

25+

Sites in more than 25 countries,  
headquartered in Aubagne France

>10,400

Employees

~19%

Sales CAGR 2011-2021

~90%

Sales share with life  
science customers

+14.9pp

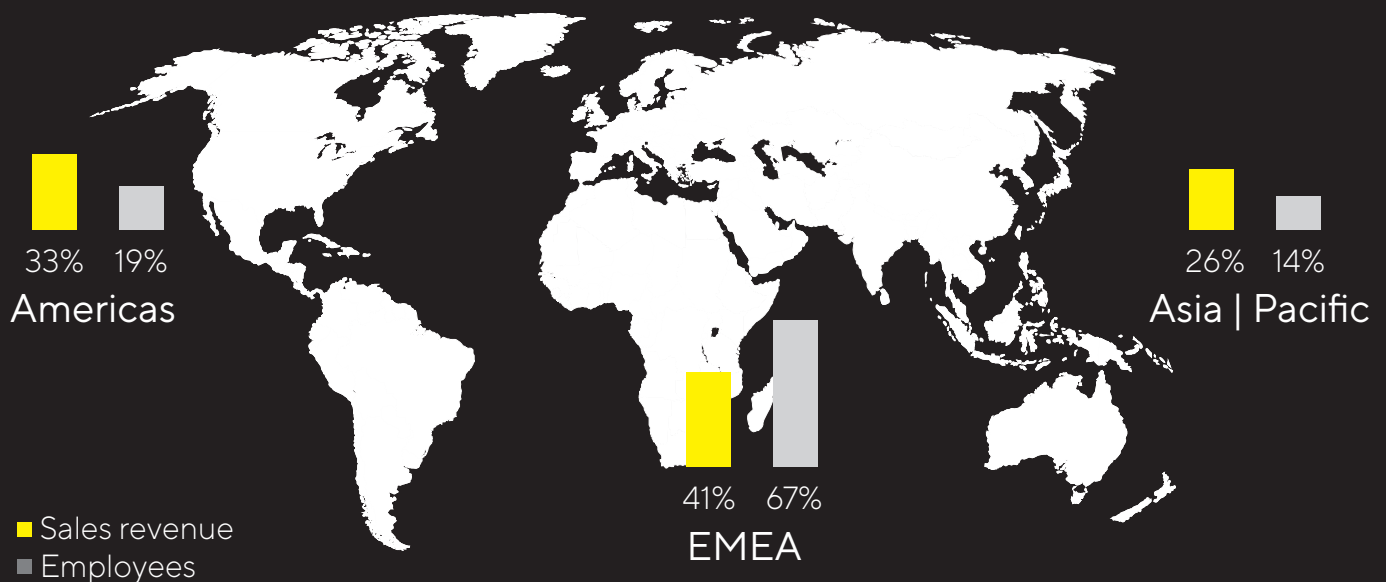
Change in underlying  
EBITDA margin 2011-2021

~€44.5bn

Sartorius Stedim Biotech S.A.  
market capitalization; listed on  
the CAC Large 60

Sales growth in constant currencies; 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 Stedim Biotech has evolved 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 18, Sartorius Stedim Biotech Group at a glance*

## Mission

At Sartorius Stedim Biotech, we empower engineers to simplify and accelerate progress in bioprocessing. In this way, we enable new and better pharmaceuticals to be manufactured and help keep medications affordable.



## 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.