at the first two management levels below the Executive

Board has considerably increased on the whole and is already at a comparably high level. To achieve the intended communications effect, the Executive Board had deliberately set ambitious targets for the first target achievement deadline by June 30, 2017, despite this short time horizon and the risk that these targets might not be completely met. Thus, this quota for the first management level was set at 25% and for the second level, at around 30%.

Although the percentage of women at these levels increased, both targets were not quite reached. At the end of June 2017, the proportion of women at the first management level was around 19%; that at the second level, around 28%. The reasons these targets were not met were the low fluctuation in the management team and, therefore, a lower number of positions to be filled, on the one hand, and consolidation of the companies acquired slightly diluted the progress achieved, on the other hand.

For the next deadline by June 30, 2022, the Executive Board resolved in the reporting year to increase the proportion of women at both levels of management below this board to around 30%.

Further Corporate Governance Practices

Risk Management

Conscientious management of commercial risks is a key principle of good corporate governance. Sartorius AG and the Group have at their disposal enterprise-wide and company-specific reporting and control systems designed to facilitate the recording, assessment and management of commercial risks. These systems are developed and adapted continuously as conditions evolve. The Executive Board informs the Supervisory Board regularly of existing risks and their development. The Audit Committee is concerned, in particular, with monitoring of the following: the accounting process including reporting; the efficacy of the internal control system; risk management and the internal auditing system; compliance; and the independent statutory audit. Details on risk management are presented in the Opportunity and Risk Report.

Transparency

Sartorius AG places great importance on disclosing consistent and complete information promptly. Information about the economic position of the Group and new developments is consequently released regularly, without delay, as it becomes known in order to inform

participants in the capital market and interested members of the public at large. The annual report, first-half financial report and quarterly reports are published within the timeframes specified for this purpose. Current developments and material events are publicized as press releases and, where appropriate, ad hoc announcements. This information is usually made available in German and English simultaneously and published via suitable media and on the internet.

The chief recurring events and publications, such as the Annual General Shareholders' Meeting, the annual report and the interim reports, are listed on a financial calendar that may be viewed at any time on the Group website.

Share Trading Activities of Supervisory and Executive Board Members

We did not receive any reports, under the applicable mandatory disclosure requirements, of any purchases or sales of shares in Sartorius AG or related financial instruments made by the members of the Executive Board and the Supervisory Board or other persons with management responsibilities or their related parties.

Supervisory Board Chairman Dr. Lothar Kappich in his capacity as executor of the estate of Horst Sartorius holds approximately 50.1% of the ordinary shares issued by the company, but otherwise no member of the Executive Board or Supervisory Board has any holding of shares or financial instruments subject to the mandatory reporting requirements that directly or indirectly exceeds 1% of the shares issued by the company.

Considering the share capital increase, by use of retained earnings, that became effective June 1, 2016, the CEO and Chairman of the Executive Board Dr. Joachim Kreuzburg holds 100,000 ordinary shares and 100,000 preference shares in the company. These were transferred to him based on a corresponding agreement arising from his employment contract of December 18, 2015, and are subject to a minimum holding period of four years. For further information on this transfer, please see the Remuneration Report on pp. 76 et seq.

Accounting and Independent Statutory Audit

The consolidated financial statements and the Group Management Report, as well as the consolidated interim financial statements and reports, are prepared in accordance with the International Financial Reporting Standards (IFRS) as they are to be applied within the

EU, and according to the commercial law regulations to be applied under Section 315a, Subsection 1, of the German Commercial Code, HGB. The annual financial statements of Sartorius AG are prepared in accordance with German commercial law, HGB. The consolidated financial statements and the annual financial statements are prepared by the Executive Board, audited by the independent auditors elected by the Annual General Shareholders' Meeting and approved by the Supervisory Board.

It has been agreed with the independent auditors that they will notify the Supervisory Board directly of any potential disqualification or bias issues and any material findings and incidents identified during the audit. This also encompasses the corporate governance reporting duties pursuant to § 161 of the German Stock Corporation Law (Aktiengesetz).

Basic Principles of Our Compliance Management System | Code of Conduct

With its compliance management system that is valid worldwide, Sartorius ensures that the members of its individual boards, executives and employees comply with all legal regulations and codes, and perform their activities in accordance with the company's internal rules and guidelines. Targeted training and awareness-raising prevent any misconduct, as well as economic damage and loss of image.

Sartorius makes every effort to ensure optimal risk management by using a combination of approaches: a preventive compliance approach designed to proactively stop any potential breaches before they occur and a repressive compliance approach intended to continuously monitor compliance with the company's rules. These processes are closely intermeshed, creating a standardized compliance management system that aims to offer the best possible protection against potential violations of rules and regulations. Sartorius has developed a Code of Conduct as a preventive component of its compliance management system and has committed to an Anti-Corruption Code. An internal system is available for reporting any suspicious circumstances involving potential compliance violations.

Further information is given in the Sustainability Report on pp. 68 et seq. and on the company's internet website at www.sartorius.com.

The Supervisory Board | The Executive Board

Remuneration Report

1. Main Features of the Remuneration Plan for the Executive Board

General Information

The full Supervisory Board is responsible for establishing the remuneration paid to members of the Executive Board of Sartorius AG. The total value of the remuneration of an Executive Board member reflects the scope of the responsibilities of the Executive Board member concerned, the Executive Board member's personal performance, the company's economic situation and sustainable progress. In addition, the extent to which this amount of remuneration is typical is considered, taking into account peer companies and the remuneration structure in place in other areas of the company and in similar companies.

Remuneration is comprised of both fixed non-performance-based components and of variable performance-based components, and is reviewed annually to ensure that it remains appropriate. The variable performance-based remuneration components consist of those to be paid annually and of multi-year components intended to have a long-term incentive. Fixed non-performance-based remuneration is paid in the year in which it is granted. For 100% target achievement, the variable annual and long-term performance-based components generally represent half of total remuneration, which excludes pension commitments under a defined benefit plan as well as fringe benefits.

Variable Performance-Based Remuneration

The portion of the variable performance-based remuneration that is paid annually is based on the following weighted components: sales revenue | order intake, underlying EBITDA and the ratio of net debt to EBITDA. Minimum target achievement is required for these components. The amount to be paid out depends on the degree to which the target is achieved, which the Supervisory Board defines by setting each individual subordinate target. A cap is provided for each variable component to be paid out. Variable performance-based remuneration is calculated upon approval of the company's annual financial statements for the respective fiscal year and not settled and paid out until the following fiscal year.

Multi-year Components as Long-term Incentives

Weighted components determined by multi-year assessment depend on the development of consolidated net profit in a multi-year period, on the one hand, and on the development of the company's share prices, on the other hand. Multi-year components providing a long-term incentive are based on a three-year average of consolidated net profit and on a four-year average of share prices, respectively. These components are paid out after two fiscal years for net profit and at the earliest after three fiscal years for share prices.

a) Consolidated Net Profit

For this subordinate target, the basis for assessment is the consolidated net profit after non-controlling interest excluding amortization (impairment of the value of intangible assets due to business combinations pursuant to IFRS 3). Target achievement for assessing annual variable remuneration in the particular year under review is based on the average taken over a period of three fiscal years, beginning with the particular year under review. To smooth the amounts to be paid out, a partial payment amounting to 50% of the target achievement for the first fiscal year of each respective average period will be effected. Any overpayments as a result of these partial payments will be offset against other remuneration components once the total target achievement has been determined after the third fiscal year of an average period. A cap for this component is provided as well.

b) Phantom Stock Plan

Through the issue of shadow shares, called phantom stock, Executive Board members are treated as if they were owners of a certain number of shares in Sartorius AG, without, however, being entitled to receive dividends. The development of the value of this phantom stock is linked with the development of the Sartorius share; both increases and decreases in the share price are taken into account. Later, the value of this phantom stock is assessed based on the share price at the time, and its equivalent is paid out, provided that the associated conditions are met. Phantom stock cannot be traded and does not entail any share subscription rights.

According to the Sartorius phantom stock plan, each Executive Board member is credited at the beginning of every year with phantom stock units valued at an

agreed monetary sum. The value of this phantom stock can be paid out only as an entire annual tranche. Payment can be requested, at the earliest, after a period of four years and no later than after eight years.

An Executive Board member is entitled to receive payment for phantom stock units only if the share price at the time of such payment request has appreciated at least 7.5% per year relative to the time the phantom stock was assigned or if the share price outperformed the TecDAX as a comparative index. The phantom stock plan rules out subsequent changes to the parameters used for comparative stock valuation. The amount to be paid is capped at a maximum of 2.5 times the share price at the time the phantom stock was assigned, based in each case on the actual annual tranche concerned.

Assignment of this phantom stock and later payment of its monetary equivalent depend on the mean value calculated from the average prices of both classes of Sartorius AG share, up to the year 2015, and as of the year 2016, on the mean value calculated from the average prices of the preference share, with said prices quoted in the closing auction of Xetra trading on the Frankfurt Stock Exchange over the last 20 days of trading of the previous year or over the last 20 days of trading prior to submission of a payment request. This serves to compensate for any short-term fluctuations in the share prices.

Payment for phantom stock is blocked for the four weeks preceding the scheduled publication date of quarterly and preliminary year-end results and for 20 days of trading on the stock exchange following the actual publication of quarterly and preliminary year-end results. These blackout periods are intended to prevent Executive Board members' profiting from their insider knowledge.

Pension Commitments

According to the company's remuneration policy, Executive Board members of Sartorius AG receive performance-related benefit commitments under a defined benefit plan when reappointed for the first time. In addition to including a basic pension, these commitments provide for the Executive Board member to make his own contribution from his variable earnings and for the company to match this contribution by a bonus amount. An Executive Board member may choose to receive such defined benefits in the form of a retirement pension for old age or as a one-time payment to cover the member's retirement pension for old age and invalidity as well as in the form of survi-

vor's benefits for the surviving spouse and children of the decedent.

Beyond such commitments, an Executive Board member is additionally entitled under a former company pension scheme to receive performance-based retirement benefits based on the salary of a German federal civil servant classified as grade 10 of salary class B for ministry officials according to the Federal Civil Service Remuneration Act (Bundesbesoldungsgesetz). Such benefits are paid in the form of a retirement pension or old age and invalidity as well as in the form of survivors' benefits for the surviving spouse and children of the decedent.

After a member has turned 65, this shall be considered the regular age limit at which this member shall automatically be entitled to receive all such benefits.

Other Remuneration Components

The remuneration system provides that the Supervisory Board of Sartorius AG at its discretion may grant an Executive Board member special compensation based on that member's exceptional performance.

Early Termination of Executive Board Duties

In the event of any early termination of Executive Board duties, the employment contracts of Executive Board members provide for severance to be capped to a maximum of two annual salaries.

Non-competition Clause

The employment contracts provide for a post-contractual non-competition clause for a duration of two years upon termination of employment with the company. In the event that this non-competition clause is not waived or is nullified, half of the remuneration last paid shall be granted for the duration of the non-competition period.

Fringe Benefits

Beyond the remuneration components stated above, the members of the Executive Board are each entitled to use a company car, reclaim expenses incurred on business travel and to be covered by accident insurance and D&O insurance as fringe benefits. D&O insurance provides for the application of a deductible or excess in the amount required by law.

Share-based Payment

As a rule, the remuneration policy for Executive Board members does not provide for the transfer of Sartorius AG shares as compensation for members. An exception to this policy rule was made in December 2014 for Dr. Kreuzburg, who was granted entitlement to receive share-based remuneration due to the third extension of his appointment as a member of the Executive Board and as its Chairman and CEO; please refer to Section 3 in this chapter.

2. Remuneration of the Executive Board Members in the Reporting Year

In 2017, the total remuneration for active service provided by all Executive Board members totaled €3,492K relative to €3,299K in 2016. Of this aggregate total, €1,892K accounted for non-performance-based components as "fixed remuneration" (2016: €1,672K) and €1,600K for variable performance-based components and multi-year components with a long-term incentive (2016: €1,627K). Furthermore, as part of the pension commitments to the Executive Board members, the pension service cost totaling €430K in the reporting year was expensed, following on €393K in the prior year.

Total Remuneration of the Executive Board Pursuant to § 314, Subsec. 1, No. 6, of the German Commercial Code (HGB)

€ in K	Executive Board (total)		Dr. Joachim Kreuzburg		Rainer Lehmann (as of March 1, 2017)		Jörg Pfirrmann (until Feb. 28, 2017)		Reinhard Vogt	
	2017	2016	2017	2016	2017	2016	2017	2016	2017	2016
Fixed remuneration	1,739	1,625	832	800	333	0	54	325	520	500
Fringe benefits ¹⁾	153	47	15	15	117	0	2	13	19	19
Fixed remuneration	1,892	1,672	847	815	450	0	56	338	539	519
Variable performance-based remuneration (1 year) ²⁾	759	850	363	418	145	0	24	170	227	262
Variable multi-year components w/ long-term incentive										
Consolidated net profit (3 years) ³⁾	375	371	184	182	0	0	79	78	112	111
Phantom stock plan (4 - 8 years) ⁴⁾	466	406	208	200	83	0	45	81	130	125
	1,600	1,627	755	800	228	0	148	329	469	498
Total remuneration	3,492	3,299	1,602	1,615	678	0	204	667	1,008	1,017

¹⁾ The amounts contributed to D&O insurance totaling €161K (2016: €171K) are not included as these refer to the executive bodies of all companies of the Sartorius Group and are not allocated to the individual insureds.

²⁾ Amount corresponds to actual target achievement

³⁾ Amount corresponds to actual target achievement of the plan in which a fiscal year ended; i.e., for 2017, consolidated net profits for 2015–2017 (2016: consolidated net profits for 2014–2016)

⁴⁾ Fair value at the time granted

As part of the remuneration component based on the consolidated net profit of three consecutive fiscal years, each Executive Board member receives a partial compensation payment of 50% of his respective target achievement for the first fiscal year under review. Once the total target achievement has been determined after the third fiscal year, final payment is then effected by deducting the particular partial payment already made. The amounts of the partial payments made in total at the end of the reporting year are shown as follows:

	2017 € in K	2016 € in K
Balance as of Jan. 1 of a fiscal year	387	311
Partial payments deducted	-165	-146
Partial payments effected	169	222
Balance as of Dec. 31 of a fiscal year	391	387

3. Disclosures on Share-based Payments

For multi-year components with a long-term incentive, the phantom stock plan must be generally classified as share-based payment just as is the share-based payment agreed to be granted to Dr. Kreuzburg in December 2014 in connection with the third extension of his appointment as a member of the Executive Board and its Chairman and CEO. Since December 18, 2015, Dr. Kreuzburg has held 100,000 ordinary shares and 100,000 preference shares. These shares transferred to him are subject to a holding period that will end on November 10, 2019. Should Dr. Kreuzburg leave the company prior to November 11, 2019, at his own request, he shall be required to transfer half of the shares granted to him back to the company.

The amount resulting since December 16, 2014, for the shares granted are to be spread as an employee benefits expense over the full vesting period and recognized as such in profit or loss. In fiscal 2017, an amount of €1,241 K was accordingly recognized as an employee benefits expense resulting from the grant of shares.

The employee benefits expense recognized in profit or loss in connection with the share-based payments is summarized as follows:

	2017 € in K	2016 € in K
Executive Board (total)		
Share-based payments	1,607	1,746
Phantom stock units	366	390
Shares granted	1,241	1,356

	2017 € in K	2016 € in K
Dr. Joachim Kreuzburg		
Share-based payments	1,404	1,548
Phantom stock units	163	192
Shares granted	1,241	1,356

	2017 € in K	2016 € in K
Rainer Lehmann (as of March 1, 2017)		
Share-based payments	65	0
Phantom stock units	65	0
Shares granted	0	0

	2017 € in K	2016 € in K
Jörg Pfirrmann (until Feb. 28, 2017)		
Share-based payments	36	78
Phantom stock units	36	78
Shares granted	0	0

	2017 € in K	2016 € in K
Reinhard Vogt		
Share-based payments	102	120
Phantom stock units	102	120
Shares granted	0	0

Disclosure of Phantom Stock Units

	Number of phantom stock units	Price on assignment in €	Fair value when granted on Jan. 1 of the particular year € in K	Fair value at year-end on Dec. 31, 2016 € in K	Fair value at year-end on Dec. 31, 2017 € in K	Paid in fiscal 2017 € in K	Change in value in fiscal 2017 € in K	Status
Dr. Joachim Kreuzburg								
Tranche for fiscal 2013	9,156	17.34	159	397	0	397	0	Paid out in 2017
Tranche for fiscal 2014	8,032	21.01	169	422	422	0	0	Not exercisable
Tranche for fiscal 2015	7,360	24.70	182	454	454	0	0	Not exercisable
Tranche for fiscal 2016	3,484	57.41	200	192	225	0	33	Not exercisable
Sum of the tranches from the previous years	28,032		710	1,465	1,101	397	33	
Tranche for fiscal 2017	2,950	70.51	208	0	163	0	-45	Not exercisable
Total sum of tranches	30,982		918	1,465	1,264	397	-12	
Rainer Lehmann								
Tranche for fiscal 2017	1,182	70.51	83	0	65	0	-18	Not exercisable
Total sum of tranches	1,182		83	0	65	0	-18	
Jörg Pfirrmann								
Tranche for fiscal 2013	3,960	17.34	69	172	0	172	0	Paid out in 2017
Tranche for fiscal 2014	3,452	21.01	73	181	181	0	0	Not exercisable
Tranche for fiscal 2015	3,140	24.70	78	194	194	0	0	Not exercisable
Tranche for fiscal 2016	1,416	57.41	81	78	91	0	13	Not exercisable
Sum of the tranches from the previous years	11,968		301	625	466	172	13	
Tranche for fiscal 2017	644	70.51	45	0	36	0	-9	Not exercisable
Total sum of tranches	12,612		346	625	502	172	4	
Reinhard Vogt								
Tranche for fiscal 2013	5,588	17.34	97	242	0	242	0	Paid out in 2017
Tranche for fiscal 2014	4,880	21.01	103	256	256	0	0	Not exercisable
Tranche for fiscal 2015	4,456	24.70	110	275	275	0	0	Not exercisable
Tranche for fiscal 2016	2,176	57.41	125	120	140	0	20	Not exercisable
Sum of the tranches from the previous years	17,100		435	893	671	242	20	
Tranche for fiscal 2017	1,844	70.51	130	0	102	0	-28	Not exercisable
Total sum of tranches	18,944		565	893	773	242	-8	

The number of phantom stock units granted as well as the particular grant prices were adjusted to the proportions following the stock split executed in 2016.

4. Pension Commitments

The retirement plan for Executive Board members provides for an old age and disability pension for Dr. Kreuzburg and for an old age pension for Messrs. Lehmann, Pfirrmann and Vogt. To cover such pensions, a benefit contribution amounting to one percent of each pensionable income and of each pensionable bonus is paid into a reinsurance policy. The benefit contribution for Dr. Kreuzburg is 10%; for Messrs. Pfirrmann and Vogt, 14% of their respective pensionable income; and for Mr. Lehmann, 9%, which equals their fixed remuneration.

If an Executive Board member elects to convert a portion of his salary to accrued retirement benefits by paying his own contribution into the reinsurance policy, Sartorius matches this by paying a corresponding additional benefit contribution on the reporting date. This amount matched by the company is 5% of the pensionable bonus earned by Dr. Kreuzburg and 7% of the same earned by Messrs. Pfirrmann and Vogt. This pensionable bonus is comprised of their respective one-year variable remuneration and of their respective multi-year remuneration based on the consolidated net profit. The amount of the retirement benefits that

Sartorius will pay later to each Executive Board member and his surviving dependents is dependent on the maturity payment of the insurance policy accrued up to the maturity date, including the policyholders' bonuses earned by the insurance company. An Executive Board member does not acquire any rights to the reinsurance policy; Sartorius shall be solely vested with such rights at all times.

Furthermore, an earlier pension agreement granted to Dr. Kreuzburg provides that he will receive a monthly pension of 70% of the basic salary of a German federal civil servant classified as grade 10 of salary class B for ministry officials according to the Federal Civil Service Remuneration Act (Bundesbesoldungsgesetz) in the respective version applicable. With each full year of service, 5% of his full pension is vested until after his full pension will have been reached after 20 years. Arrangements for pensions of surviving dependents basically provide for a widow's pension of 60% and an orphan's pension for each child amounting to 20% of his pension.

The projected pension payments, the present value of pension obligations and service cost are shown in the following table:

€ in K	Projected pension payment p.a.	Present value of the obligation (IFRS)		Service cost (IFRS)	
		Dec. 31, 2017	Dec. 31, 2016	2017	2016
Dr. Joachim Kreuzburg	238	2,989	2,741	258	233
Rainer Lehmann	26	27	0	30	0
Jörg Pfirrmann	84	442	375	43	64
Reinhard Vogt	40	604	475	99	96
	388	4,062	3,591	430	393

5. Disclosures Required by the German Corporate Governance Code (DCGK)

The following table shows the benefits granted for the year 2017, including the fringe benefits and the attainable maximum and minimum remuneration for the variable remuneration components in line with the requirements of the DCGK of lit. 4.2.5 of February 2017:

	Dr. Joachim Kreuzburg				Rainer Lehmann (as of March 1, 2017)				Jörg Pfirrmann (until Feb. 28, 2017)				Reinhard Vogt			
	2017 (min)	2017 (max)	2017	2016	2017 (min)	2017 (max)	2017	2016	2017 (min)	2017 (max)	2017	2016	2017 (min)	2017 (max)	2017	2016
Benefits granted € in K																
Fixed remuneration	832	832	832	800	333	333	333	0	54	54	54	325	520	520	520	500
Fringe benefits	15	15	15	15	117	117	117	0	2	2	2	13	19	19	19	19
Total non-performance-based remuneration	847	847	847	815	450	450	450	0	56	56	56	338	539	539	539	519
Variable performance-based remuneration (1 year)	0	499	416	400	0	200	167	0	0	32	27	163	0	312	260	250
Variable multi-year components w/ long-term incentive																
Consolidated net profit (2017 - 2019)	0	250	208		0	100	83		0	54	45		0	156	130	
Consolidated net profit (2016 - 2018)				200				0				81				125
Phantom stock plan 2017 (holding period 2017 - 2020)	0	520	208		0	208	83		0	113	45		0	325	130	
Phantom stock plan 2016 (holding period 2016 - 2019)				200				0				81				125
	847	2,116	1,679	1,615	450	958	783	0	56	255	173	663	539	1,332	1,059	1,019
Post-employment benefits	258	258	258	233	30	30	30	0	43	43	43	64	99	99	99	96
Total remuneration	1,105	2,374	1,937	1,848	480	988	813	0	99	298	216	727	638	1,431	1,158	1,115

The inflows of the various remuneration components in the reporting year are shown in the following table:

	Dr. Joachim Kreuzburg		Rainer Lehmann (as of March 1, 2017)		Jörg Pfirrmann (until Feb. 28, 2017)		Reinhard Vogt	
	2017	2016	2017	2016	2017	2016	2017	2016
Benefits received € in K								
Fixed remuneration	832	800	333	0	54	325	520	500
Fringe benefits ¹⁾	15	15	117	0	2	13	19	19
Total non-performance-based remuneration	847	815	450	0	56	338	539	519
Variable performance-based remuneration (1 year) ¹⁾	363	418	145	0	24	170	227	262
Variable multi-year components w/ long-term incentive								
Consolidated net profit (2015 - 2017) ¹⁾	184		0		79		112	
Consolidated net profit (2014 - 2016) ¹⁾		182		0		78		111
Phantom stock plan 2013 ²⁾	397		0		172		242	
Phantom stock plan 2012 ²⁾		365		0		160		223
	1,791	1,780	595	0	331	746	1,120	1,115
Post-employment benefits	258	233	30	0	43	64	99	96
Total remuneration	2,049	2,013	625	0	374	810	1,219	1,211

¹⁾ Amount equal to actual target achievement

²⁾ Paid out in the fiscal year

6. Main Features of the Remuneration Plan for the Supervisory Board

The remuneration for Supervisory Board members is defined in the Articles of Association of Sartorius AG and comprises fixed remuneration, meeting attendance fees and reimbursement of out-of-pocket expenses. Members serving as chairperson and vice chairperson of the Supervisory Board receive higher fixed remuneration.

Members and chairpersons of Supervisory Board committees are entitled to receive additional annual fixed amounts and meeting attendance fees and reimbursement of their out-of-pocket expenses. These amounts do not apply in relation to the Nomination Committee or to the committee pursuant to Section 27, Subsection 3, of the German Codetermination Law (MitBestG).

7. Remuneration of the Supervisory Board Members

	2017 € in K	2016 € in K
Remuneration for the Supervisory Board Members		
Total remuneration	922	957
Fixed remuneration	587	600
Compensation for committee work	80	80
Meeting attendance fee	165	184
Total remuneration for the Sartorius Stedim Biotech subgroup	90	93
Remuneration from Sartorius Stedim Biotech GmbH, Göttingen	10	38
Remuneration from Sartorius Stedim Biotech S.A., Aubagne	80	55

	2017 € in K	2016 € in K
Dr. Lothar Kappich (Chairman as of July 19, 2017)		
Total remuneration	142	49
Fixed remuneration	76	40
Compensation for committee work	13	0
Meeting attendance fee	15	9
Vergütung von der Sartorius Stedim Biotech S.A., Aubagne	38	0

	2017 € in K	2016 € in K
Manfred Zaffke¹⁾ (Vice Chairman)		
Total remuneration	118	125
Fixed remuneration	81	80
Compensation for committee work	16	16
Meeting attendance fee	21	29

	2017 € in K	2016 € in K
Annette Becker¹⁾		
Total remuneration	51	49
Fixed remuneration	40	40
Meeting attendance fee	11	9

	2017 € in K	2016 € in K
Uwe Bretthauer¹⁾		
Total remuneration	77	85
Fixed remuneration	40	40
Compensation for committee work	16	16
Meeting attendance fee	21	29

	2017 € in K	2016 € in K
Michael Dohrmann¹⁾		
Total remuneration	51	49
Fixed remuneration	40	40
Meeting attendance fee	11	9

	2017 € in K	2016 € in K
Dr. Daniela Favocchia (as of April 6, 2017)		
Total remuneration	36	0
Fixed remuneration	30	0
Meeting attendance fee	6	0

	2017 € in K	2016 € in K
Petra Kirchhoff		
Total remuneration	51	49
Fixed remuneration	40	40
Meeting attendance fee	11	9

	2017 € in K	2016 € in K
Karoline Kleinschmidt¹⁾		
Total remuneration	51	49
Fixed remuneration	40	40
Meeting attendance fee	11	9

	2017 € in K	2016 € in K
Dr. Guido Oelkers (as of November 11, 2017)		
Total remuneration	8	0
Fixed remuneration	6	0
Meeting attendance fee	2	0

	2017 € in K	2016 € in K
Ilke Hildegard Panzer (as of April 6, 2017)		
Total remuneration	36	0
Fixed remuneration	30	0
Meeting attendance fee	6	0

	2017 € in K	2016 € in K
Prof. Dr. Thomas Scheper		
Total remuneration	51	48
Fixed remuneration	40	40
Meeting attendance fee	11	8

	2017 € in K	2016 € in K
Prof. Dr. Klaus Trützschler		
Total remuneration	78	71
Fixed remuneration	40	40
Compensation for committee work	20	16
Meeting attendance fee	18	15

1) The employee representatives declared that they donate their Supervisory Board remuneration to the foundation Hans-Böckler-Stiftung according to the guidelines of the German Trade Union Association.

Beyond their Supervisory Board remuneration, the employee representatives who are employees within the Sartorius Group receive compensation that is not related to their service on the Supervisory Board.

Former Supervisory Board Members

	2017 € in K	2016 € in K
Prof. Dr. Dres. h.c. Arnold Picot (Chairman until July 9, 2017)		
Total remuneration	139	266
Fixed remuneration	62	120
Compensation for committee work	13	24
Meeting attendance fee	12	29
Total remuneration for the Sartorius Stedim Biotech subgroup	52	93
Remuneration from Sartorius Stedim Biotech GmbH, Göttingen	10	38
Remuneration from Sartorius Stedim Biotech S.A., Aubagne	42	55

	2017 € in K	2016 € in K
Dr. Dirk Basting (until April 6, 2017)		
Total remuneration	14	49
Fixed remuneration	11	40
Meeting attendance fee	3	9

	2017 € in K	2016 € in K
Prof. Dr. Gerd Krieger (until April 6, 2017)		
Total remuneration	19	68
Fixed remuneration	11	40
Compensation for committee work	2	8
Meeting attendance fee	6	20

8. Remuneration of Former Managing Directors

	2017 € in K	2016 € in K
Remuneration of Former Managing Directors		
Remuneration of former managing directors and members of the Executive Board as well as their surviving dependents	871	498
Retirement benefits and pension obligations to former managing directors and members of the Executive Board as well as their surviving dependents	8,098	7,485

Any circumstances beyond the disclosures made above and required to be reported according to Section 289, Subsection 4, and Section 315, Subsection 4, of the German Commercial Code "HGB" do not exist or are unknown.