

Consolidated Financial Statements and Notes

04

Statement of Profit or Loss and Other Comprehensive Income

	Notes	2018 12 months € in K	2017 12 months € in K
Sales revenue	[9]	1,212,152	1,081,033
Cost of sales		-582,589	-526,208
Gross profit on sales		629,563	554,825
Selling and distribution costs		-215,208	-195,241
Research and development costs		-60,616	-53,210
General administrative expenses		-67,004	-61,685
Other operating income and expenses	[11]	13,500	-22,987
Earnings before interest and taxes (EBIT)		300,234	221,702
Financial income	[12]	5,349	9,500
Financial expenses	[12]	-21,035	-10,589
Financial result		-15,685	-1,089
Profit before tax		284,549	220,613
Income taxes	[13]	-74,590	-56,849
Net profit for the period		209,959	163,763
Attributable to:			
Equity holders of Sartorius Stedim Biotech		208,052	161,088
Non-controlling interest	[22]	1,907	2,676
Earnings per share (€)	[15]	2.26	1.75
Diluted earnings per share (€)	[15]	2.26	1.75

Other Comprehensive Income

	Notes	2018 12 months € in K	2017 12 months € in K
Net profit for the period		209,959	163,763
Cash flow hedges	[36]	-11,547	26,299
of which effective portion of changes in fair value		-20,016	26,901
of which reclassified to profit or loss		8,469	-602
Income tax on cash flow hedges	[19]	3,463	-7,889
Foreign currency translation differences		6,433	-27,034
Items that are or may be reclassified subsequently to profit or loss		-1,651	-8,624
Remeasurements of the net defined benefit liabilities	[23]	469	-667
Income tax on remeasurements of the net defined benefits liabilities	[19]	-160	481
Items that will not be reclassified to profit or loss		309	-186
Other comprehensive income after tax		-1,342	-8,810
Total comprehensive income		208,617	154,953
Attributable to:			
Equity holders of Sartorius Stedim Biotech		206,673	152,335
Non-controlling interest		1,943	2,619

Statement of Financial Position

	Notes	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Non-current assets			
Goodwill	[16]	384,695	386,045
Other Intangible Assets	[16]	177,492	172,906
Property, plant and equipment	[17]	435,980	331,295
Financial Assets		5,644	11,808
Other Assets		632	18
Deferred tax assets	[19]	14,490	10,989
		1,018,932	913,060
Current assets			
Inventories	[20]	252,002	185,604
Trade receivables	[28]	220,231	211,616
Other financial assets	[29]	22,036	20,049
Current tax assets		12,950	24,465
Other assets		21,334	16,561
Cash and cash equivalents	[27]	23,975	32,552
		552,529	490,845
Total assets		1,571,461	1,403,905
Equity			
Equity attributable to SSB S.A. shareholders		1,036,398	872,028
Issued capital	[21]	18,436	18,436
Capital reserves		231,526	231,526
Retained earnings (including net profit)		786,436	622,066
Non-controlling interest		8,476	7,426
		1,044,874	879,454
Non-current liabilities			
Pension provisions	[23]	35,595	35,446
Other provisions	[24]	2,877	3,103
Loans and borrowings	[30]	43,125	46,250
Finance lease liabilities		14,976	15,829
Other financial liabilities	[31]	8,889	30,392
Deferred tax liabilities	[19]	39,150	42,987
		144,612	174,007
Current liabilities			
Provisions	[24]	12,283	9,558
Trade payables	[32]	154,568	117,959
Loans and borrowings	[30]	89,817	95,933
Finance lease liabilities		1,717	1,658
Other financial liabilities	[33]	14,312	34,964
Employee benefits		39,335	31,895
Current tax liabilities		33,070	31,791
Other liabilities		36,873	26,686
		381,975	350,444
Total equity and liabilities		1,571,461	1,403,905

Statement of Cash Flows

	Notes	2018 12 months € in K	2017 12 months € in K
Profit before tax		284,549	220,613
Financial result	[12]	15,685	1,089
Earnings before interest and taxes (EBIT)		300,234	221,702
Depreciation amortization of fixed assets	[16][17]	60,914	50,678
Increase decrease in provisions	[23][24]	2,361	655
Income taxes paid	[13]	-65,537	-49,248
Other non-cash items		-32,909	3,489
Gross cash flows from operating activities		265,062	227,276
Increase decrease in receivables and other assets	[28][29]	-21,604	-46,316
Increase decrease in inventories	[20]	-64,962	-21,091
Increase decrease in liabilities	[31][32][33]	48,793	14,819
Net cash flow from operating activities		227,289	174,689
Acquisitions of intangible and tangible assets	[16][17]	-176,540	-126,826
Net cash flow from investing activities		-176,540	-126,826
Payments for acquisitions of consolidated subsidiaries and other business operations; net of cash acquired	[8]	0	-68,100
Net cash flow from investing activities and acquisitions		-176,540	-194,926
Interest received	[12]	2,569	290
Interest paid and other financial charges	[12]	-8,995	-1,914
Dividends paid to:			
- Shareholders of Sartorius Stedim Biotech SA		-42,403	-38,713
- Non-controlling interest		-783	-704
Gross cash flows from financing activities		-49,612	-41,041
Proceeds from loans and borrowings raised	[6][30]	12,829	60,994
Repayments of loans and borrowings	[6][30]	-22,811	-3,333
Net cash flow from financing activities		-59,594	16,620
Net increase decrease in cash and cash equivalents		-8,845	-3,617
Cash and cash equivalents at the beginning of the period		32,552	34,756
Net effect of currency translation on cash and cash equivalents		268	1,414
Cash and cash equivalents at the end of the period		23,975	32,552

The Notes to the Consolidated Financial Statements are an integral part of these statements.

Statement of Changes in Equity

€ in K	Issued capital	Capital reserves	Hedging reserves	Pension reserves	Retained earnings	Foreign currency translation reserves	Group equity	Non-controlling interest	Total equity
Balance at Jan. 1, 2017	18,083	231,526	-6,961	-10,983	490,275	36,064	758,004	5,551	763,555
Net profit for the period	0	0	0	0	161,088	0	161,088	2,676	163,763
Cash flow hedges	0	0	26,299	0	0	0	26,299	0	26,299
Remeasurements of the net defined benefit liabilities	0	0	0	-667	0	0	-667	0	-667
Foreign currency translation differences	0	0	0	0	0	-26,977	-26,977	-57	-27,034
Deferred taxes	0	0	-7,889	481	0	0	-7,408	0	-7,408
Other comprehensive income for the period	0	0	18,410	-186	0	-26,977	-8,753	-57	-8,810
Total comprehensive income	0	0	18,410	-186	161,088	-26,977	152,335	2,619	154,953
Dividends	0	0	0	0	-38,713	0	-38,713	-704	-39,417
Changes in non-controlling interest	0	0	0	0	40	0	40	-40	0
Other changes	353	0	0	0	9	0	362	0	362
Balance at Dec. 31, 2017	18,436	231,526	11,449	-11,169	612,699	9,087	872,027	7,426	879,453
Adjustment on adoption of IFRS 9					321		321		321
Balance at Jan. 1, 2018	18,436	231,526	11,449	-11,169	613,020	9,087	872,349	7,426	879,775
Net profit for the period	0	0	0	0	208,052	0	208,052	1,907	209,959
Cash flow hedges	0	0	-11,547	0	0	0	-11,547	0	-11,547
Remeasurements of the net defined benefit liabilities	0	0	0	469	0	0	469	0	469
Foreign currency translation differences	0	0	0	0	0	6,396	6,396	37	6,433
Deferred taxes	0	0	3,463	-160	0	0	3,303	0	3,303
Other comprehensive income for the period	0	0	-8,084	309	0	6,396	-1,379	37	-1,342
Total comprehensive income	0	0	-8,084	309	208,052	6,396	206,673	1,943	208,617
Dividends	0	0	0	0	-42,403	0	-42,403	-783	-43,186
Other changes		0	0	0	-221	0	-221	-110	-331
Balance at December 31, 2018	18,436	231,526	3,365	-10,860	778,448	15,483	1,036,398	8,476	1,044,874

Notes to the Financial Statements

1. General Information

Sartorius Stedim Biotech is a leading international supplier of products and services that enable the biopharmaceutical industry to develop and manufacture drugs safely and efficiently. As a total solutions provider, Sartorius Stedim Biotech offers a portfolio covering nearly all steps of biopharmaceutical manufacture. The company focuses on single-use technologies and value-added services to meet the rapidly changing technology requirements of the industry it serves. With its own manufacturing and R&D sites in Europe, North America and Asia and an international network of sales companies, Sartorius Stedim Biotech has a global reach.

Headquartered in Aubagne, France, Sartorius Stedim Biotech S.A. is listed on the Euronext Paris (ISIN code: FR 0000053266).

Sartorius Stedim Biotech S.A.'s ultimate parent company is Sartorius AG, headquartered in Goettingen, Germany, and listed at several German stock exchanges (ISIN codes: 0007165607 ordinary shares, 0007165631 preference shares).

In compliance with the European Regulation 1606/2002 of July 19, 2002, requiring listed companies to use International Accounting Standards, the consolidated financial statements of the Sartorius Stedim Biotech Group for the year ended December 31, 2018, are compliant with the Standards and Interpretations IFRS and IFRIC of the IASB as adopted by the European Union, which are available at the following site:

https://ec.europa.eu/commission/index_fr.

The consolidated financial statements are prepared in euros. Unless otherwise specified, all amounts are disclosed in thousands of euros (abbreviated as € in K). In some cases, the sum of the figures given in this report may not precisely equal the stated totals and percentages may not be exact due to rounding.

These consolidated financial statements were approved by the Board of Directors on February 7, 2019 and will be submitted for approval by the Shareholders' Meeting on March 26, 2019.

2. Effects of New Financial Reporting Standards

The following major new accounting rules were applicable for the first time to the present consolidated financial statements of the Group:

– IFRS 9, Financial Instruments

This standard changes the accounting for financial instruments. For detailed information on these changes and the initial application of IFRS 9, see note 26.

– IFRS 15, Revenue from Contracts with Customers

– Clarifications to IFRS 15, Revenue from Contracts with Customers

This standard defines when and in which amount revenue is recognized. For detailed information on the related changes and the initial application of IFRS 15, see note 9.

The following new accounting rules were applicable for the first time to the present financial statements and had no material impact on the presentation of the company's financial position and financial performance:

– Annual Improvements to IFRSs 2014 - 2016 Cycle, Amendments to IFRS 1 and IAS 28

The amendments are related to the first-time application of IFRS (IFRS 1) as well as the option for venture capital organizations and similar organizations to measure investments in associates or joint ventures at fair value through profit or loss (IAS 28).

– Amendments to IFRS 2, Classification and Measurement of Share-based Payment Transactions

The amendments include clarifications with regard to the accounting for share-based payment transactions.

– Amendments to IAS 40, Transfers of Investment Property

The amendments intend to clarify the guidance regarding transfers to, or from, investment properties.

– IFRIC 22, Foreign Currency Transactions and Advance Consideration

The interpretation addresses a question in relation to IAS 21. It clarifies which exchange rate is used for foreign currency transactions that involve advance payments received or made.

The following standards, interpretations and amendments were not yet applied to the consolidated financial statements of the reporting year as they had not yet been adopted by the EU or their application was not mandatory for 2018:

Standard Interpretation	Title	Applicable for financial years from ¹⁾	Endorsement by the EU Commission
IFRS 14	Regulatory Deferral Accounts	January 1, 2016	No
IFRS 16	Leases	January 1, 2019	Yes
IFRIC 23	Uncertainty over Income Tax Treatments	January 1, 2019	Yes
Amendments to IFRS 9	Prepayment Features with Negative Compensation	January 1, 2019	Yes
Amendments to IFRS 28	Long-term Interests in Associates and Joint Ventures	January 1, 2019	No
Amendments to IFRS 3, IFRS 11, IAS 12 and IAS 23	Annual Improvements to IFRSs 2015 - 2017 Cycle (issued in Dec. 2017)	January 1, 2019	No
Amendments to IAS 19	Plan Amendment, Curtailment or Settlement	January 1, 2019	No
Amendments to IAS 1 and IAS 8	Definition of Material	January 1, 2020	No
Various Standards / Framework	Amendments to References to the Conceptual Framework in IFRS Standards	January 1, 2020	No
Amendments to IFRS 3	Definition of a Business	January 1, 2020	No
IFRS 17	Insurance Contracts	January 1, 2021	No
Amendments to IFRS 10 and IAS 28	Sale or Contribution of Assets between an Investor and its Associate or Joint Venture	n/a	No

¹⁾ These are required to be applied once they are endorsed by the EU Commission. The dates mentioned above are those required by the standard themselves (IASB effective dates).

IFRS 16, Leases

The Group did not yet apply IFRS 16, Leases. The application of the standard is mandatory for periods from 2019 onwards. IFRS 16 introduces a standardized accounting model according to which leases are generally to be recognized on the lessee's balance sheet.

A lessee recognizes a right-of-use asset representing its right to use a lease asset, as well as a liability resulting from the lease, which represents its obligation to make lease payments. There are exemptions for short-term leases and leases of low-value assets. Accounting for the lessor is comparable to that of the current standard; i.e., lessors continue to classify leases as financial or operating leases.

The Group has analyzed the effects of this new standard on the consolidated financial statements during the year 2018. The Group did not apply the standard early. The Group does not plan to apply the standard retrospectively in accordance with IAS 8, but rather plans to recognize the effect from initially applying IFRS 16 in retained earnings at the date of

initial application. The Group will use the exemptions for short-term leases and leases of low-value assets and recognize the corresponding lease payments as an expense generally on a straight-line basis over the particular lease term.

IFRS 16 will lead to an increase in fixed assets and financial liabilities. Based on its present level of knowledge, the Group does not expect any significant impact overall on its most important key figures, such as equity ratio or underlying EBITDA.

For example, on the basis of a differentiated analysis of the operating lease commitments (see also note 18) conducted in the course of the preparation of the present consolidated financial statements, total assets would increase by approx. €35 million as of December 31, 2018. This would result in a reduction of the equity ratio by approx. 1.5%. On this basis, the EBITDA margin would increase by about 0.5 - 1% for the year 2018.

3. Significant Accounting Policies

Basis of Preparation

The consolidated financial statements of the Group are based on the principle of the historical cost of acquisition, construction or production, with the exception of the items reflected at fair value, such as derivative financial instruments.

Consolidation

The consolidated financial statements of the Sartorius Stedim Biotech Group include the annual financial statements of all companies, which are controlled directly or indirectly by Sartorius Stedim Biotech S.A. In terms of IFRS 10, Consolidated Financial Statements, the Group Sartorius Stedim Biotech controls an entity when it is exposed to, or has rights to, variable returns from its involvement with the entity and has the ability to affect those returns through its power over the entity.

Such entities are included in the consolidated financial statements from the time when Sartorius Stedim Biotech S.A. or its subsidiaries obtain such control until the date on which control ceases.

Subsidiaries have been included on the basis of their annual financial statements for the same reporting period as the parent company, using uniform Group recognition and measurement methods.

All intragroup assets and liabilities, equity, income, expenses and cash flows relating to transactions between members of the Group are eliminated on consolidation.

Business Combinations

Business combinations are accounted according to the acquisition method. The identifiable acquired assets and assumed liabilities are generally recorded at fair value on the date of acquisition.

For significant acquisitions, the purchase price allocation is carried out with assistance from independent third-party valuation specialists. The valuations are based on the information available at the acquisition date.

The Group determines goodwill at the acquisition date as:

- the fair value of the consideration transferred; and
- the amount recognized for any non-controlling interest in the acquiree; and
- if the business combination is carried out in stages, the fair value of any previously held equity interest in the acquiree; less
- the net recognized amount for the identifiable assets acquired and liabilities assumed.

When the difference is negative, the purchase gain is recognized immediately in income. Expenses directly related to business combinations are recorded in the profit or loss as they are incurred.

Foreign Currency Transactions

The presentation currency of the consolidated financial statements of the Sartorius Stedim Biotech Group is the euro (financial statements presented in thousands of euros). In the financial statements of each company, transactions denominated in foreign currencies have been translated into the functional currency of the subsidiary at the exchange rate applicable on the date of the transaction. Monetary assets and debts denominated in a foreign currency have been translated at the exchange rate on the balance sheet date. Rate gains and losses have been recognized in profit or loss for the period.

Translation of financial statements prepared in foreign currencies

Subsidiaries' financial statements prepared in foreign currencies have been translated pursuant to IAS 21, The Effects of Changes in Foreign Exchange Rates, in accordance with the concept of a functional currency. Foreign subsidiaries have been regarded as independent subdivisions of the Sartorius Stedim Biotech Group. The assets (including goodwill) and liabilities of the entities that have a functional currency different from the presentation currency are translated at the exchange rate prevailing at the balance sheet date. The incomes, expenses, and cash flows of these entities have been translated using the average rate for the year, to the extent that this rate represents an approximate value of exchange rates used as of the date of the transaction in the absence of significant fluctuations. Resulting translation differences are recognized in other comprehensive income.

For long-term loans for which settlement is neither planned nor likely in the foreseeable future, the Group applies the principle of "net investment in a foreign operation." Exchange differences resulting from these loans are recognized in other comprehensive income in accordance with IAS 21.32.

The exchange rates for major currencies against the euro were considered as follows:

For 1 €	Year-end exchange rates		Average exchange rates	
	2018	2017	2018	2017
USD	1.14500	1.19930	1.18129	1.12955
GBP	0.89453	0.88723	0.88465	0.87670
JPY	125.85000	135.01000	130.40242	126.70218
CHF	1.12690	1.17020	1.15517	1.11173
INR	79.72980	76.60550	80.69389	73.52751
KRW	1277.93000	1279.61000	1299.12464	1276.62397
CNY	7.87510	7.80440	7.80666	7.62790

4. Use of Judgments and Estimates

During the preparation of consolidated financial statements, management uses estimates and assumptions based on their best knowledge of the current and future situation. However, actual results may differ from these estimates. These estimates and assumptions are revised on a regular basis, and the impact of changes in estimates is recognized prospectively.

In addition, Group management exercises its judgment in defining the accounting treatment of specific transactions when the existing Standards and Interpretations do not specifically treat the accounting problems concerned.

Assumptions and estimates primarily concern the following topics:

Business Combinations

The accounting for business combinations requires that the assets acquired and liabilities assumed be recorded at their respective fair values on the date the Group obtains control. The application of the acquisition method requires estimates and assumptions to be made, especially concerning the fair values of the acquired intangible assets, property, plant and equipment and the liabilities assumed at the acquisition date, and the useful lives of the acquired intangible assets, property, plant and equipment.

These measurements are based to a large extent on anticipated cash flows. If actual cash flows vary from those used in calculating fair values, this may materially affect the Group's future results of operations.

Impairment of Assets

The book values (carrying amounts) of property, plant and equipment and intangible assets are subject to impairment testing if there is an indication of impairment and at least once a year for assets with an indefinite useful life or not yet available for use in accordance with IAS 36, Impairment of Assets. When an asset is tested, the recoverable amount of the asset is estimated. The recoverable amount of an asset or a cash-generating unit (CGU) is the higher of its fair value – less costs to sell the asset or its CGU – and its value in use. In the event the individual asset's recoverable amount cannot be estimated, the recoverable amount of the asset's CGU is estimated.

If the estimated recoverable amount of an asset (or a CGU) goes below its book value (carrying amount), this carrying amount is reduced to the recoverable amount (impairment allocated in priority to goodwill).

If the causes of the asset impairment are removed, the book value of the asset (or the CGU) is credited to the newly estimated recoverable amount. However, the book value increase is limited to the value that the asset (or CGU) would have had if no asset impairment loss had been recognized in previous financial years.

The calculation of the value in use is generally based on discounted cash flow methods using cash flow projections up to five years. These projections take into account past experience and represent management's best estimate about future sales revenue and cost developments. Cash flows after the planning period are extrapolated using individual growth rates. Key assumptions on which management has based its determination of the value in use include estimated growth rates, weighted average cost of capital and tax rates. These estimates can have a material impact on the respective values and ultimately the amount of any impairment.

Intangible Assets

The capitalization of self-constructed intangible assets also includes a significant level of judgment, e.g. the evaluation of feasibility of a development project, the expected market prospects and the determination of useful lives.

Employee Benefits – Pension Provisions

Obligations for pension and other post-employment benefits are determined in accordance with actuarial valuations. These valuations rely on key assumptions including discount rates, expected salary increases and mortality rates. The discount rate assumptions are determined by reference to yields on high-quality corporate bonds of appropriate duration and currency at the end of the reporting period.

Due to changing market and economic conditions the underlying key assumptions may differ from actual developments and may lead to significant changes in pension and other post-employment benefit obligations.

Such differences are recognized in other comprehensive income in the period in which they occur. For a sensitivity analysis, see note 23, Pension and Employee Benefits Provisions.

Provisions, Contingent Liabilities and Contingent Assets

Provisions are recognized for legal or constructive obligations that exist as of the balance sheet date. To determine the amount of the obligations, certain estimates and assumptions have to be applied, including the determination of the probability and the amount of future outflows of resources. Typically, significant estimates are involved in the determination of provisions related to onerous contracts, warranty costs, asset retirement obligations and legal proceedings.

Income Taxes

The Group operates in various tax jurisdictions and therefore has to determine tax positions under respective local tax laws and tax authorities' views which can be complex and subject to different interpretations of taxpayers and local tax authorities. Deferred tax assets have to be recognized for all deductible temporary differences and unused tax losses to the extent that it is probable that future taxable profit will be available against which the deductible temporary differences and unused tax losses can be utilized. As future developments are uncertain and partly beyond management's control,

assumptions are necessary to estimate future taxable profits as well as the period in which deferred tax assets will be recovered.

Estimates are revised in the period in which there is sufficient evidence to revise the assumption. If management considers it probable that all or a portion of a deferred tax asset cannot be realized, the corresponding amount is not recorded as an asset.

Fair Value Measurement

A number of the Group's accounting policies and disclosures may require the measurement of fair values, for both financial and non-financial assets and liabilities, including Level 3 fair values (unobservable inputs).

If third party information, such as broker quotes or pricing services, is used to measure fair values, then management assesses the evidence obtained from the third parties to support the conclusion that these valuations meet the requirements of IFRS, including the level in the fair value hierarchy in which the valuations should be classified.

When measuring the fair value of an asset or a liability, the Group uses observable market data as far as possible. If the inputs used to measure the fair value of an asset or a liability fall into different levels of the fair value hierarchy, then the fair value measurement is categorized in its entirety in the same level of the fair value hierarchy as the lowest level input that is significant to the entire measurement.

5. Operating Segments

According to IFRS 8, Operating Segments the identification of reportable operating segments is based on the "management approach"; i.e. the segments are defined analogously to the internal financial reporting of an entity. Therefore, an area of activity is to be considered an operating segment if its business activities may result in revenues and expenses, its operating results are regularly reviewed by the entity's chief operating decision maker (= the Executive Members of the Board of Directors) and discrete financial information is available in its internal reporting. Internal control and reporting within Sartorius Stedim Biotech is based on the approach of operating as a "total solution provider" for its customers. Accordingly, there is only one single segment to be identified for Sartorius Stedim Biotech, driven by the product and customer perspective: Biopharm.

The key performance indicator of the operating segment of the Sartorius Stedim Biotech Group is the so-called "underlying EBITDA", as the board monitors this performance measure at a consolidated level and they believe this measure is relevant to an understanding of the Group's financial performance.

EBITDA corresponds to earnings before interest, taxes, depreciation and amortization; "underlying EBITDA" means EBITDA adjusted for extraordinary items. In this connection, extraordinary items are expenses and income that are of an exceptional or a one-time nature and accordingly distort the sustainable profitability of a segment and have a material impact on the net worth, financial position and earnings of the Group. Examples of such items are restructuring

expenses, large Group projects as well as proceeds or losses from the disposal, sale or other transfer of financial assets or of property, plant and equipment, provided that these are not of a recurrent nature.

Underlying EBITDA is not a defined performance measure in IFRS. The Group's definition of underlying EBITDA may not be comparable with similarly titled performance measures and disclosures by other entities.

Segment assets and segment liabilities are not analyzed on a regular basis to the chief operating decision maker and are therefore not part of the segment report.

€ in K	Biopharm			Group		
	2018	2017	Change	2018	2017	Change
Sales revenue	1,212,152	1,081,033	12%	1,212,152	1,081,033	12%
Underlying EBITDA	342,430	294,912	16%	342,430	294,912	16%
as a % of sales revenue	28.2%	27.3%		28.2%	27.3%	
EBIT	300,234	221,702	35%	300,234	221,702	35%
as a % of sales revenue	24.8%	20.5%		24.8%	20.5%	
Acquisitions of intangible and tangible assets	176,945	136,734	29%	176,945	136,734	29%
as a % of sales revenue	14.6%	12.6%		14.6%	12.6%	

Reconciliation of Segment Profit or Loss:

€ in K	2018	2017
Underlying EBITDA of the segment	342,430	294,912
Depreciation and amortization	-54,918	-50,645
Extraordinary effects	12,722	-22,564
EBIT	300,234	221,702
Financial result	-15,685	-1,089
Profit before tax	284,549	220,613

Supplementary Information by Region

To provide additional information required by IFRS 8, the table below presents the supplementary information by geographical region. The key figures for non-current assets of the geographical areas refer to the company location, whereas sales revenue is reported according to the customer's location.

The non-current assets correspond to property, plant and equipment as well as to intangible assets (including goodwill) of the Group affiliates that are to be allocated to these various regions.

The amount of sales revenue with a single customer does not exceed 5% of the consolidated sales revenue (2018 and 2017).

€ in K	Sales revenue			Non-current assets
	2018	2017	2018	2017
EMEA	508,241	460,664	816,639	756,242
thereof Germany	144,602	130,454	340,599	296,567
thereof France	54,977	56,955	331,371	305,508
Americas	422,625	368,817	159,484	113,722
thereof USA	396,509	339,319	159,484	113,722
Asia Pacific	281,286	251,552	22,044	20,281
thereof China	83,904	66,978	1,474	1,526
thereof South Korea	72,824	76,585	7,502	7,109
Group	1,212,152	1,081,033	998,167	890,245

6. Statement of Cash Flows

The statement of cash flows shows the impact of cash inflows and outflows on the cash and cash equivalents of the Group. The cash flows are classified by operating, investing and financing activities according to IAS 7 (Statement of Cash Flows).

In this context cash equivalents are assets that can be converted into cash within a short maturity (generally less than three months). The amount considered in the

statement of cash flows is equal to the amount of cash and cash equivalents in the statement of financial position.

The impact resulting from the modified agreement with Lonza in the field of cell culture media posted in 2018 had no impact on the statement of cash flows (see note 11).

The following table summarizes the development of the liabilities arising from financing activities during the reporting period:

	Balance at Dec. 31, 2016 € in K	cash flows	Currency effects	other non-cash changes	Balance at Dec. 31, 2017 € in K
Loans and borrowings	84,052	58,347	-215	0	142,184
Finance lease liabilities	18,270	-686	-173	76	17,487
Liability for phantom units in connection with the AllPure acquisition	5,833	0	-706	155	5,282
Total financial liabilities from financing activities	108,155	57,661	-1,094	231	164,952

	Balance at Dec. 31, 2017 € in K	cash flows	Currency effects	other non-cash changes	Balance at Dec. 31, 2018 € in K
Loans and borrowings	142,184	-9,229	-12	0	132,943
Finance lease liabilities	17,487	-754	-40	0	16,693
Liability for phantom units in connection with the AllPure acquisition	5,282	0	250	3,206	8,739
Total financial liabilities from financing activities	164,953	-9,982	198	3,206	158,375

7. Scope of Consolidation

The 2018 financial statements of the following entities:

- TAP Biosystems (PHC) Ltd., UK
- TAP Biosystems Ltd., UK
- Distribo GmbH, Germany

were not included in the scope of consolidation, because the figures were of minor importance for assessing the financial position of the Group.

The sales revenue and total assets of the non-consolidated companies are below 1% of the Group figures.

The financial statements of the following companies have been included in the Group financial statements:

	Ownership in %
EMEA	
Sartorius Stedim Biotech S.A., Aubagne, France	Parent company
Sartorius Stedim Belgium N.V., Brussels, Belgium	100
Sartorius Stedim Nordics Oy, Helsinki, Finland	100
Sartorius Stedim Biotech GmbH, Goettingen, Germany	100
Sartorius Stedim Plastics GmbH, Goettingen, Germany	100
Sartorius Stedim North America Holding GmbH, Goettingen, Germany	100
Sartorius Stedim Systems GmbH, Guxhagen, Germany	100
Sartorius Stedim Cellca GmbH, Laupheim, Germany	100
Sartorius Stedim UK Ltd., Epsom, UK	100
Sartorius Stedim BioOutsource Ltd., Glasgow, UK	100
Sartorius Stedim Lab Ltd., Stonehouse, UK	100
TAP Biosystems Group Ltd., Royston, UK	100
TAP ESOP Management Ltd., Royston, UK	100
The Automation Partnership Cambridge Ltd., Royston, UK	100
Sartorius Stedim FMT S.A.S., Aubagne, France	100
Sartorius Stedim France S.A.S., Aubagne, France	100
Sartorius Stedim Aseptics S.A., Lourdes, France	100
Sartorius Stedim Ireland Ltd., Dublin, Ireland	100
Sartorius Stedim Italy S.p.A., Florence, Italy	100
Sartorius Stedim Netherlands B.V., Amersfoort, Netherlands	100
Sartorius Stedim Austria GmbH, Vienna, Austria	100
Sartorius Stedim Poland sp. z.o.o., Kostrzyn, Poland	100
Sartorius Stedim RUS, St. Petersburg, Russia	100
Sartorius Stedim Data Analytics AB, Umeå, Sweden	100
Sartorius Stedim Switzerland AG, Tagelswangen, Switzerland	100
Sartorius Stedim Spain S.A., Madrid, Spain	100
Sartorius Stedim Hungaria Kft., Budapest, Hungary	100
Sartorius Stedim Bioprocess S.A.R.L., M'Hamdia, Tunisia	100
Americas	
Sartorius Stedim Filters Inc., Yauco, Puerto Rico	100
Sartorius Stedim North America Inc., Wilmington, USA	100
Asia Pacific	
Sartorius Stedim Australia Pty. Ltd., Dandenong South, Victoria, Australia	100
Sartorius Stedim Biotech (Beijing) Co. Ltd., Beijing, China	100
Sartorius Stedim (Shanghai) Trading Co. Ltd., Shanghai, China	100
Sartorius Stedim India Pvt. Ltd., Bangalore, India	100
Sartorius Stedim Japan K.K., Tokyo, Japan	100
Sartorius Korea Biotech Co. Ltd., Seoul, South Korea	69
Sartorius Stedim Malaysia Sdn. Bhd., Kuala Lumpur, Malaysia	100
Sartorius Stedim Singapore Pte. Ltd., Singapore, Singapore	100

There are no associates or joint ventures included in the scope of consolidation, all companies are consolidated in full. The ownership rate equals the share in voting rights.

8. Business Combinations

None.

Notes to the Statement of Profit or Loss

9. Sales Revenue

Initial Application of IFRS 15, Revenue from Contracts with Customers

IFRS 15, Revenue from Contracts with Customers, defines a comprehensive model to determine when to recognize revenue and in which amount. It replaces existing guidelines for recognition of revenue, including IAS 18, Revenue; IAS 11, Construction Contracts; and IFRIC 13, Customer Loyalty Programmes.

The Group applied IFRS 15 for the first time as of January 1, 2018 on the basis of the modified retrospective method. Accordingly, the comparative period is presented in line with previous standards. Any effects of the initial application of IFRS 15 are recorded as an adjustment to the opening balance of retained earnings at the date of initial application. Furthermore, upon initial adoption, the Group has been applying IFRS 15 only to contracts that are not considered completed contracts at the date of initial application. Besides, the Group used the practical expedients offered regarding contract modifications that occurred prior to the date of initial application of IFRS 15 and did not retrospectively restate such contracts. As the impact of the new standard on the Group's consolidated financial statements is low, the use of this practical expedient is expected not to have a material impact.

There are no material effects as a result of the application of the new standard. However, as of the date of initial application, there are certain effects with regard to construction contracts in the company's project business. Sales revenue of approximately €5 million, which was recognized in the past, is not recognized over time according to IFRS 15, but rather at a point in time. This results from new criteria introduced by IFRS 15 for revenue recognition over time for contracts on the construction of customer-specific goods. The effects resulted in a balance sheet extension by about €2 million at the date of initial application. There were no effects to be recognized in equity. The effects of applying the modified retrospective method on the opening statement of financial position as of January 1, 2018 are shown in the following table.

Effects from the application of IFRS 15 on the opening balances of the statement of financial position as of Jan. 1, 2018	Carrying amount Dec. 31, 2017 € in mn	Adjustments on adoption of IFRS 15 € in mn	Carrying amount Jan. 1, 2018 € in mn ³⁾
Inventories	185.6	4.1	189.7
Amounts due from customers for contract work ¹⁾	7.0	-1.9	5.0
Trade receivables	204.6		204.6
Trade receivables	211.6	-1.9	209.7
Trade payables payments received for orders ²⁾	40.7	2.2	42.9
Trade payables	77.2		77.2
Trade payables	118.0	2.2	120.2

¹⁾ Contract assets according to IFRS 15.

²⁾ Contract liabilities according to IFRS 15.

³⁾ Carrying amount without adjustments for the new credit impairment model according to IFRS 9.

If the Group had applied the previous standards, IAS 18 and IAS 11, respectively, during the reporting period, sales revenue as well as cost of sales would have been higher by about €0.6 million while there would be no impact on the net result. Total assets and total liabilities would have been lower by about €2 million. The effects result from the revised criteria

for revenue recognition over time according to which revenue is now recognized at a point in time rather than over time for one project. The following table presents the impact on the consolidated statement of financial position as of December 31, 2018 that the continued application of the previous standards would have had in comparison with IFRS 15.

Effects from the application of IFRS 15 on the consolidated statement of financial position as of December 31, 2018	Carrying amount December 31, 2018 (IFRS 15) € in mn	Adjustments on adoption of IFRS 15 € in mn	Carrying amount December 31, 2018 (IAS 18, IAS 11) € in mn
Inventories	252.0	-5.1	246.9
Amounts due from customers for contract work ¹⁾	3.9	2.9	6.8
Trade receivables	216.3		216.3
Trade receivables	220.2	2.9	223.1
Trade payables payments received for orders ²⁾	51.0	-2.2	48.8
Trade payables	103.6		103.6
Trade payables	154.6	-2.2	152.4

¹⁾ Contract assets according to IFRS 15.

²⁾ Contract liabilities according to IFRS 15.

Other Disclosures

The revenues from contracts with customers according to IFRS 15 are disaggregated into geographical regions (see Segment Report, note 5). For details on the impairments on trade receivables and contract assets recognized in the reporting period see note 38.

The following table presents the opening balances as of January 1, 2018, and the balances as of December 31, 2018 for the Group's contract liabilities. The balances of trade receivables and contract assets are presented in note 28.

Line item in statement of financial position	Carrying amount as of January 1, 2018 € in k	Carrying amount as of December 31, 2018 € in k
Deferred revenue	10,153	16,235
Payments received on account of orders	42,949	51,000
Contract liabilities (total)	53,101	67,234

There were no extraordinary changes in the carrying amounts of the contract liabilities and contract assets in the reporting period. Revenue in the amount of 41,182 k euros was recognized in the reporting period that was included in the contract liability balance at the beginning of the reporting period.

The Group produces and sells instruments and consumables for customers in the Biopharm segment.

The Group satisfies its performance obligations depending on the goods to be transferred and the promised services. The far majority of the revenues from sales of products is recognized at a point in time when the customer obtains control over the goods. This is typically the case when the significant risks and rewards of ownership of the goods are transferred to the customer. Therefore, the point in time may vary

depending on the agreement with the individual customer.

For complex products that require installation at the customer's site, revenue is recognized upon formal customer acceptance. To a low extent, revenue is recognized over time in the customer-specific project business. In these cases, revenue is recognized according to the project progress which is measured based on the percentage of costs to date compared to the total estimated contract costs. The amount of actual costs incurred to date reflects the progress and the transfer of control to the customer appropriately as the Group has a right to a reimbursement of cost to date plus an appropriate margin, if the project is cancelled by the customer without cause.

Revenue from services is generally recognized when the services are performed or have been performed. When the services are performed continuously over a period of time, the Group recognizes the related revenue over time. In this case, revenue is generally recognized pro rata temporis in relation to the total contract period. Product sales are typically accompanied by the legally required warranty. Any material extended warranties are accounted for as separate performance obligations.

According to the general payment terms, customer payments are due in the short-term, typically within 30 days. To some extent, the Group obtains advance payments, e.g. to avoid credit risks. Therefore, the Group regularly has contract liabilities (payments received on account of orders). In addition, the Group recognizes contract liabilities in connection with service contracts (deferred revenues) when customers pay in advance.

There are no material effects from contracts with significant financing components. The Group uses the practical expedient regarding the existence of a significant financing component. This means that a financing component is only taken into consideration when the length of time between the transfer of goods or services and the receipt of consideration is expected to exceed one year and the effect is material. As of December 31, 2018, the Group has refund liabilities in the amount of 5,755k euros as a result from incentive agreements with customers.

The aggregate amount of the transaction price allocated to the performance obligations that are unsatisfied (or partially unsatisfied) at the end of the reporting period (orders on hand) amounts to 489.0million euros. The Group expects that these unsatisfied performance obligations will be satisfied to a large extent in 2019.

10. Functional Costs

The statement of profit or loss has been presented according to the "cost of sales format", i.e. expenses have been allocated to the relevant functions production, sales & marketing, research & development and general administration.

Expenses relating to cross-functional initiatives or projects are assigned to the respective functional costs based on an appropriate allocation principle.

The caption "cost of sales" includes the costs of products sold and the acquisition costs of merchandise sold. In addition to directly attributable expenses, such as raw materials and supplies, employee benefits expense and energy expenses, cost of sales also includes overhead, which can be allocated to the manufacturing area, and the corresponding depreciation and amortization.

The selling and distribution costs pertain, in particular, to the costs of the sales and marketing function, distribution and market research.

Research and development costs comprise the costs of research and product and process development, unless they are recognized as assets.

The item "general administrative expenses" mainly includes employee benefits expense and the cost of materials of the general administrative area.

All profit and loss items that cannot be allocated to one of the mentioned functional areas are recognized as other income and expenses. This includes essentially effects from translation of transactions in foreign currencies, sale of fixed assets, allowances on trade receivables and reorganization and other non-recurring expenses. Income from grants related to income is recognized as other income, when there is reasonable assurance that the conditions attached to the grants are complied with and the grants will be received. They are recognized systematically as income over the period in which the related costs are recorded.

Operating expenses by nature are presented in the Profit or Loss Statement by nature in note 14.

The material expense and personnel cost are as follows:

Raw Materials and Supplies

This caption consists of the following:

	2018 12 months € in K	2017 12 months € in K
Purchases consumed	237,889	235,762
Cost of purchased services	54,079	48,353
Total	291,968	284,114

Personnel Cost

This caption can be broken down as follows:

	2018 12 months € in K	2017 12 months € in K
Wages and salaries	274,115	244,438
Social security	59,877	54,481
Expenses for retirement benefits and pensions	6,861	8,088
Total	340,853	307,007

11. Other Operating Income and Expenses

	2018 12 months € in K	2017 12 months € in K
Currency translation gains	10,392	7,652
Extraordinary income	35,248	0
Income from the decrease in allowances for bad debts	475	780
Income from release of provisions and liabilities	353	973
Income from grants	4,235	8,013
Other income	1,921	1,139
Other operating income	52,624	18,557
Currency translation losses	-10,403	-14,762
Extraordinary expenses	-22,526	-22,564
Allowances for bad debts	-2,041	-801
Other expenses	-4,153	-3,416
Other operating expenses	-39,124	-41,543
Total other operating income and expenses	13,500	-22,987

The item reported as income from grants discloses the grants for expenses (essentially related to research and development projects).

Extraordinary items amounted to €12.7million (net) (previous year: €-22.6million). The extraordinary income in 2018 refers mainly to the derecognition of the liability for the acquisition of the cell culture media business as a consequence of the modified cooperation agreement with Lonza; extraordinary expenses essentially cover one-time expenses for strategic Group projects and integration and acquisition-related items.

12. Financial Result

	2018 12 months € in K	2017 12 months € in K
Interest and similar income	385	274
- of which from affiliated companies	126	18
Income from derivative financial instruments	2,798	801
Other financial income	2,167	8,425
Financial income	5,349	9,500
Interest and similar expenses	-3,344	-3,834
- of which from affiliated companies	-1,378	-1,814
Expenses for derivative financial instruments	-5,779	-1,706
Interest expense for pensions	-649	-585
Other financial expenses	-11,263	-4,465
Financial expenses	-21,035	-10,589
Total	-15,685	-1,089

The other financial income in 2017 (other financial expenses in 2018) includes mainly foreign exchange gains (losses) in connection with bank deposits, loans denominated in foreign currencies as well as the Lonza liability. Furthermore, in 2018 an increase in the liability for the phantom units in AllPure has been posted with an amount of approx. €3 mn (see note 31).

The interest expenses to affiliated companies are in connection with the loan granted by the Group's ultimate parent Sartorius AG (see also chapter 41).

13. Income Taxes

	2018 12 months € in K	2017 12 months € in K
Current income taxes	-78,330	-56,173
Deferred taxes	3,740	-676
Total	-74,590	-56,849

Current income taxes are determined based on the respective local taxable income of the period and local tax rules. In addition, current income taxes include adjustments for uncertain tax payments or tax refunds for periods not yet assessed. Changes in deferred tax assets and liabilities are included in income taxes except for changes recognized in other comprehensive income or equity.

Income taxes in France are calculated at 33.33% of the estimated taxable profit for the year. For Germany, a rate of approx. 30% was applied to the taxable income. Income generated outside France and Germany is taxed at rates applicable in the corresponding country.

Considering the French and German average tax rates and the impact of other tax legislations, the expected tax rate for the Sartorius Stedim Biotech Group is roughly 26% (29% in 2017). The reduction of 3% in the expected tax rate is mainly in connection with the reduced corporate tax rate in the US. The following table describes the difference between the expected tax expense and the income tax expenses reported for the particular financial year.

	2018 12 months € in K	2017 12 months € in K
Expected tax rate	26%	29%
Expected tax expense	-73,983	-63,978
Differences from the Group average income tax rate	9,771	8,638
Permanent differences	-11,172	-1,551
Tax-free income and other tax exemptions	1,222	1,262
Taxes for previous years	604	-853
Withholding and similar taxes	-404	-596
Other	-628	228
Total	-74,590	-56,849
Effective tax rate	-26.2%	-25.8%

14. Profit or Loss Statement by Nature

	2018 12 months € in K	2017 12 months € in K
Sales revenue	1,212,152	1,081,033
Purchases consumed	-237,889	-235,762
Cost of purchased services	-54,079	-48,353
Personnel costs	-340,853	-307,007
Amortization and depreciation	-60,914	-50,678
Other operating costs	-218,183	-217,531
Subtotal	-911,918	-859,331
Operating profit (EBIT)	300,234	221,702
Financial income expenses	-15,685	-1,089
Income tax	-74,590	-56,849
Non-controlling interest	-1,907	-2,676
Net profit after non-controlling interest	208,052	161,088

15. Earnings per Share

According to IAS 33, the earnings per share must be determined as follows: The basic earnings per share (basic EPS) are calculated on the basis of the weighted average number of ordinary shares during the period.

	2018	2017
Net profit after tax (€ in K)	209,959	163,763
Group net profit after tax (€ in K)	208,052	161,088
Earnings per share (€)	2.26	1.75
Diluted earnings per share (€)	2.26	1.75
Number of shares (statutory level)	92,180,190	92,180,190
Treasury shares	-3,416	0
Weighted average number of shares used in earnings per share calculation	92,176,774	92,180,190
Weighted average number of shares used in diluted earnings per share calculation	92,176,774	92,180,190

Notes to the Individual Balance Sheet Items

16. Goodwill and Other Intangible Assets

Goodwill

	Goodwill € in K
Gross book values at Jan. 1, 2017	344,777
Currency translation	-4,791
Business combinations	46,059
Gross book values at Dec. 31, 2017	386,045
Impairment losses at Jan. 1, 2017	0
Currency translation	0
Impairment losses	0
Impairment losses at Dec. 31, 2017	0
Net book values at Dec. 31, 2017	386,045
	Goodwill € in K
Gross book values at Jan. 1, 2018	386,045
Currency translation	-1,350
Business combinations	0
Gross book values at Dec. 31, 2018	384,695
Impairment losses at Jan. 1, 2018	0
Currency translation	0
Impairment losses	0
Impairment losses at Dec. 31, 2018	0
Net book values at Dec. 31, 2018	384,695

The caption reported as goodwill in the amount of €384,695K is the capitalized difference in assets resulting from business combinations. According to IAS 36, goodwill acquired in a business combination may not be amortized, but rather, must be tested annually for impairment and as soon as there is any indication of asset impairment. The increase recorded in 2017 concerns the acquisition of Umetrics.

For the purpose of impairment testing, goodwill must be allocated to each of the acquirer's cash-generating units (CGUs) that are expected to benefit from the synergies of the combination. The CGU represents the lowest level within the entity at which goodwill is monitored for internal management purposes and may not be larger than a segment. Sartorius Stedim Biotech

Group follows the strategy to be a total solution provider for its customers. Because of the various interdependencies within the business, the lowest level at which goodwill is monitored is that of the Biopharm segment. Therefore, the acquired goodwill is allocated to this group of CGU's.

As in 2017, the impairment test conducted for 2018 measures the recoverable amount on the basis of the value in use of the particular cash-generating unit (Biopharm segment). The cash flow forecasts consider previous experiences and are generally based on Group management's forecasts for a period of four years. The calculations were based on a terminal year growth rate of 2.5% for the years after 2022. This rate is derived from market expectations, which forecast significant growth rates for the targeted biopharmaceutical market. The major growth driver for the Sartorius Stedim Biotech Group will be the aging and increase in population and the improved access to drugs in the emerging markets as well as the ongoing paradigm shift from reusable products to single-use products utilized in bio manufacturing by the biopharmaceutical industry.

The discount rates applied correspond to the weighted capital cost rates and were recognized as follows:

	2018		2017	
	Before tax	After tax	Before tax	After tax
Biopharm segment	8.9%	7.2%	8.5%	6.8%

In 2018, our impairment test did not result in recognition of impairment losses. In this context, various sensitivity analyses based on realistic variations of the assumptions disclosed above did not result in an impairment either. The following variations would theoretically represent the "break-even point":

	2018	2017
Discount rates	22.9%	20.4%
Terminal growth rate	-29.6%	-28.4%
Cash flows	-78.5%	-80.2%

Intangible Assets

	Patents, licenses and similar rights € in K	Brand name € in K	Customer relationships € in K	Capitalized development costs € in K	Payments on account € in K	Total € in K
Gross book values at Jan. 1, 2017	74,664	11,061	121,351	68,054	36	275,165
Currency translation	-2,817	-47	-1,006	-387	0	-4,257
Business combinations	11,614	892	14,486	0	0	26,992
Acquisitions	5,767	0	469	20,788	0	27,024
Disposals	-34	0	0	-2,685	0	-2,720
Transfers	-143	0	0	155	0	12
Gross book values at Dec. 31, 2017	89,051	11,906	135,300	85,924	36	322,216
Amortization and impairment losses at Jan. 1, 2017	-27,813	-68	-67,710	-35,556	0	-131,147
Currency translation	1,298	5	410	139	0	1,853
Amortization and impairment losses	-7,306	-98	-10,072	-5,176	0	-22,652
Disposals	34	0	0	2,607	0	2,641
Transfers	-5	0	0	0	0	-5
Amortization and impairment losses at Dec. 31, 2017	-33,792	-160	-77,372	-37,986	0	-149,310
Net book values at Dec. 31, 2017	55,259	11,746	57,928	47,937	36	172,906

	Patents, licenses and similar rights € in K	Brand name € in K	Customer relationships € in K	Capitalized development costs € in K	Payments on account € in K	Total € in K
Gross book values at Jan. 1, 2018	89,051	11,906	135,300	85,924	36	322,216
Currency translation	246	-32	172	-52	1	335
Business combinations	0	0	0	0	0	0
Acquisitions	20,443	0	0	22,839	19	43,301
Disposals	-11,010	0	-11,804	-1,089	0	-23,902
Transfers	363	0	0	0	0	363
Gross book values at Dec. 31, 2018	99,094	11,874	123,669	107,622	55	342,313
Amortization and impairment losses at Jan. 1, 2018	-33,792	-160	-77,372	-37,986	0	-149,310
Currency translation	-444	1	-41	28	0	-455
Amortization and impairment losses	-11,519	-152	-13,991	-4,341	0	-30,004
Disposals	8,183	0	5,680	1,089	0	14,952
Transfers	-3	0	0	0	0	-3
Amortization and impairment losses at Dec. 31, 2018	-37,576	-311	-85,724	-41,210	0	-164,821
Net book values at Dec. 31, 2018	61,518	11,563	37,944	66,412	55	177,492

Intangible assets acquired are recorded at cost less the accumulated, regular amortization that is calculated according to the straight-line method and any impairment loss. The useful life of an intangible asset is the period during which the Group expects to use the asset.

Amortization of intangible assets is based on the following estimated useful lives:

Software	2 to 5 years
Capitalized R&D expenses	4 to 6 years
Customer relations and technologies	5 to 15 years
Brand name	5 years to indefinite

Costs incurred within the scope of the development of new products and methods were capitalized as internally generated intangible assets if the following criteria were met:

- The technical feasibility of completing the intangible assets so that it will be available for use or sale;
- The intention to complete the intangible asset and use or sell it;
- The ability to use or sell the intangible asset;
- The demonstration of how the intangible asset will generate probable future economic benefits;
- The availability of adequate technical, financial and other resources to complete the development and to use or sell the intangible asset;
- The ability to measure reliably the expenditure attributable to the intangible asset during its development.

The capitalized development costs essentially cover the costs that were allocated to the staff involved in R&D, raw materials and supplies, outside services and directly attributable overhead. Intangible assets generated internally are amortized on a straight line basis over their useful lives, which generally do not exceed six years.

If an internally generated intangible asset may not be recognized, the development costs are expensed in the period in which they are incurred. Costs for research activities are reported as expenses in the period in which they are incurred.

The Stedim brand name acquired in 2007 is considered to have an indefinite useful life and is therefore not amortized. There is no foreseeable limit to the period over which the brand name is expected to generate net cash inflows for the Group. The brand name is tested annually for impairment at the level of the "Biopharm segment" cash-generating unit (CGU).

In 2018, the development costs of €22,839K were recognized as assets (€20,788 K in 2017).

Amortization of intangible assets is allocated to the corresponding functions in the statement of profit or loss. For capitalized development costs, amortization is disclosed in the line "cost of sales".

As a consequence of the modified agreement with Lonza in the cell culture media business, the related intangible assets (technology and customer relationships) have been recognized as disposals (see note 11).

In 2018, impairments were recognized in the amount of €5.8million. The impairments relate mainly to technologies and customer relationships.

17. Property, Plant and Equipment

	Land, buildings and improvements € in K	Technical machinery and equipment € in K	Factory and office equipment and other equipment € in K	Payments on account and construction in progress € in K	Total € in K
Gross book values at Jan. 1, 2017	171,308	152,063	81,784	36,704	441,860
Currency translation	-4,127	-2,165	-2,873	-4,742	-13,907
Business combinations	0	0	141	0	141
Acquisitions	2,317	7,438	10,995	88,958	109,708
Disposals	-454	-2,683	-3,153	-1,797	-8,087
Transfers	7,494	7,209	1,710	-15,944	469
Gross book values at Dec. 31, 2017	176,538	161,862	88,603	103,179	530,183
Depreciation at Jan. 1, 2017	-52,038	-78,460	-49,899	1	-180,396
Currency translation	1,033	764	2,246	0	4,042
Depreciation	-6,947	-12,687	-8,393	0	-28,026
Disposals	436	2,553	2,978	0	5,967
Transfers	-405	190	-260	0	-475
Depreciation at Dec. 31, 2017	-57,921	-87,641	-53,327	1	-198,888
Net book values at Dec. 31, 2017	118,617	74,222	35,275	103,181	331,295

	Land, buildings and improvements € in K	Technical machinery and equipment € in K	Factory and office equipment and other equipment € in K	Payments on account and construction in progress € in K	Total € in K
Gross book values at Jan. 1, 2018	176,538	161,862	88,603	103,179	530,183
Currency translation	386	-302	445	3,709	4,238
Business combinations	0	0	0	0	0
Acquisitions	20,018	11,247	9,137	93,242	133,644
Disposals	-3,643	-7,316	-9,959	-299	-21,218
Transfers	6,035	11,899	852	-19,113	-326
Gross book values at Dec. 31, 2018	199,335	177,390	89,078	180,718	646,521
Depreciation at Jan. 1, 2018	-57,921	-87,641	-53,327	1	-198,888
Currency translation	35	-242	-133	0	-341
Depreciation	-8,490	-13,443	-8,980	0	-30,913
Disposals	3,320	6,712	9,257	0	19,290
Transfers	-11	72	250	0	311
Depreciation at Dec. 31, 2018	-63,067	-94,542	-52,933	1	-210,541
Net book values at Dec. 31, 2018	136,268	82,848	36,146	180,719	435,980

The "Property, plant and equipment" caption is recorded at cost, and related assets are depreciated over their estimated useful life using the straight line method. Tangible assets are subject to impairment tests whenever there are indicators of impairment.

Depreciation of fixed assets is based on the following periods of useful life:

Buildings	15 to 50 years
Machinery	5 to 15 years
Factory and office equipment	3 to 13 years

Depreciation is included in the statement of profit or loss according to use of the assets in the cost of sales, selling and distribution costs, research and development costs, general administrative expenses and other operating expenses.

Borrowing costs are expensed as incurred unless they are directly attributable to the acquisition, construction or production of a qualifying asset and are therefore part of the cost of that asset. A qualifying asset is defined as an asset that takes a substantial period of time (six to twelve months) to get ready for its intended use.

Grants related to assets are deducted from the cost of the related asset.

In 2018, as for fiscal 2017, there were no significant impairment losses to be recognized with regard to property, plant and equipment.

18. Leases

A lease is an arrangement whereby the lessor conveys to the lessee in return for a payment or series of payments the right to use an asset for an agreed period of time. According to IAS 17 a lease is classified as either an operating or a finance lease. A finance lease is a lease that transfers substantially all the risks and rewards incidental to ownership of an asset. All other leases are designated as operating leases.

When the Group is a lessee in a finance lease, the amount equal to the fair value of the leased property, or if lower, the present value of the minimum lease payments is recognized as an asset on the balance sheet and simultaneously recognized as a financial liability. The minimum lease payments essentially consist of the finance charge and the reduction of the outstanding liability, which are measured according to the effective interest method. A leased asset is depreciated on a straight-line basis over the period of its expected useful life or over the shorter lease term.

For an operating lease, the lease instalments to be paid by the lessee are recognized as expenses over the lease term and the lease payments received by the lessor are recognized as income, respectively. The leased asset continues to be recognized on the lessor's balance sheet as property, plant and equipment.

Capitalized property, plant and equipment for assets held under finance leases amount to €14,599 K (2017: €15,672). The cost of acquisition of these assets totals €19,631 K (2017: €19,682 K).

In 2018 rental payments amounting to €14.0 million (2017: €11.3 million) were made for assets leased under operating leases.

For the new lease accounting rules according to IFRS 16 (applicable from 2019 on) please refer to note 2.

The Group's financial obligations from operating lease commitments (future minimum lease payments under non-cancellable leases), are as follows:

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Operate leases		
- due within one year	11,434	8,682
- due within 2 to 5 years	22,145	13,900
- due thereafter	8,813	4,817

19. Deferred Tax

	Dec. 31, 2018 € in K	Deferred Tax Assets		Deferred Tax Liabilities	
		Dec. 31, 2017 € in K	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K	Dec. 31, 2018 € in K
Intangible assets	1,066	783	37,746	36,630	
Tangible assets	0	0	5,865	6,553	
Inventory	10,236	9,143	0	0	
Receivables and other current assets	230	189	640	5,105	
Provisions	6,384	6,649	0	0	
Liabilities	5,391	2,181	1,866	3,140	
Gross amount	23,307	18,945	46,117	51,428	
Carry forward of taxable losses	0	2,334	0	0	
Tax on undistributed earnings of subsidiaries	0	0	1,850	1,848	
Offset	-8,817	-10,289	-8,817	-10,289	
Net amount	14,490	10,989	39,150	42,987	
Change	3,501	235	3,837	-14,207	
thereof recognized in profit or loss	3,450	292	291	-968	

Deferred tax assets or liabilities are determined based on temporary differences between the carrying amounts and the tax basis of assets and liabilities (except in special cases provided by IAS 12) including loss carry forwards and tax credits. Measurement is based on the tax rates expected to be effective in the period in which an asset is realized or a liability is settled.

For this purpose, the tax rates and tax rules are used which have been enacted or substantively enacted at the reporting date. Deferred tax assets are recognized for deductible temporary differences and tax losses and unused tax credits only to the extent that it is probable that the Group will have future taxable income against which they can be charged.

Deferred Tax Assets

On the balance sheet date, the Group had unused tax loss amounts carried forward of €11.6 million to be deducted from future taxable profits (€18.2 million in 2017). A deferred tax amount was reported on approx. €0.0 million of these losses (€9.8 million in 2017). Concerning the remaining losses to be carried forward, no deferred tax amounts were recognized because of the lack of visibility of future taxable profits.

Deferred tax assets in the amount of €0.0 million (€2.3 million) relate to companies that reported losses in this year under review or in the earlier reporting year.

Deferred Tax Liabilities

The deferred tax liabilities in connection with intangible assets refer to assets acquired in business combinations and consequently are mainly linked to customer relationships and technologies.

The Group did not record deferred tax liabilities on approx. €658 million (€480 million) in cumulative undistributed earnings of subsidiaries because these earnings are intended to be reinvested in these operations. When the dividends are paid out, an amount of 5% of the dividends will be taxed under the French and German taxation rules and, if applicable, with withholding tax. Furthermore, additional income tax consequences could arise in the case of an intermediate holding company.

In fiscal 2018, as in the previous years, the tax effect from hedging instruments, and the deferred tax assets from the recognition of the remeasurements of defined benefit liabilities (assets) were recognized in other comprehensive income. The income taxes recognized in other comprehensive income are disclosed as follows in the table:

€ in K	2018	2017
Cash flow hedges	3,463	-7,889
Remeasurements of the net defined benefit obligations	-160	481
Total	3,303	-7,408

20. Inventories

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Raw materials and supplies	68,318	60,130
Work in progress	71,985	53,205
Finished goods and merchandise	108,520	68,008
Payments on account	3,179	4,261
Total	252,002	185,604

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Gross amount inventories	269,013	199,824
Write-downs	-17,011	-14,221
Net Amount Inventories	252,002	185,604

Raw materials and supplies, including merchandise, are reported under "Inventories" at average cost. In principle, finished goods and work in progress are reported at cost of conversion. This cost includes direct costs, which can be allocated to these materials, and the appropriate portion of production and materials handling overhead, general administrative expenses and fixed assets at normal depreciation and/or amortization rates, based on the normal production capacity, provided that these expenses are caused by production.

Inventories must be valued at the lower amount of cost and the net realizable value. The net realizable value represents the estimated selling price in the ordinary course of business less the estimated costs of completion and the estimated costs necessary for marketing, sales and distribution. Where inventory risks exist, such as the risk of reduced shelf life as a result of storage periods or limited usability, inventories are marked down accordingly.

21. Issued Capital

Sartorius Stedim Biotech S.A. share capital consists of 92,180,190 shares with a par value of 0.20 €.

As of December 31, 2017, and December 31, 2018, there were no dilutive instruments. Shares registered

in the name of the same owner for at least four years benefit from a double voting right.

	Dec. 31, 2018	Dec. 31, 2017
Number of shares at the beginning of the period	92,180,190	92,180,190
Number of shares at the end of the period	92,180,190	92,180,190
Nominal value per share (in €)	0.20	0.20
Issued capital amount (€ in K)	18,436	18,436

Dividends

The Board of Directors will submit a proposal to the Annual General Shareholders' Meeting for payment of a dividend for the year ended December 31, 2018, as follows: payment of a net dividend of €0.57 per share (2017: €0.46), i.e., a total disbursement of 52,540,761.00€ (excluding treasury shares; 2017: 42,402,887.00€).

22. Non-Controlling Interest

The non-controlling interest recognized in the statement of financial position amounting to €8,476K relate to the subsidiary Sartorius Korea Biotech. The Group's interest in Sartorius Korea Biotech is 69%, the remaining 31% are subject to an exercisable call option.

The purchase price for this non-controlling interest is variable and depends on the future performance of the entity.

Key Figures

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Sartorius Korea Biotech Co. Ltd.		
Sales revenue	63,111	61,477
Net result	6,150	8,631
Total assets	36,225	34,399
Attributed profit or loss	1,907	2,676

There are no significant restrictions on the Group's ability to access or use the assets or settle the liabilities of the mentioned entity.

23. Pension and Employee Benefits Provisions

Pension Obligations

Pension provisions and similar obligations are recognized in the consolidated financial statements of Sartorius Stedim Biotech Group in accordance with actuarial principles. IAS 19, Employee Benefits, stipulates the Projected Unit Credit Method as the method of measurement. In addition to known pensions and life expectancies, this expected cash value method takes into account future salary and pension increases.

All remeasurements of the net defined benefit liability are recognized in other comprehensive income (pension reserves) in accordance with the standard IAS 19.

Defined Contribution Plans

Most of the Sartorius Stedim Biotech Group companies make payments under defined contributions plans, primarily related to government-run pension plans. In 2018, the total expense recognized for the defined contribution plans amounted to €20,807 K (2017: €19,527 K).

Defined Benefit Plans

The remeasurements of defined benefit liabilities (asset) are shown in other comprehensive income according to the standard IAS 19. The actuarial losses, which were transferred to the pension reserves, essentially resulted from a change in the discount rate and totaled €469 K (€-667 K in 2017).

An amount of €24,441 K relates in particular to pension provisions for retirement pension plans in Germany. These provisions totaled €25,068 K in 2017 and primarily relate to direct commitments under defined benefit pension plans. Under these commitments, the employees earn benefits for each year of service rendered to the company. The benefits earned depend on the salary level and the age of the respective employees. The pension benefits are generally not funded with assets.

The assumed discount rates reflect the interest rates payable on the reporting date for high-quality corporate bonds with matching maturities and denominated in the relevant currencies (mainly Euro). If such corporate bonds are not available with matching long-term maturities or are insufficiently

available, their matching interest rates are determined by extrapolation.

Measurement of the post-employment benefit obligations is based on the following actuarial assumptions:

For Germany:

in %	Dec. 31, 2018	Dec. 31, 2017
Discount rate	1.81	1.75
Future salary increases	3.00	3.00
Future pension increases	2.00	2.00

With regard to the assumptions for mortality and disability the tables "Richttafeln (RT) 2018 G" by Klaus Heubeck were applied.

For France:

in %	Dec. 31, 2018	Dec. 31, 2017
Discount rate	1.80	1.70
Future salary increases	2.00	2.50
Future pension increases	2.00	2.00

The amounts reported in the statement of profit or loss and other comprehensive income consist of the following:

	2018 € in K	2017 € in K
Current service cost	-1,978	-1,856
Past service cost	485	408
Net interest expenses	-531	-509
Components of defined benefit costs recognized in profit or loss	-2,023	-1,958
Return on plan assets (excl. interest)	56	57
Remeasurements	413	-723
Components of defined benefit costs recognized in other comprehensive income	469	-667
Total	-1,554	-2,624

In the statement of profit or loss, the current service cost is disclosed according to the assignment of employees to the respective functions.

The amount included in the consolidated statement of financial position arising from the Group's obligation in respect of defined benefit plans is as follows:

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Present value of the obligations	46,459	43,752
Fair value of the plan assets	-10,865	-8,306
Net Liability	35,595	35,446

The present value of the defined benefit obligation developed as follows:

	2018 € in K	2017 € in K
Present value of the obligations as of Jan. 1	43,752	40,841
Current service cost	1,978	1,856
Past service cost	-485	-408
Interest cost	649	585
Remeasurements	-548	745
Foreign currency translation differences	386	-725
Retirement benefits paid in the reporting year	-1,011	-852
Employee contributions	261	220
Contributions by plan participants	1,466	1,550
Other changes	14	-60
Present value of the obligations as of Dec. 31	46,459	43,752

The remeasurements of defined benefit liability (asset) can be allocated as follows:

	2018 € in K	2017 € in K
Experience adjustments	367	850
Changes in demographic assumptions	306	632
Changes in financial assumptions	-1,221	-737
Total	-548	745

Plan Assets

	2018 € in K	2017 € in K
Plan assets as of Jan. 1	8,306	6,622
Interest income	118	75
Return on plan assets (excl. interest)	56	57
Remeasurements	-136	22
Group contribution & payments	-887	-709
Foreign currency translation differences	279	-481
Employee contributions	261	220
Employer contributions	1,402	894
Contributions by plan participants	1,466	1,606
Other changes	0	0
Plan assets as of Dec. 31	10,865	8,306

Composition of Plan Assets

The plan assets do primarily refer to insurance contracts in Germany and Switzerland, there are no major equity or debt investments included. The subsidiary in South Korea has deposited an amount of €2.4 million (€1.8 million in 2017) to local banks as cash and cash equivalents.

Sensitivity Analysis

An increase|decrease of the actuarial assumptions would have the following impacts on the defined benefit obligations (a positive sign (+) means an increase of the obligation):

2017:

	€ in K	
Demographic assumptions		
Life expectancy	+1 year	-1 year
Effect	874	-870
Financial assumptions		
Discount rate	+100 bps	-100 bps
Effect	-6,105	7,848
Future salary increases	+50 bps	-50 bps
Effect	865	-803
Future pension increases	+25 bps	-25 bps
Effect	984	-937

2018:

€ in K		
Demographic assumptions		
Life expectancy	+1 year	-1 year
Effect	1,053	-1,052
Financial assumptions		
Discount rate	+100 bps	-100 bps
Effect	-6,389	8,171
Future salary increases	+50 bps	-50 bps
Effect	938	-861
Future pension increases	+25 bps	-25 bps
Effect	1,020	-972

The sensitivity analysis presented above might not be representative of the actual change in the defined benefit obligation as it is unlikely that the change in assumptions would occur in isolation of one another. Furthermore, the present value of the defined benefit obligation has been calculated using the same method that was applied in calculating the defined benefit obligation liability recognized in the statement of financial position (projected unit credit method).

Maturity Analysis

The undiscounted cash flows from defined benefits obligations can be allocated to maturities as follows:

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
<1 year	1,664	1,477
1-5 years	8,336	7,616
6-10 years	12,430	11,564
>10 years	73,605	58,288
Total	96,035	78,944

The weighted average duration of the defined benefit obligations is 16.6 years (2017: 17.6 years).

24. Other Provisions

A provision is recognized when a present legal or constructive obligation to third parties arising from past events has been incurred, an outflow of resources is probable and the amount of the obligation can be reasonably estimated. The amount recognized as a provision represents the best estimate of the obligation at the closing date.

Restructuring provisions are set up in connection with programs that materially change the scope of business performed by a segment or business unit or the manner in which business is conducted. In most cases, restructuring expenses include termination benefits and compensation payments due to the termination of agreements with suppliers and dealers, including leasing contracts. Restructuring provisions are recognized when the Group has a detailed formal plan that has either commenced implementation or been announced.

Other Non-current Provisions

	Payments to employees on early retirement plan € in K	Other € in K	Total € in K
Balance at Jan. 1, 2017	2,331	752	3,083
Currency translation	0	-11	-11
Consumption	-1,035	-16	-1,051
Additions	877	205	1,082
Balance at Dec. 31, 2017	2,173	930	3,103

	Payments to employees on early retirement plan € in K	Other € in K	Total € in K
Balance at Jan. 1, 2018	2,173	930	3,103
Currency translation	0	9	9
Consumption	-1,124	-41	-1,165
Additions	824	105	929
Balance at Dec. 31, 2018	1,873	1,004	2,877

The non-current provisions comprise mainly provisions for partial retirement and employee anniversary bonuses (included in the item "other"). These obligations arise mainly in German Group companies. The partial retirement plans allow employees to work part-time for 3 - 5 years before their actual retirement.

According to IAS 19 these obligations are treated as severance payments to be earned in future periods and

are therefore recognized in profit or loss over the respective period of service. Actuarial gains and losses, as well as past service costs, on these obligations are recognized as income or expense.

Non-current provisions are reported at their present value on the reporting date. The discount rate for employees on the early retirement plan is 0.1% (2017: 0.0%).

Current Provisions

During financial 2017 and 2018, the current provisions developed as follows:

	Warranties € in K	Other € in K	Total € in K
Balance at Jan. 1, 2017	4,914	4,367	9,281
Currency translation	-56	-73	-129
Consumption	-1,362	-1,095	-2,457
Release	-837	-1,539	-2,377
Additions	2,165	3,651	5,816
Other changes	0	-577	-577
Balance at Dec. 31, 2017	4,824	4,735	9,558

	Warranties € in K	Other € in K	Total € in K
Balance at Jan. 1, 2018	4,824	4,735	9,558
Currency translation	4	-14	-9
Consumption	-934	-1,045	-1,980
Release	-2,550	-649	-3,199
Additions	5,021	2,892	7,913
Other changes	0	0	0
Balance at Dec. 31, 2018	6,364	5,919	12,283

Warranty provisions contain expenses for replacement deliveries and repairs. Specific risks are recognized when the occurrence is more likely than not. General warranty risks are considered on the basis of experiences in the past. The other provisions contain mainly onerous contracts and uncertain liabilities to employees.

25. Other Financial Obligations | Contingent Assets and Liabilities

As was the case in the previous years there are no significant contingent liabilities or contingent assets to be reported.

Financial Instruments | Financial Risks

A financial instrument is any contract that gives rise to a financial asset of one entity and a financial liability or equity instrument of another entity. The following sections give an overview of the impact of financial instruments on the financial statements of the Sartorius Stedim Biotech Group and provide additional information on the balance sheet items, which contain financial instruments.

26. Initial Application of IFRS 9, Financial Instruments: Transition Effects and Accounting Policies

The initial date of application of IFRS 9, Financial Instruments is January 1, 2018. The Group applies IFRS 9 retrospectively without restating the prior-year figures. The comparative period is presented in line with previous rules (IAS 39). The aggregate effects from the transition to IFRS 9 are accounted for by way of an adjustment to the opening balance of retained earnings. The effect from the initial application of IFRS 9 results from the adjustment of the credit impairment approach for trade receivables and amounts to €0.3 million. As a result, this increased retained earnings from €612.7 million as of December 31, 2017 to €613.0 million as of January 1, 2018.

IFRS 9 replaces the existing guidelines in IAS 39, Financial Instruments: recognition and measurement. IFRS 9 includes revised guidelines for classification and measurement of financial instruments, including a new model of expected credit losses for the calculation of impairments of financial assets, as well as the new general guidelines for hedge accounting. This standard also adopts the guidance of IAS 39 for recognition and derecognition of financial instruments.

Under IFRS 9, the new classification and measurement approach for financial assets reflects both the entity's business model (held-to-collect, held-to-collect-and-sell, other) within the scope of which assets are held and the contractual cash flow characteristics (so-called "SPPI" criterion, solely payments of principal and interest). For each class of the Group's financial assets, the following table provides an overview of the categories according to IAS 39 and IFRS 9, as well as the respective carrying amounts at the date of initial application of IFRS 9.

Carrying amounts of financial assets at the date of initial application of IFRS 9	Category acc. to IAS 39	Carrying amount acc. to IAS 39 Dec. 31, 2017 € in mn	Reclassification € in mn	Remeasure-ment € in mn	Carrying amount acc. to IFRS 9 Jan. 1, 2018 € in mn	Category acc. to IFRS 9
Cash and cash equivalents	Loans and receivables	32.6			32.6	Measured at amortized cost
Receivables and other assets	Loans and receivables	12.8			12.8	Measured at amortized cost
Derivative financial instruments in hedge relationships	n/a	7.2			7.2	n/a
Other financial assets (current)		20.0			20.0	
Trade receivables	Loans and receivables	204.6		0.3	205.0	Measured at amortized cost
Amounts due from customers for contract work	n/a	7.0		0.0	7.0	n/a
Trade receivables		211.6		0.3	211.9	
Financial assets	Loans and receivables	2.7	0.7		3.4	Measured at amortized cost
Financial assets	Available for sale	1.6	-1.0		0.6	Debt instruments at fair value through profit or loss
Financial assets	Available for sale	0.1			0.1	Equity instruments at fair value through profit or loss
Financial assets (investments in non-consolidated subsidiaries)	n/a	0.1			0.1	n/a (investments in non-consolidated subsidiaries)
Derivative financial instruments in hedge relationships	n/a	7.3			7.3	n/a
Financial assets (non-current)		11.8	-0.4		11.5	

There were no effects on the Group's financial liabilities. The few reclassifications of financial assets were determined on the basis of individual assessments of the financial instruments, especially with regard to the contractual cash flow characteristics. For equity instruments that existed as of the date of initial application of the Standard and that were not held for trading, the Group decided to recognize future changes in the fair value of those instruments in profit or loss. This choice is generally to be made on an instrument-by-instrument basis upon initial recognition of the instrument. Reclassification at the date of initial application did not lead to measurement effects. The measurement effects presented result from the adjustment of the impairment approach for trade receivables. There were no reclassifications of financial instruments during the reporting period.

With regard to the impairment of financial assets, IFRS 9 replaces the incurred-loss model by the expected-loss model. Financial assets are generally regarded as credit-impaired when there are objective

indications that cast doubt about the full collection of the cash flows of the respective financial assets. With regard to the financial assets of the Group, the simplified approach which is applied to trade receivables is especially of relevance.

At the date of initial application of IFRS 9, the allowance for expected credit losses amounted to €0.1 million. This amount includes the expected credit loss in relation to contract assets according to IFRS 15 that do not contain a significant financing component. In the course of the transition to IFRS 9, the Group also analyzed and revised its approach for the recognition of incurred losses, given its low level of historical losses. At the date of initial application of IFRS 9, the total effect from the adjustments to the impairment approach for trade receivables amounted to €0.3 million and increased the Group's retained earnings as of January 1, 2018.

The following table illustrates the effects from the adjusted impairment model for trade receivables:

Valuation allowances at the date of initial application of IFRS 9	Category acc. to IAS 39	Loss allowances acc. to IAS 39 Dec. 31, 2017 € in mn	Remeasurement due to implementation of IFRS 9 € in mn	Loss allowances acc. to IFRS 9 Jan. 1, 2018 € in mn	Category acc. to IFRS 9
Trade receivables	Loans and receivables	-2.4	0.3	-2.1	Measured at amortized cost
Amounts due from customers for contract work	n/a	0.0	0.0	0.0	n/a
Trade receivables		-2.4	0.3	-2.1	

Besides trade receivables, cash and cash equivalents are the most material financial assets on the Group's statement of financial position as of the date of initial application of IFRS 9 and as of the reporting date, December 31, 2018. No impairment is recognized for these financial assets due to materiality considerations.

For the remaining financial assets that are measured at amortized cost, no impairment is recognized as of the date of initial application of IFRS 9 and as of December 31, 2018, for the 12-month expected credit losses, given the Group's immaterial historical losses.

Derivatives are measured at fair value determined according to the mark-to-market method in which recognized mathematical methods are used. The fair values are based on the market data available at the time the value of these derivatives is calculated and reflect the estimates of the market conditions at the end of the year. Those instruments that are not designated as hedging instruments and for which no hedge accounting is applied, are categorized as held for trading. Changes of the fair values of derivative financial instruments are either recognized in profit or loss or, in case of hedging relationships, in other comprehensive income.

The Group applies the new hedge accounting rules of IFRS 9 prospectively. The Group uses forward transactions to hedge cash flow risks that result from changes in foreign exchange rates in relation to sales of products and the procurement of materials, and designates only the spot element of the hedging instrument. No transition effects resulted from the application of the new hedge accounting requirements.

Financial Assets

Financial assets of the Group mainly include cash and cash equivalents, trade and loan receivables and derivative financial instruments with a positive fair value.

27. Cash and Cash Equivalents

The Group considers all highly liquid investments with less than three months maturity from the date of acquisition to be cash equivalents. This mainly includes checks, cash on hand and deposits in banks. Cash and cash equivalents are measured at cost. For purposes of the consolidated cash flow statement, cash and cash equivalents include cash and cash equivalents as defined above. As of December 31, 2018, cash and cash equivalents amounted to €23,975 K (2017: €32,552 K).

28. Current Trade Receivables | Other Receivables

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Trade receivables from third parties	204,931	192,866
Amounts due from customers for contract work*	3,902	6,967
Receivables from subsidiaries of the Sartorius AG Group	11,399	11,782
Trade receivables	220,231	211,616

*Contract assets according to IFRS 15.

The book values of trade receivables and other receivables are representative of their fair value considering the maturity date and the credit risks. The contract assets are recognized in connection with customer-specific construction contracts that meet the requirements for revenue recognition over time according to IFRS 15 (see section 9). The "Receivables from subsidiaries of the Sartorius AG Group" item refers to other companies of the Sartorius Group (please refer to section 41). Impairment losses on trade and other receivables are recognized using separate allowance accounts. For details on the determination of the impairment allowances see section 38.

30. Loans and Borrowings

	Balance at Dec. 31, 2018 € in K	of which current Dec. 31, 2018 € in K	Balance at Dec. 31, 2017 € in K	of which current Dec. 31, 2017 € in K
Liabilities to banks	38,278	35,153	29,367	23,117
Loans from Sartorius AG	94,501	54,501	109,481	69,481
Other loans from Sartorius Group companies	163	163	3,335	3,335
Total loans and borrowings	132,942	89,817	142,183	95,933

Sartorius Stedim Biotech Group has signed a loan agreement with its parent company Sartorius AG which secures the financing of the Sartorius Stedim Biotech Group over the long term. The credit volume of this agreement is approx. 300 million euros and the interest rate is variable with a credit margin based on arms'-length principles.

29. Other Financial Assets

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Derivative financial instruments	2,824	7,207
Other financial assets	19,212	12,842
Current financial assets	22,036	20,049

The amount shown as derivative financial instruments represents the fair value of foreign currency hedging instruments, mainly forward contracts (for details refer to section 35).

Other financial assets are measured at amortized cost using the effective interest method less any impairment losses. The item "Other financial assets" includes loan receivables to other entities of the Sartorius AG Group in the amount of €2,588 K (2017: €1,542 K).

Financial Liabilities

Financial liabilities of the Group mainly comprise loans borrowed from Sartorius AG, trade payables, finance lease liabilities and derivative financial instruments with a negative fair value. The financial liabilities other than derivative financial instruments are measured at amortized cost.

The non-current loans and borrowings do not include the liabilities in connection with acquisitions which are presented in the caption "other non-current liabilities".

31. Other Non-current Liabilities

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Other liabilities	8,889	30,392
Total	8,889	30,392

The other non-current liabilities include the liability for phantom units that was incurred in connection with the acquisition of the non-controlling interests in the company AllPure Technologies, LLC amounting to €8,739K. The purchase price depends on the future sales and is due 2022 at the latest. As a result of the positive development of the AllPure business, the expected payments are determined by considering future revenue at an annual growth rate of about 20% on average. Due to this change in expectations regarding revenue growth, the liability has been increased by about €3 million. The effect is recognized in profit or loss. An increase (decrease) of the sales revenue by 10% in each of the following years would lead to an increase (decrease) of the liability by €0.9 million (€0.9 million).

In the previous reporting period, the other non-current liabilities included the non-current part of the liability for the remaining purchase price for the cell culture media business of the company Lonza. In connection with the change in cooperation between the Group and Lonza, the Group was released from the obligation to pay the remaining purchase price in the reporting period. Therefore, the related liability was released in 2018 and the effects on profit or loss are recognized in other operating income (see note 11).

Trade Payables

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Payments received on account of orders*	51,000	40,749
Trade payables to third parties	86,187	69,900
Payables to participations	6	2
Payables to subsidiaries of the Sartorius AG Group	17,375	7,309
Total	154,568	117,959

*Contract liabilities according to IFRS 15.

32. Other Current Financial Liabilities

	Dec. 31, 2018 € in K	Dec. 31, 2017 € in K
Derivative financial instruments	1,346	1,928
Other liabilities	12,966	33,036
Total	14,312	34,964

The derivative financial instruments refer to the fair values of foreign currency hedging transactions such as forward contracts (mainly related to the US\$).

The reduction in other financial liabilities is linked to the release of the liability for the acquisition of the cell culture media business (see note 11).

33. Carrying Amounts and Fair Values of Financial Instruments according to Categories

The following tables show the carrying amounts and fair values of financial assets and liabilities by category of financial instrument as of December 31, 2018, under IFRS 9 and as of December 31, 2017, under IAS 39:

	Category acc. to IFRS 9	Carrying amount Dec. 31, 2018 € in K	Fair value Dec. 31, 2018 € in K
Investments in non-consolidated subsidiaries	n/a	109	109
Financial assets	Equity instruments at fair value through profit or loss	50	50
Financial assets	Debt instruments at fair value through profit or loss	671	671
Financial assets	Measured at amortized cost	4,814	4,814
Financial assets (non-current)		5,644	5,644
Amounts due from customers for contract work	n/a	3,902	3,902
Trade receivables	Measured at amortized cost	216,330	216,330
Trade receivables		220,231	220,231
Receivables and other assets	Measured at amortized cost	19,212	19,212
Derivative financial instruments designated as hedging instruments*	n/a	2,824	2,824
Other financial assets (current)		22,036	22,036
Cash and cash equivalents	Measured at amortized cost	23,975	23,975
Loans and borrowings	Financial liabilities at cost	132,942	133,175
Trade payables	Financial liabilities at cost	103,568	103,568
Trade payables payments received for orders	n/a	51,000	51,000
Trade payables		154,568	154,568
Derivative financial instruments	Held for trading	1,346	1,346
Other financial liabilities	Financial liabilities at cost	21,855	21,326
Other financial liabilities		23,201	22,672

*The amounts include the non-designated part of the contracts.

	Category acc. to IAS 39	Carrying amount Dec. 31, 2017 € in K	Fair value Dec. 31, 2017 € in K
Financial Assets	Available for sale	1,813	1,813
Financial Assets	Loans and receivables	2,734	2,734
Derivative financial instruments in hedging relationship	Hedging Instruments*	7,260	7,260
Non-current financial assets		11,808	11,808
Trade receivables	Loans and receivables	211,616	211,616
Financial Assets	Loans and receivables	12,842	12,842
Derivative financial instruments in hedging relationship	Hedging Instruments*	7,207	7,207
Other financial assets		20,049	20,049
Cash and cash equivalents	Loans and receivables	32,552	32,552
Loans and borrowings	Financial liabilities at cost	142,183	142,671
Finance lease liabilities	IAS 17	17,487	28,086
Trade payables	Financial liabilities at cost	77,211	77,211
Trade payables	n/a	40,749	40,749
Trade payables		117,959	117,959
Derivative financial instruments	Held for trading	552	552
Derivative financial instruments in hedging relationship	Hedging Instruments*	1,376	1,376
Other financial liabilities	Financial liabilities at cost	63,428	67,719
Other financial liabilities		65,356	69,647

*The amounts include the non-designated part of the contracts.

The fair values of the financial instruments were determined on the basis of the market information available on the reporting date and are to be allocated to one of the three levels of the fair value hierarchy in accordance with IFRS 13.

Level 1 financial instruments are measured on the basis of prices quoted on active markets for identical assets and liabilities. In Level 2, financial instruments are measured on the basis of input factors which are derivable from observable market data or on the basis of market prices for similar instruments. Level 3 financial instruments are measured on the basis of input factors that cannot be derived from observable market data.

The financial instruments to be recognized at fair value on the reporting date are mainly derivatives in the form of forward contracts. They were measured on the basis of their quoted exchange rates and market yield curves (Level 2).

The fair values to be disclosed for financial liabilities recognized at amortized cost, especially liabilities to banks and finance leases, were measured on the basis of the market interest rate, taking the current indicative credit spreads into account (Level 2).

The fair values of the remaining financial assets and liabilities to be disclosed approximate the carrying amounts on account of their predominantly short-term maturity. The maximum credit loss risk is reflected by the carrying amounts of the financial assets recognized in the Statement of Financial Position.

The Group recognizes transfers between the levels of the fair value hierarchies at the end of the reporting period during which a change has occurred. In the current reporting period, there were no transfers between the levels.

34. Net Gains and Losses from Financial Instruments

The net gains and losses of the various categories of financial instruments are presented in the following table:

	2018 12 months € in K
Categories according to IFRS 9	
Financial assets measured at amortized cost	2,771
Financial assets and liabilities measured at fair value through profit or loss	552
Financial liabilities measured at amortized cost	-12,733

	2017 12 months € in K
Categories according to IAS 39	
Loans and receivables	-4,121
Financial assets and liabilities held for trading	2,358
Financial liabilities at cost	1,606

The net result from financial assets measured at amortized cost mainly includes the effects of currency translation and changes in allowances.

The net result from financial assets and liabilities held for trading predominantly comprises changes in the fair value of derivative financial instruments as well as interest income and interest expenses for these financial instruments.

The net result from liabilities measured at amortized cost mainly comprises the effects of foreign currency translation.

Total interest income and expenses for financial assets and liabilities that are not measured at fair value through profit or loss were as follows:

	2018 12 months € in K	2017 12 months € in K
Interest income	463	282
Interest expenses	-4,563	-5,454

Capital and Financial Risk Management

Capital Risk Management

In the Sartorius Stedim Biotech Group, capital is managed in order to maximize earnings of those participating in the company by optimizing the ratio of equity to liabilities. Furthermore, we ensure that all Group companies operate under the premise of the going-concern principle.

The financial liabilities detailed above are regarded as managed capital and, furthermore, so are the cash and cash equivalents as well as equity capital.

Goals of Financial Risk Management

The Treasury Department of the Sartorius Stedim Biotech Group is centrally focused in Sartorius Corporate Administration GmbH, a subsidiary of Sartorius AG. This centralized Treasury Department performs services for all companies of the Sartorius Group, including the Sartorius Stedim Biotech Group, and coordinates access to national and international financial markets. In addition, the Treasury Department monitors and controls financial risks by internal risk reporting, which analyzes risks according to their degree and scope. Essentially, these risks entail currency, interest rate and liquidity risks as well as credit risks.

The Sartorius Stedim Biotech Group strives to minimize the impact of currency and interest rate risks using appropriate derivative financial instruments. Hedging transactions and their controlling are carried out by different staff members. Moreover, the Group's Internal Auditing Department regularly monitors the use of such financial instruments. Trading with derivative financial instruments is done for hedging purposes only.

35. Management of Exchange Rate Risks and Hedge Accounting

The Group is exposed to currency risks as more than one third of sales revenue is generated in U.S. dollars or currencies linked to the U.S. dollar and, to a lesser extent, in other foreign currencies. At the same time, the Group is able to compensate the major part of the revenues denominated in foreign currencies with costs incurred in the same currencies due its global production network. The share of revenues generated in foreign currencies that exceeds such costs, the so-called net currency exposure, is hedged with derivative

financial instruments to a large extent (generally 70% to 80%). The Group generally follows a rolling hedging strategy of up to 12 months in advance. Also, the hedging measures are reviewed at regular intervals in order to adapt them to currency fluctuations.

For currency hedging, forward contracts are used. Forward contracts secure the right, and simultaneously create the obligation, to sell an established foreign currency amount on the exercise date at a specific exchange rate against the euro, independently of the exchange rate actually valid on this date. The profit or loss resulting from the difference between the current and the previously established exchange rate is

recognized as income or expense in the statement of profit or loss.

At the balance sheet date, forward contracts have been carried out in an amount of \$137 million (2017: \$190 million) to hedge against the risk of fluctuation in the EUR|USD exchange rate. This amount covers roughly half of the expected net exposure for the U.S. dollar within the period of 12 months. Furthermore, other foreign currencies have been hedged in smaller volumes.

The following table shows the forward transactions as of the balance sheet date:

Dec. 31, 2017	Currency	Volume	Maturity	Fair value € in K
Forward contract	USD	110,000	2018	6,988
	USD	80,000	2019	7,260
	USD	190,000		14,248
Forward contract	JPY	700,000	2018	72
	JPY	700,000		72
Forward contract	CHF	24,500	2018	-1,375
	CHF	24,500		-1,375
Forward contract	CAD	8,800	2018	148
	CAD	8,800		148
Dec. 31, 2018	Currency	Volume	Maturity	Fair value € in K
Forward contract	USD	50,000	2019	-53
	USD	50,000	2019	296
	USD	21,000	2019	712
	USD	16,000	2019	644
	USD	137,000		1,599
Forward contract	JPY	400,000	2019	-138
	JPY	200,000	2019	-67
	JPY	200,000	2019	-51
	JPY	250,000	2019	-52
	JPY	1,050,000		-308
Forward contract	CHF	3,000	2019	4
	CHF	3,000		4
Forward contract	GBP	3,500	2019	18
	GBP	3,500		18
Forward contract	CAD	2,000	2019	32
	CAD	4,000	2019	90
	CAD	2,000	2019	43
	CAD	8,000		165

Derivative financial instruments are measured at the time of acquisition at cost and at fair value on subsequent balance sheet dates. The changes in value of the derivative financial instruments are generally recognized in the statement of profit or loss on the reporting date. If the derivative financial instruments serve to hedge against cash flow risk arising from exchange rate risks and a qualified hedging relationship exists based on the criteria of IFRS 9, the valuation adjustments of the effective portion of the instrument are recognized in other comprehensive income (cumulative amount in 2018: €4.8 million; 2017: €16.4 million). Only the spot element of the forward contracts used to hedge the cash flow risks is designated as hedging instrument. The amounts recognized in equity are included in the profit or loss in the period in which the hedged transactions affect profit or loss. The changes of the hedging reserves are explained below as well as in the statement of changes

in equity. The non-designated or ineffective portion of the hedging instruments is recognized within the financial result in profit or loss.

The economic relationship between hedging instrument and hedged item and the effectiveness of the hedge relationship is determined based on consistency of the significant contractual features of the transactions ("Critical Terms Match"). In this regard, the Group performs a qualitative assessment. Hedge ineffectiveness may possibly arise when the timing of future transactions deviates from the original assumptions or the credit risk of the counterparties of the hedging instrument changes.

The following table presents the effects of the hedging instruments related to exchange rate risks on the financial position and performance of the Group:

Currency	Carrying amount (asset) as of Dec. 31, 2018	Carrying amount (liability) as of Dec. 31, 2018	Hedge ratio	Change in value of hedging instruments	Change in value of hedged item	Nominal amount	Maturity: 1-6 months	Maturity: 7-12	Average exercise price
	€ in K	€ in K		€ in K	€ in K	in K of respective currency			
USD	5,554	-661	100%	4,893	4,893	137,000	100,000	37,000	1.14
CHF	5	0	100%	5	5	3,000	3,000	0	1.13
CAD	194	0	100%	194	194	8,000	6,000	2,000	1.52
JPY	0	-302	100%	-302	-302	1,050,000	600,000	450,000	130.33
GBP	16	0	100%	16	16	3,500	3,500	0	0.90

Hedging instruments that have a positive fair value are shown in the line item "financial assets (non-current)" or "other financial assets (current)" in the statement of financial position. Hedging instruments that have a negative fair value are shown in the line item "other financial liabilities (non-current)" or "other financial liabilities (current)" in the statement of financial position.

The amounts that are recognized in the reporting period in connection with the cash flow hedges in other comprehensive income as well as those amounts that were reclassified from other comprehensive income to profit or loss (in the line item "sales revenue") are presented in the statement of other comprehensive income and the statement of changes in equity.

If the U.S. dollar would have depreciated 10% against the euro, the equity would have increased by €10.7 million (2017: increase by €13.2 million), the

impact on the result would have been +€1.6 million (2017: €0.1 million). Vice versa, if the U.S. dollar would have appreciated 10% against the euro, the resulting impact on the result would have been -€2.0 million (2017: -€0.1 million) and on the other comprehensive income -€13.1 million (2017: -€16.1 million).

36. Interest Risk Management

Sartorius Stedim Biotech is mainly financed through its parent company Sartorius AG. This major loan is taken out at variable interest rates; therefore the Group continues to be exposed to interest rate risks. To control the interest risk, an appropriate ratio between fixed and variable loans is maintained. As of December 31, 2018, the Group has no open interest rate derivative contracts to hedge the risk of increasing interest rates.

As of Dec. 31, 2018 the raised loans with variable interest rates amount to approx. €95 million.

If the market interest rate had been 1.0 percentage point higher, the interest expenses in the statement of profit or loss would have been €0.95 million (2017: €0.8 million) higher.

With regard to a decrease in interest rates a base interest rate of 0% has been considered. The impact on the financial result would have been €0.0 million (2017: –€0.2 million).

37. Liquidity Risk Management

The maturity of the financial liabilities excluding derivative financial instruments shows the following pattern:

	Carrying amount Dec. 31, 2017 € in K	Cash Flow Dec. 31, 2017 € in K	< 1 year € in K	1 - 5 years € in K	> 5 years € in K
Loans and borrowings	142,183	142,723	96,224	46,499	0
Finance Leases	17,487	32,825	1,900	7,558	23,368
Trade payables	77,211	77,211	77,211	0	0
Other liabilities (excluding derivatives)	63,428	68,965	33,036	35,780	150
Financial Liabilities	300,309	321,724	208,370	89,837	23,518

	Carrying amount Dec. 31, 2018 € in K	Cash Flow Dec. 31, 2018 € in K	< 1 year € in K	1 - 5 years € in K	> 5 years € in K
Loans and borrowings	132,942	133,191	89,997	43,194	0
Finance Leases	16,693	33,100	1,988	9,374	21,738
Trade payables	103,568	103,568	103,568	0	0
Other liabilities (excluding derivatives)	21,855	22,208	17,515	4,542	150
Financial Liabilities	275,059	292,068	213,069	57,111	21,888

The cash flows shown in the above tables include the undiscounted expected payments in connection with the respective financial liabilities including the associated interest payments based on the interest rates as of the balance sheet date.

The loans and borrowings include the loan raised from the parent company Sartorius AG. The other liabilities include the liability for the phantom units in AllPure.

The following tables illustrate the liquidity analysis for derivative financial instruments based on undiscounted cash flows:

	Carrying amount Dec. 31, 2017 € in K	Cash Flow Dec. 31, 2017 € in K	< 1 year € in K	1 - 5 years € in K	> 5 years € in K
Gross fulfilment					
Forward contracts	1,376	1,375	1,375	0	0
Payment obligation		22,308	22,308	0	0
Payment claim		-20,932	-20,932	0	0
Net fulfilment					
Interest rate swaps	552	552	552	0	0
Derivatives	1,928	1,927	1,927	0	0

	Carrying amount Dec. 31, 2018 € in K	Cash Flow Dec. 31, 2018 € in K	< 1 year € in K	1 - 5 years € in K	> 5 years € in K
Gross fulfilment					
Forward contracts	1,344	1,344	1,344	0	0
Payment obligation		42,177	42,177		
Payment claim		-40,833	-40,833		
Derivatives	1,344	1,344	1,344	0	0

The Group controls liquidity risks by maintaining credit lines and additional facilities with banks, by continuously tracking the forecasted and actual cash flows and by managing the maturity profiles of financial assets and liabilities. It is not expected that

the cash outflows will occur at significantly different times or at significantly different amounts.

The table below provides an overview of the credit lines available on the reporting date:

	Credit line at Dec. 31, 2017	< 1 year € in K	1 - 5 years € in K	> 5 years € in K	Interest rate	Credit line used at Dec. 31, 2017	Credit line unused as of Dec. 31, 2017
Loan from Sartorius AG	300,000	0	300,000	0	Variable	109,481	190,519
Bilateral loans	9,375	2,344	7,031	0	fix	9,375	0
Bilateral credit line	31,335	31,335	0	0	Variable	19,235	12,100
Total	340,710	33,679	307,031	0		138,091	202,619

	Credit line at Dec. 31, 2018	< 1 year € in K	1 - 5 years € in K	> 5 years € in K	Interest rate	Credit line used at Dec. 31, 2018	Credit line unused as of Dec. 31, 2018
Loan from Sartorius AG	309,601	0	309,601	0	Variable	94,500	215,101
Bilateral loans	6,250	3,125	3,125	0	fix	6,250	0
Bilateral credit line	23,163	23,163	0	0	Variable	16,750	6,413
Total	339,014	26,288	312,726	0		117,500	221,514

38. Credit Risk Management

Credit risk is the risk of financial loss to the Group if a counterparty to a financial instrument fails to meet its contractual obligations. Credit risk arises principally from cash and cash equivalents and trade receivables. In addition to that, the Group is exposed to credit risks arising from derivative financial instruments with positive fair values and, to a low extent, from contractual cash flows from debt securities.

Credit risk is controlled centrally for the Group by the Treasury Management unit. For counterparties such as banks and financial institutions the creditworthiness is continuously monitored in order to recognize increases in credit risks at an early stage. If no new information is obtained, the Group assumes that the related financial assets still have a low credit risk.

Customers are assigned risk limits that principally depend on the business volume, past experience and the financial position of the customer. Compliance with the limits is regularly reviewed by the management responsible. In some cases, the Group receives advance payments in order to avoid credit risks. There are no significant concentrations of credit risks from individual customers or regions.

For some trade receivables the Group possibly has collateral such as guarantees that can be used within the contractual agreements in case the counterparty does not meet its contractual payment obligations.

Impairment of Trade Receivables and Contract Assets

The new impairment model of IFRS 9, recognition of expected credit losses, is of particular relevance for the Group's trade receivables and contract assets according to IFRS 15. The Group applies the simplified approach according to IFRS 9 to trade receivables and contract assets. Accordingly, lifetime expected credit losses are recognized for these assets. The starting point of the new impairment model is an analysis of the actual historical credit loss rates. These are adjusted, taking into consideration forward-looking information and the effects of current changes in the macroeconomic environment, if significant. Due to the immaterial level of historical credit losses the Group currently determines the expected credit losses for the Group's portfolio of trade receivables as a whole. However, historical loss rates are analyzed regularly in more detail in order to apply different loss rates to different portfolios, where appropriate.

The contract assets are related to projects for typical customers of the Group. Therefore, it is assumed that the loss rates applied to trade receivables are appropriate approximations for the loss rates of the contract assets. Accordingly, there is no further differentiation between trade receivables and contract assets.

On this basis, the allowances for trade receivables and contract assets were determined as follows as of December 31, 2018 and as of the date of initial application of IFRS 9:

December 31, 2018 € in K	Not due	1 – 30 days overdue	31 – 60 days overdue	61 – 90 days overdue	More than 90 days overdue	Total
Gross carrying amount of trade receivables	133,707	39,781	17,502	6,115	22,843	219,947
Gross carrying amount of contract assets	3,902	0	0	0	0	3,902
Impairment loss allowance	77	24	49	280	3,188	3,617
January 1, 2018 € in K	Not due	1 – 30 days overdue	31 – 60 days overdue	61 – 90 days overdue	More than 90 days overdue	Total
Gross carrying amount of trade receivables	149,266	27,460	10,757	4,180	15,419	207,083
Gross carrying amount of contract assets	6,967	0	0	0	0	6,967
Impairment loss allowance	78	18	108	4	1,903	2,111

The expected credit losses are determined based on a loss rate of 0.05%. In addition, impairments are determined on the basis of individual assessments. Days overdue are one essential criterion in this context. A default is generally presumed when there is no

reasonable expectation of recovering a financial asset. In such a case, the respective receivables are derecognized.

The movements in the allowance for impairment in respect of trade receivables and contract assets in the reporting period are presented below:

	€ in K
Balance at January 1, 2018 under IAS 39	-2,433
Adjustment on initial application of IFRS 9	321
Balance at January 1, 2018 under IFRS 9	-2,111
Net remeasurement of loss allowance recognized in profit or loss in the reporting period	-2,041
Derecognition and consumption	39
Recoveries of amounts previously impaired	474
Foreign currency translation differences	23
Balance at December 31, 2018 under IFRS 9	-3,617

Impairment of Other Financial Assets

Besides trade receivables, cash and cash equivalents are the most material financial assets on the Group's statement of financial position as of the reporting date, December, 2018. The expected credit losses are monitored at regular intervals. Due to the high creditworthiness of the counterparties and the short maturities, the impairment which would have to be recognized for these financial assets is immaterial. Therefore, no impairment is recognized for cash and cash equivalents.

For the other financial assets measured at amortized cost no impairment is recognized as of December 31, 2018 for the twelve months of expected credit losses due to immaterial historical credit losses. In the event of a significant increase in credit risk, which is generally presumed when a payment is more than 30 days past due, the lifetime expected credit losses are

recognized for the respective financial asset. A default is generally presumed when there is no reasonable expectation of recovering a financial asset. This is generally presumed when payments are more than 90 days past due. As of the reporting date, there are no indications of increases in credit risk to a material extent. The carrying amounts of the financial assets reflect the maximum credit loss for these assets at the end of the reporting period.

39. Other Risks Associated with Financial Instruments

As of the reporting date, the Sartorius Stedim Biotech Group has not been exposed to the risk of volatility in share prices.

40. Share-based Payments

On the level of Sartorius Stedim Biotech's majority shareholder Sartorius AG, share-based payments exist in the form of so-called phantom stock units. Under this plan the respective board member is granted a certain number of phantom stocks each year that represent an agreed amount of money. The exercise of these stocks is not possible before four years and is depending on certain requirements regarding the performance of the Sartorius AG shares.

When the stocks are paid out the amount is based on the share price at the exercise date. The payment is capped at an amount of 2.5 times the share price at the time these virtual options were granted. For further details please refer to the Remuneration Report.

The fair value of the phantom stock units is disclosed as follows:

	Number of phantom stock units	Subscription price in €	Fair value when granted on Jan. 1 of the particular year € in K	Fair value at year-end on Dec. 31, 2018 € in K	Paid out € in K	Exercisable
Tranche of phantom stock units for 2014	8,032	21.01	169	0	422	
Tranche of phantom stock units for 2015	7,360	24.70	182	454	0	no
Tranche of phantom stock units for 2016	3,484	57.41	200	359	0	no
Tranche of phantom stock units for 2017	2,950	70.51	208	281	0	no
Tranche of phantom stock units for 2018	2,685	80.32	216	239	0	no
Total	24,511		974	1,333	422	

Other Disclosures

The consolidated financial statements were prepared on a going concern basis.

Material Events after the Reporting Date

No material events occurred after the reporting date.

Number of Employees

The average workforce employed during the year 2018 was 5,412 (4,991 in 2017).

41. Related Parties

General

The majority shareholder of Sartorius Stedim Biotech S.A. is Sartorius AG, which holds a controlling stake in the company of 74.3% in equity capital – and 84.5% of the voting rights. The Sartorius Group itself is organized in two divisions: Bioprocess Solutions (mainly run by the Sartorius Stedim Biotech Group) and Lab Products & Services (mainly run by the other companies of Sartorius Group). This structure leads to the fact that the Group holds two subsidiaries in most of the countries and these companies partially share space, staff and other resources. Furthermore, the German group companies carry out various central functions and accordingly deliver services to the worldwide entities (e.g. IT support). The company Sartorius Corporate Administration GmbH, a 100% subsidiary of Sartorius AG has incorporated numerous Group functions like Group Finance, HR, IT, Investor Relations and Legal. These services are charged within the Group and to a significant extent also to Sartorius Stedim Biotech.

The described structures lead to various relations and transactions with related parties. Transactions between Sartorius Stedim Biotech S.A. and its subsidiaries (presented in Note 7), which are related parties of the company, have been eliminated on consolidation and are not disclosed under this Note. Details of transactions between the Group and other related parties, belonging to the Sartorius Group, are disclosed below.

Sales, Purchases and Commissions

In certain business areas members of the Sartorius Group act as contract manufacturers for the Sartorius Stedim Biotech Group and vice versa. The respective transactions are carried out at arms' length principles and are disclosed in the table below as "sales revenue" and "purchases".

	Sales revenue 2018 € in K	Purchases 2018 € in K
Related parties of Sartorius Group	72,221	7,061
	Sales revenue 2017 € in K	Purchases 2017 € in K
Related parties of Sartorius Group	60,292	5,085

Certain product groups of the Sartorius Stedim Biotech portfolio are sold through the sales force of other Sartorius entities. For the arranging of the sale the Sartorius Stedim Biotech Group has paid commissions in the amount of €0.4 million (€1.1 million in 2017). These commissions are typically calculated as a percentage of the generated sales revenue.

Management Fees and Other Shareholder Costs

One of Sartorius Stedim Biotech S.A.'s board members is also member of the Sartorius AG Executive Board and is paid by the German parent company. For his services for Sartorius Stedim Biotech a portion of his remuneration is charged to Sartorius Stedim Biotech S.A. (€0.7 million in 2018; €0.8 million in 2017) and charged to Sartorius Stedim Biotech GmbH (€0.7 million in 2018; €0.8 million in 2017).

Other shareholder functions like Group Financial Reporting, Compliance and Investor Relations are performed by the above mentioned Sartorius Corporate Administration GmbH in Germany. These services have been charged to Sartorius Stedim Biotech S.A. in the amount of €0.8 million (2017: €0.8 million).

Shareholder Loan

As described in note 30 the Sartorius Stedim Biotech Group has raised a loan from its parent company Sartorius AG with a credit volume of approx. €300 million and a current utilization of approx. €95 million (2017: €109 million). The interest charged is based on a variable interest rate plus an arms'-length credit margin.

Administration Charges and Shared Costs

As described above the companies in most of the countries share certain functions and costs. The underlying contracts include mainly subleases for office space and central administrative functions, such as accounting and controlling, human resource

management and IT. In this respect, the relevant companies charge rent, salaries, social security costs and other expenses for such services, as well as a pro-rated profit margin for the services they provide.

The most significant contract in this context is the one between Sartorius Stedim Biotech GmbH, Germany, and Sartorius Corporate Administration GmbH. This company provides all central service and administrative functions to Sartorius Stedim Biotech GmbH and other Group companies. The calculation for services fees typically includes a surcharge of 3% on total costs. 3% is a surcharge compliant with arm's length principles for routine tasks, following OECD and EU guidelines. In 2018, services for approx. €53.4 million were provided to Sartorius Stedim Biotech GmbH (€40.1 million in 2017). This amount covers the following functions:

- Marketing Communication, e-Business, Business Development
- Environment, Health & Security, Factory Maintenance
- Finance, Human Resources, Information Technology
- Central Services & General Organization.

During the reporting period, the Sartorius Stedim Biotech Group acquired a perpetual right to use the Sartorius global process template for some of its entities for a total amount of €19.9 million. This transaction enhances the independence of the Sartorius Stedim Biotech Group from Sartorius Corporate Administration GmbH and will lead to a corresponding reduction in future IT service charges.

Compensation of Key Management Personnel

In 2017 and 2018, the Executive Board Management received the following remuneration:

	Total € in K	Short-term benefits € in K	Post-employment benefits € in K	Other long-term benefits € in K	Termination benefits € in K	Share-based payments € in K
2018¹⁾	2,522	1,333	257	189	0	743
2017 ¹⁾	3,056	1,210	258	184	0	1,404

¹⁾ For more information please refer to the chapter Corporate Governance (See pages 57 to 87)

Statutory Auditors' Report on the Consolidated Financial Statements

This is a translation into English of the statutory auditors' report on the consolidated financial statements of the Company issued in French and it is provided solely for the convenience of English speaking users.

This statutory auditors' report includes information required by European regulation and French law, such as information about the appointment of the statutory auditors or verification of the information concerning the Group presented in the management report.

This report should be read in conjunction with, and construed in accordance with, French law and professional auditing standards applicable in France

For the year ended 31 December 2018

To the Shareholders of SARTORIUS STEDIM BIOTECH S.A.,

Opinion

In compliance with the assignment entrusted to us by your shareholders' meetings, we have audited the accompanying consolidated financial statements of SARTORIUS STEDIM BIOTECH S.A. for the year ended 31 December 2018.

In our opinion, the consolidated financial statements give a true and fair view of the assets and liabilities and of the financial position of the Group as at 31 December 2018 and of the results of its operations for the year then ended in accordance with International Financial Reporting Standards as adopted by the European Union.

The audit opinion expressed above is consistent with our report to the Audit Committee.

Basis for Opinion

Audit Framework

We conducted our audit in accordance with professional standards applicable in France. We believe that the audit evidence we have obtained is sufficient and appropriate to provide a basis for our opinion.

Our responsibilities under those standards are further described in the Statutory Auditors' Responsibilities for the Audit of the Consolidated Financial Statements section of our report.

Independence

We conducted our audit engagement in compliance with independence rules applicable to us, for the period from 1 January 2018 to the date of our report and specifically we did not provide any prohibited non-audit services referred to in Article 5(1) of Regulation (EU) No 537/2014 or in the French Code of ethics (code de déontologie) for statutory auditors.

Justification of Assessments – Key Audit Matters

In accordance with the requirements of Articles L.823 - 9 and R.823 - 7 of the French Commercial Code (code de commerce) relating to the justification of our assessments, we inform you of the key audit matters relating to risks of material misstatement that, in our professional judgment, were of most significance in our audit of the consolidated financial statements of the current period, as well as how we addressed those risks.

These matters were addressed in the context of our audit of the consolidated financial statements as a whole, and in forming our opinion thereon, and we do not provide a separate opinion on specific items of the consolidated financial statements.

Goodwill valuation – Impairment test

Identified risk

As of 31 December 31 2018, goodwill amounts to 384.7 M€, or 25% of total consolidated assets.

As described in Note 5 to the consolidated financial statements, SARTORIUS STEDIM BIOTECH is an "integrated solution provider" for its customers, and as a result there is only one operating segment from a product and customer perspective: "Biopharma ". In addition, as indicated in Note 16 to the consolidated financial statements, because of the interdependence of the market in which your group operates, the

lowest level at which goodwill can be allocated is the Biopharma segment. The goodwill has therefore been fully allocated to the Cash Generating Units (C.G.U.) corresponding to the Biopharma segment.

Goodwill is subject to an annual impairment test and whenever there are indicators of impairment in accordance with the methods and assumptions described in Notes 3 and 16 to the consolidated financial statements. In particular, in view of what has been described above, the impairment test is carried out at the level of the Biopharma segment.

We considered that the determination of the value of goodwill is a key point of our audit given its significance in the consolidated financial statements of your group, and because the determination of the recoverable amount taken into account in the impairment test on the basis of the value in use of the C.G.U. requires the use of estimates and assumptions (including future cash flows, perpetual revenue growth rates and the discount rate) that require significant management judgment.

Responses obtained during our audit

We obtained the impairment test from the C.G.U. group corresponding to the Biopharma segment as well as the forecasts underlying the calculation (4-year plan).

We reviewed the compliance of the company's methodology in accordance with applicable accounting standards.

We also performed a critical review of how the company has implemented this methodology, including the following procedures:

- Assess the reasonableness of the key assumptions used to determine the cash flow of the Biopharma segment as well as that used for the perpetual growth rate;
- Assess, with the support of our valuation specialists, the discount rate used by management. We compared this rate with our own estimates and analysed its various constituent elements;
- Check the arithmetical accuracy of the impairment test performed by your group.

We have also obtained and assessed the sensitivity analyses carried out by management, as shown in Note 16 of the notes to the consolidated financial statements. As a result, we were able to verify that

only an extremely large change in the main assumptions could lead to the recognition of an impairment of goodwill.

Lastly, we verified the appropriateness of the information provided in Notes 3 and 16 to the consolidated financial statements.

Specific verifications

As required by French law and regulations, we have also verified in accordance with professional standards applicable in France the information pertaining to the Group presented in the management report of the Board of Directors.

We have no matters to report as to its fair presentation and its consistency with the consolidated financial statements

Report on Other Legal and Regulatory Requirements

Appointment of the Statutory Auditors

We were appointed as statutory auditors of SARTORIUS STEDIM BIOTECH S.A. by the annual general meeting held on 7 April 2015 for KPMG S.A. and on 19 May 2006 for Deloitte & Associés.

As at 31 December 2018, KPMG S.A. was in its 4th year of the audit mandate without interruption and Deloitte & Associés was in its 13th year.

Responsibilities of Management and Those Charged with Governance for the Consolidated Financial Statements

Management is responsible for the preparation and fair presentation of the consolidated financial statements in accordance with International Financial Reporting Standards as adopted by the European Union and for such internal control as management determines is necessary to enable the preparation of consolidated financial statements that are free from material misstatement, whether due to fraud or error.

In preparing the consolidated financial statements, management is responsible for assessing the Company's ability to continue as a going concern, disclosing, as applicable, matters related to going concern and using the going concern basis of

accounting unless it is expected to liquidate the Company or to cease operations.

The Audit Committee is responsible for monitoring the financial reporting process and the effectiveness of internal control and risks management systems and where applicable, its internal audit, regarding the accounting and financial reporting procedures.

The consolidated financial statements were approved by the Board of Directors.

Statutory Auditors' Responsibilities for the Audit of the Consolidated Financial Statements

Objectives and audit approach

Our role is to issue a report on the consolidated financial statements. Our objective is to obtain reasonable assurance about whether the consolidated financial statements as a whole are free from material misstatement. Reasonable assurance is a high level of assurance, but is not a guarantee that an audit conducted in accordance with professional standards will always detect a material misstatement when it exists. Misstatements can arise from fraud or error and are considered material if, individually or in the aggregate, they could reasonably be expected to influence the economic decisions of users taken on the basis of these consolidated financial statements.

As specified in Article L.823-10-1 of the French Commercial Code (code de commerce), our statutory audit does not include assurance on the viability of the Company or the quality of management of the affairs of the Company.

As part of an audit conducted in accordance with professional standards applicable in France, the statutory auditor exercises professional judgment throughout the audit and furthermore:

- Identifies and assesses the risks of material misstatement of the consolidated financial statements, whether due to fraud or error, designs and performs audit procedures responsive to those risks, and obtains audit evidence considered to be sufficient and appropriate to provide a basis for his opinion. The risk of not detecting a material misstatement resulting from fraud is higher than for one resulting from error, as fraud may involve collusion, forgery, intentional omissions, misrepresentations, or the override of internal control. Obtains an understanding of internal control relevant to the audit in order to design audit

procedures that are appropriate in the circumstances, but not for the purpose of expressing an opinion on the effectiveness of the internal control.

- Obtains an understanding of internal control relevant to the audit in order to design audit procedures that are appropriate in the circumstances, but not for the purpose of expressing an opinion on the effectiveness of the internal control.
- Evaluates the appropriateness of accounting policies used and the reasonableness of accounting estimates and related disclosures made by management in the consolidated financial statements.
- Assesses the appropriateness of management's use of the going concern basis of accounting and, based on the audit evidence obtained, whether a material uncertainty exists related to events or conditions that may cast significant doubt on the Company's ability to continue as a going concern. This assessment is based on the audit evidence obtained up to the date of his audit report. However, future events or conditions may cause the Company to cease to continue as a going concern. If the statutory auditor concludes that a material uncertainty exists, there is a requirement to draw attention in the audit report to the related disclosures in the consolidated financial statements or, if such disclosures are not provided or inadequate, to modify the opinion expressed therein.
- Evaluates the overall presentation of the consolidated financial statements and assesses whether these statements represent the underlying transactions and events in a manner that achieves fair presentation.
- Obtains sufficient appropriate audit evidence regarding the financial information of the entities or business activities within the Group to express an opinion on the consolidated financial statements. The statutory auditor is responsible for the direction, supervision and performance of the audit of the consolidated financial statements and for the opinion expressed on these consolidated financial statements.

Report to the Audit Committee

We submit a report to the Audit Committee which includes in particular a description of the scope of the audit and the audit program implemented, as well as the results of our audit. We also report, if any, significant deficiencies in internal control regarding

the accounting and financial reporting procedures that we have identified.

Our report to the Audit Committee includes the risks of material misstatement that, in our professional judgment, were of most significance in the audit of the consolidated financial statements of the current period and which are therefore the key audit matters, that we are required to describe in this audit report.

We also provide the Audit Committee with the declaration provided for in Article 6 of Regulation (EU) N° 537/2014, confirming our independence within the meaning of the rules applicable in France such as they are set in particular by Articles L.822-10 to L.822-14 of the French Commercial Code (code de commerce) and in the French Code of Ethics (code de déontologie) for statutory auditors. Where appropriate, we discuss with the Audit Committee the risks that may reasonably be thought to bear on our independence, and the related safeguards.

Marseille, 8 February 2019

The Statutory Auditors

French original signed by

KPMG Audit
A division of KPMG S.A

Deloitte & Associés

John Evans

Vincent Gros

