

1870

Founded by Florenz Sartorius,  
headquartered in Göttingen, Germany

60+

Production and sales sites worldwide

>14,600

Employees

~16%

Sales CAGR 2013–2023

+7.8pp

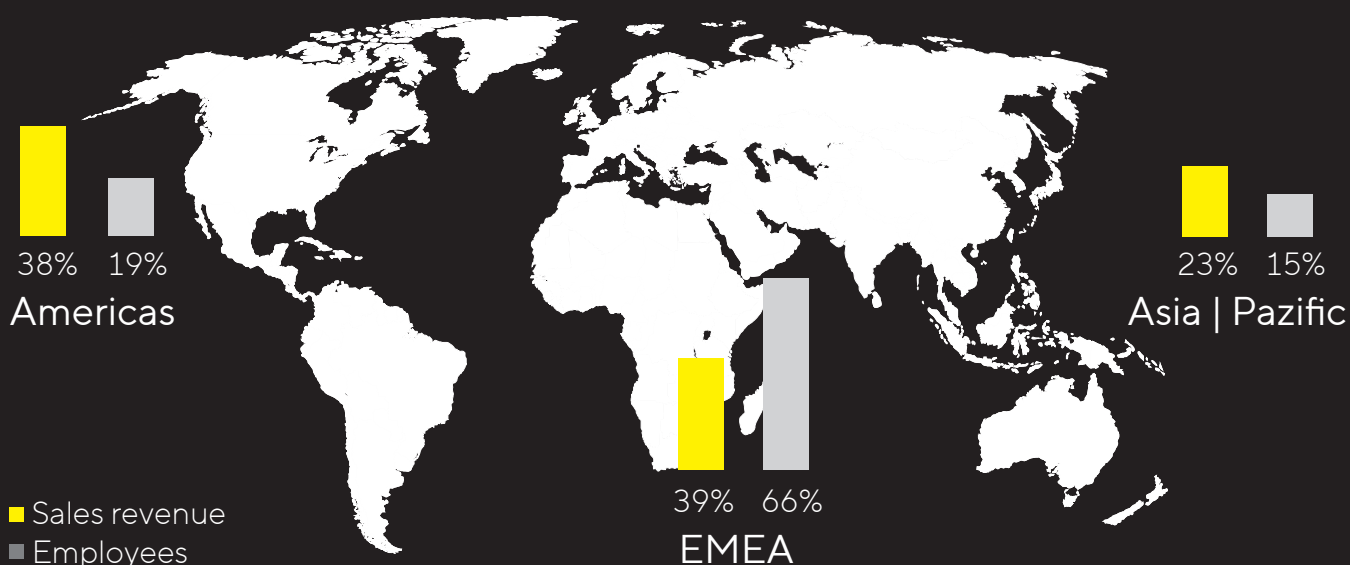
Change in underlying  
EBITDA margin 2013–2023

~€20.5bn

Sartorius AG market capitalization;  
listed on the DAX and TecDAX

Sales growth for continued operations, underlying = excluding extraordinary items

## Strong Presence in All Major Biopharma Markets



# Innovative Solutions for Better Medications

With its pioneering spirit and a profound understanding of customer requirements, Sartorius has evolved throughout its 150-year history into a key partner for biopharmaceutical research and the industry. Our goal is to make complex and expensive development of biotech medicines and their production safer and more efficient. We cover the entire value-added chain of the biopharmaceutical industry and help with our products and services to ensure that novel therapies and vaccines reach the market faster and are accessible to more people worldwide.

*See page 25, Sartorius Group at a Glance*

## Mission

We empower scientists and engineers to simplify and accelerate progress in life science and bioprocessing, enabling the development of new and better therapies and more affordable medicine.



## Vision

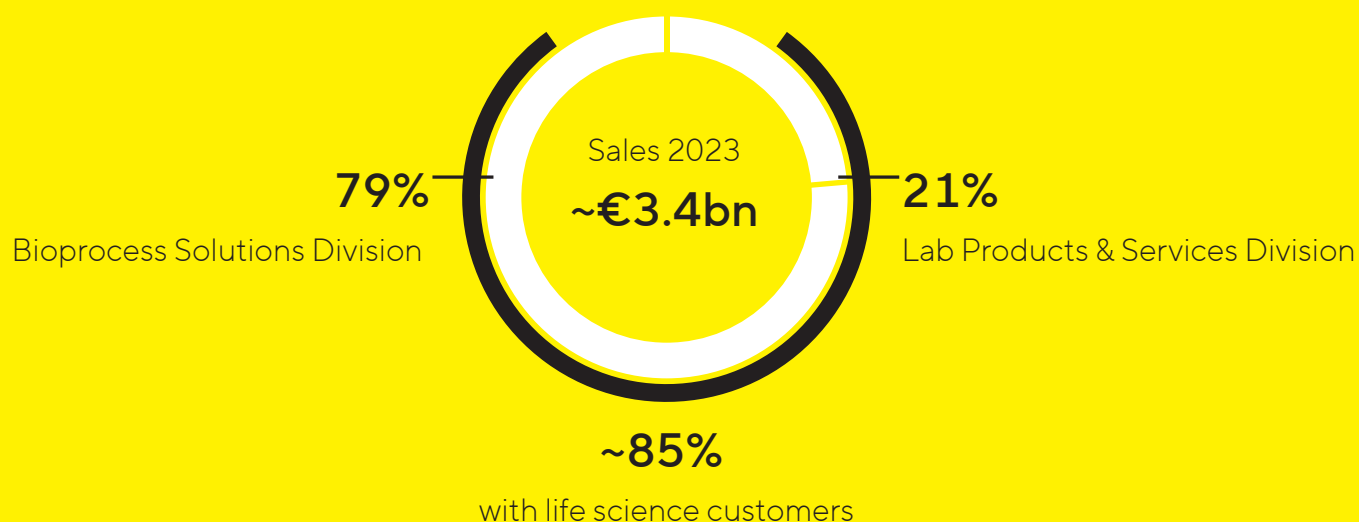
We are a magnet and dynamic platform for pioneers and leading experts in our field. We bring creative minds together for a common goal: technological breakthroughs that lead to better health for more people.

# Bioprocess Solutions

In the Bioprocess Solutions Division, Sartorius offers a broad product portfolio that covers all steps in the production of a biopharmaceutical. The company has held leading market positions for years in its core technologies, such as filtration, fermentation, cell cultivation, and fluid management.



We Operate in Two Divisions with a Clear Focus on the Life Science Industry



## Lab Products & Services

The Lab Products & Services Division offers laboratories in the pharmaceutical and biopharmaceutical industries as well as at academic research institutes innovative solutions for bioanalytics, in addition to premium laboratory products, consumables, and services. Sartorius is among the market leaders in laboratory balances, pipettes, and lab consumables.