

SARTORIUS

Working at Sartorius

Where Your Growth and Ideas Simplify Progress



This is Sartorius

Many diseases, such as cancer, dementia or inflammatory diseases, are still incurable. Hundreds of millions of people lack access to affordable and safe medicines. Our biopharma customers are working hard to develop new drugs to improve people's lives.

Sartorius is a leading international partner to biopharmaceutical research and industry. With our innovative technologies, we aim to simplify progress in the development and production of new therapeutics. So that more people will have access to better medicine.

More than 14,000 employees at over 60 production and sales locations worldwide are committed to this vision.

As a growing global life science company, we are looking for professionals, graduates and students who share our ambition and want to make a contribution in a vibrant environment while advancing their careers.

Join our global team and grow with us.



- 3 This is Sartorius
- 4 Sartorius in Numbers
- 6 What Drives Us
- 7 Corporate Values
- 8 Simplifying Medical Progress
- 10 At Home Worldwide
- 12 Where Your Growth and Ideas Simplify Progress
- 14 Your Benefits at Sartorius
- 15 Grow With Us

Sartorius in Numbers

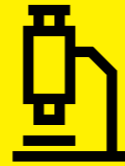
Sartorius, founded in 1870 in Göttingen, Germany, has evolved into a leading international partner to biopharmaceutical research and industry.

Back then and today, our innovative solutions are helping our customers to accelerate research work and manufacture biotech medications, vaccines and cell and gene therapies more safely, rapidly and sustainably.

~3.5 bn
Euros Sales Revenue in 2025



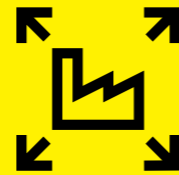
~85%
Sales With Life Science Customers



60+
Locations Worldwide



440 mn
Euros Investment in 2025



2045
Ambition for Climate Neutrality



1870
Year of Foundation



A Diverse and Ambitious Team

We believe that diverse perspectives are crucial to better understanding our customers and finding innovative solutions. That's why we foster diversity in our global team.



Explore more about people and diversity.
www.sartorius.com/diversity

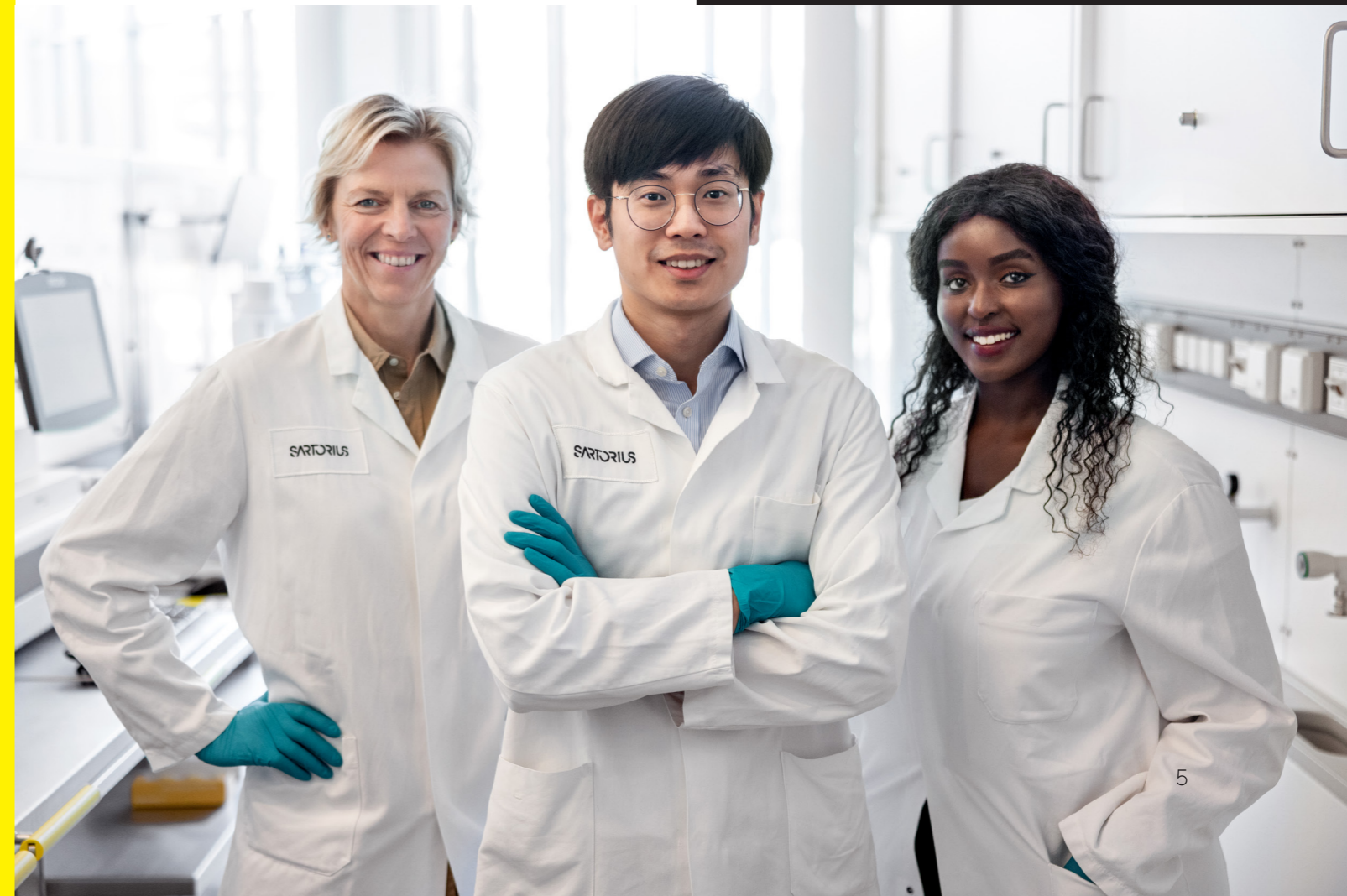
+14,000 Employees

100+ Nationalities




41 Average Age

39% Share of Women

76% Internally Filled Management Positions



What Drives Us

Our Purpose	Better health for more people.	
Our Mission	We empower biopharma customers to accelerate drug development and enhance manufacturing efficiency for better therapies and more affordable medicine.	
Our Promise	Simplifying Progress	



Corporate Values

Sustainability, openness and enjoyment are the values that shape our work with colleagues, customers and business partners alike. By incorporating them into our thinking and actions on a daily basis, we create a positive corporate culture, which is an important source of long-term economic success, lasting motivation and above-average performance.



Sustainability

Growing profitably and acting responsibly



Openness

A source of change and progress



Enjoyment

A working environment with freedom and appreciation

Simplifying Medical Progress

With our product portfolio, we focus on the biopharmaceutical industry. By offering products and services covering the entire value chain, Sartorius is part of the solution in the fight against diseases worldwide. We support customers in increasing efficiency in drug development and production.

Bioprocess Solutions

With its broad product portfolio covering filtration, fermentation, membrane chromatography and fluid management, Sartorius helps ensure that biotech drugs and vaccines can be produced safely and efficiently.

Cell Cultivation, Media, Reagents and Cell Culture Components



Fluid Management



Filtration



Purification



Bioprocessing Portfolio



Lab Portfolio

Lab Instruments and Consumables



Bioanalytical Instruments



Lab Products & Services

Sartorius aims to speed up the time-consuming and expensive development process for biopharmaceuticals with innovative solutions for bioanalytics and premium lab products, consumables and services.

At Home Worldwide

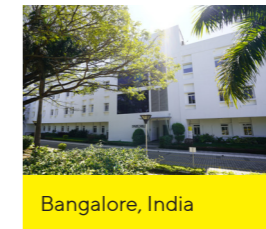
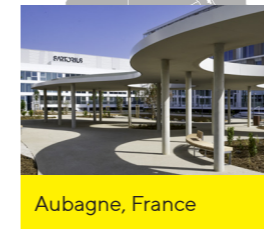
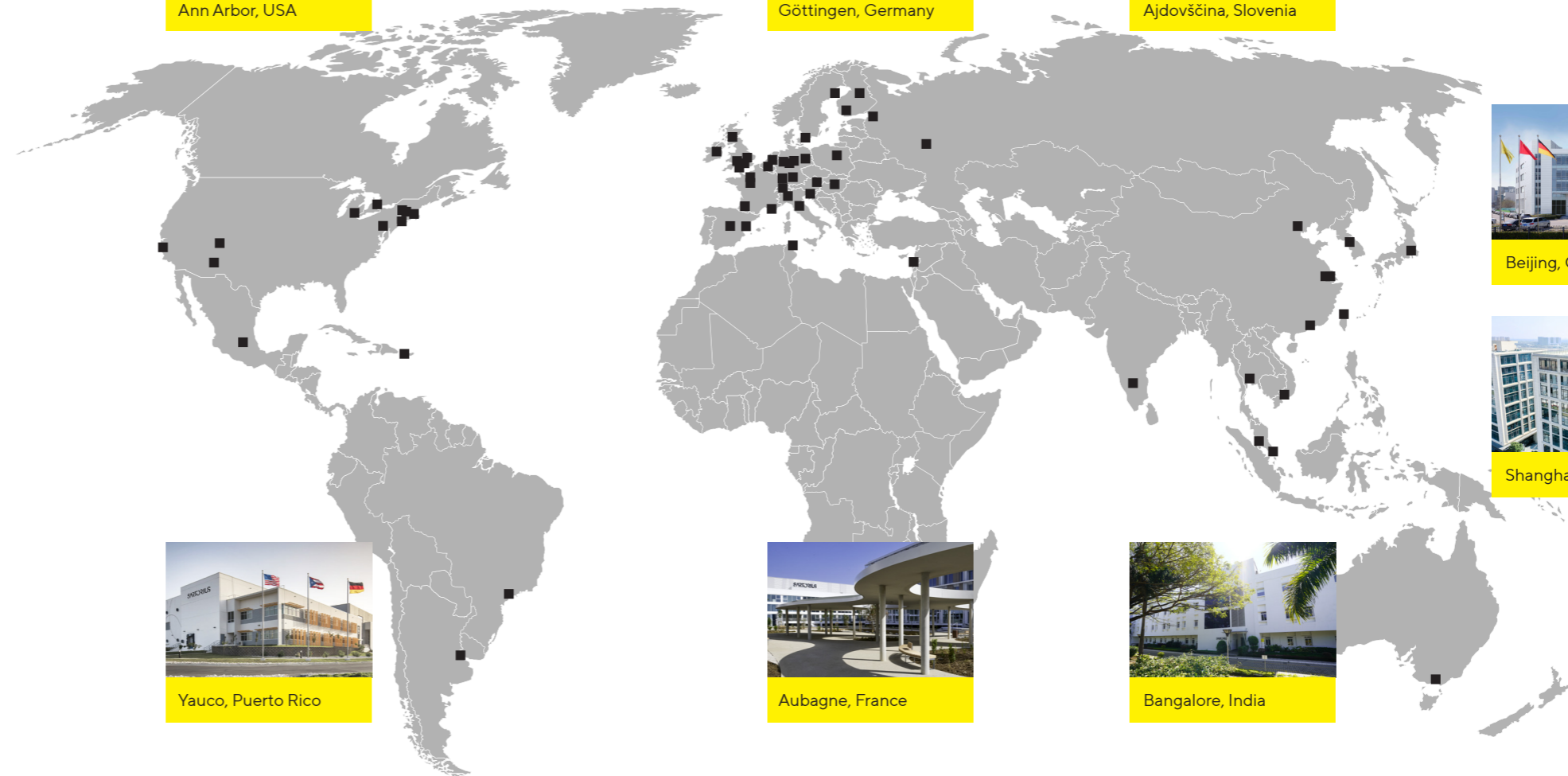
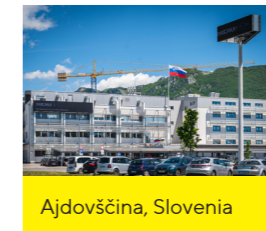
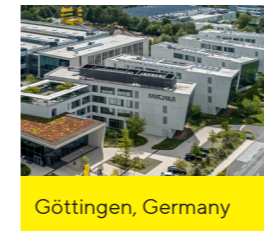
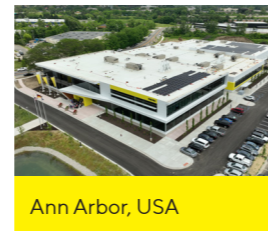
Sartorius focuses its business activities on the major biopharma markets, earning more than 85% of its sales revenue outside of Germany, the company's country of origin.

That is why we feel equally at home in the company's hometown of Göttingen as in any of the 60+ countries, where our production facilities, product development activities, customer and testing laboratories as well as sales and service teams are present.

Therefore, we are looking for people who think beyond borders and want to proactively strive to develop these markets further along with us for our international teams.



Find a current list of the Sartorius locations worldwide here



Where Your Growth and Ideas Simplify Progress

Sartorius unites life science enthusiasts, doers, risk takers, creative minds, and, above all, partners for our customers. They all share one ambition: Accelerating the development and production of innovative therapies – for better health for more people. And it all starts with the ones who move forward. Like you.

Be Part of the Solution

Sartorius is part of the solution in overcoming global diseases such as cancer and dementia. Our technologies help researchers and manufacturers to move faster from barrier to breakthrough, simplifying progress in the development and production of new therapeutics. For our people, this purpose is both a commitment and a chance. If you bring technical expertise, innovative thinking, and ambition, at Sartorius you will be part of the solution.

Grow With Us

At Sartorius, growth is more than a metric. We focus on dynamic markets that will continue to grow for decades. This creates opportunities for those who seize them and take risks to move forward. Depending on goals and changing priorities in life, you can find new roles, and discover new areas of expertise – driving your own development as well as that of our business. That's how personal growth at Sartorius is not a cherry on top. It's the bread and butter.

Add Your Spark

Ambitious pioneers and creative minds at Sartorius thrive in a unique blend of entrepreneurial spirit and global scope: Your ideas could be the spark that leads to the next game-changing discovery. Yet we know progress does not come from a single view, but from a multitude of perspectives that challenge one another and refine solutions. While great minds think alike, at Sartorius, greater minds think together – and build on each other. Those who boldly contribute new ideas prosper in a culture that encourages each individual to realize their ambitions.



Valérie Morillon

Leads the sales team for lab essentials in North and South Europe

“After several years in another company, I wanted to use my experience to serve customers and employees in a growing business such as Sartorius. I hoped that in a growing company, I would find an innovative culture and exciting challenges. Looking back, I can say I definitely found this in Sartorius.”



Muhamed Karalić

Head of Operations for several bioprocess production sites in Europe and the U.S.

“I started as a clerk at Sartorius in 2013 – today, I am responsible for bioprocess operations in countries across Europe and the U.S. This was only possible because Sartorius gave me the opportunity to continuously develop – both professionally and personally. I am proud to be part of a company involved in the manufacturing of products that are used in the fight against diseases worldwide.”



Susan Xu

Responsible for supply chain management in China

“Over my more than 25 years at Sartorius, I’ve often stepped outside my comfort zone and embraced change – and that’s where real growth happened. I started as an order processing specialist, and today, I’m leading supply chain management for bioprocessing solutions and lab products in China. At Sartorius, every challenge along the way is an opportunity for personal and professional growth.”



Read more career stories on the Sartorius blog.
www.sartorius.com/blog



Your Benefits at Sartorius

We are flexible, so you can give your best. Depending on region, job requirements and local policies, Sartorius offers benefits, such as:



Career Development | Actively advance your career at one of the fastest growing global life science companies with diverse development opportunities—from technical to geographical.



Well-Being | Benefit from a wide range of health support: from safe working environment, to healthcare insurance, to preventive medical check-ups, to sport activities or even gyms at some sites—and more.



Work-Life Balance | Enjoy the flexibility your life needs: Flexible hours and home office options (depending on position and region) as well as support with childcare services and elder care at some sites.



Smart Working Environment | Enjoy working in state-of-the-art facilities with the latest technology and equipment that supports global collaboration and encourages your creativity.



Open and Highly Collaborative Culture | Engage in our diverse, global team and experience the freedom to contribute through agile working methods, short decision-making paths and mutual support.



Regional Benefits | Discover additional advantages depending on local policies: From company restaurants, to subsidies for public transport, to bicycle and car leasing, to student loan tuition refinance service and many more.

Grow With Us

The company's dynamic growth creates opportunities to develop personally and professionally for those who seize them. We aim to fill most management positions from within our own ranks and provide development initiatives from the very beginning:

Onboarding | Global and site-specific onboarding programs aim to achieve your social, cultural and professional integration.

Learning | Our learning platform offers a broad range of internal and external trainings, with both digital and in-person formats: from product and application trainings, to soft skill and leadership development, to IT, software and languages programs.

Development | To foster personal and professional development, we are committed to annual performance reviews, talent identification initiatives, leadership trainings as well as coaching and mentoring offers.

Feedback | Drive your personal development through annual performance reviews and co-create your working environment through our regular global employee surveys.



It all starts
with the ones
who move
forward.
Like you.

Visit www.sartorius.com/careers
or find us on LinkedIn.

Simplifying Progress