

# SARTORIUS

## Working at Sartorius

Candidates' Guide



# 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 life sciences research and the biopharmaceutical 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 13,500 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 Mission & Vision
- 7 Corporate Values
- 8 Simplifying Progress in Life Sciences and Bioprocessing
- 10 At Home Worldwide
- 12 Why Work At Sartorius
- 14 Benefit From Attractive Working Conditions
- 15 Advance Your Career





# Sartorius in Numbers

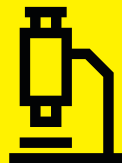
Sartorius, founded in 1870 in Göttingen, Germany and listed on the German DAX since 2021, has evolved into a leading international partner to life sciences research and the biopharmaceutical industry.

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.

~3.4 bn  
Euros Sales Revenue 2024



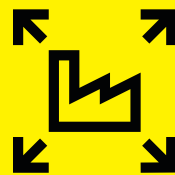
~85%  
Sales With Life Science Customers



60+  
Locations Worldwide



410 mn  
Euros Investment in 2024



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](http://www.sartorius.com/diversity)

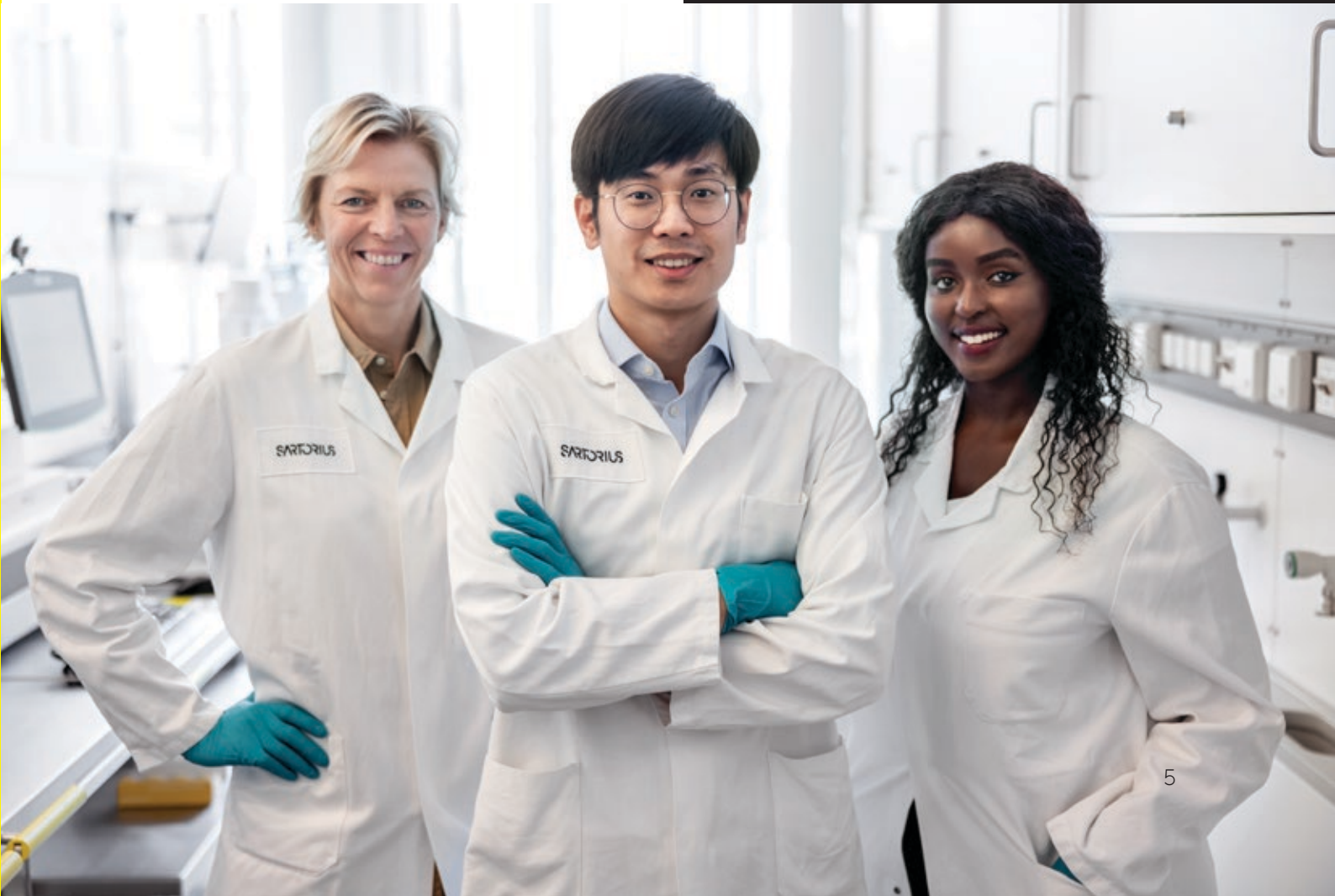
+13,500  
Employees

110+  
Nationalities

41  
Average Age

39%  
Share of Women

52%  
Internally Filled Positions



# Mission

At Sartorius, 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.

# 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 towards all stakeholders



## Openness

A source of change and progress



## Enjoyment

A working environment with freedom and appreciation



# Simplifying Progress in Life Sciences and Bioprocessing

With its two divisions, Bioprocess Solutions and Lab Products & Services, Sartorius focuses on the biopharmaceutical industry. Offering solutions covering the entire value chain, the company supports its 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



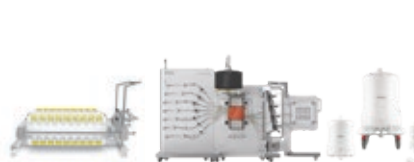
Fluid Management



Filtration



Purification



## Bioprocess Solutions Division

Molecule Development



Cell Line and Process

Development



Upstream and Downstream Production

Quality Control and Testing

## Lab Products & Services Division

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](#)



Ann Arbor, USA



Göttingen, Germany



Ajdovščina, Slovenia



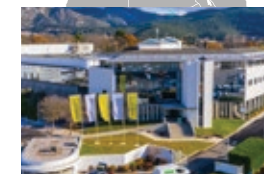
Beijing, China



Shanghai, China



Yauco, Puerto Rico



Aubagne, France



Bangalore, India



# Why Work at Sartorius

Sartorius unites life science enthusiasts, doers, risk takers, creative minds, and, above all, partners for our customers. They all share one ambition: translating scientific discoveries into medicine faster and more efficiently. As a globally listed company, Sartorius has retained the mindset that made it a pioneer in its industry as a start-up more than 150 years ago: customer focus, can-do attitude, hands-on approach, and each individual’s ambition to contribute.

### Be Part of the Solution

Sartorius is part of the solution in overcoming global diseases such as cancer or dementia. For our people, this purpose is both a commitment and a chance. If you are ready to contribute technical expertise, innovative thinking, ambition and a high level of persistence, at Sartorius you will find the chance to make an impact.

### Grow With Us

Sartorius is a growing global life science company, focusing on dynamic markets that will continue to grow for decades. This creates opportunities for those who seize them. We entrust our people with responsibility and in return expect the courage to take risks to move forward. Depending on personal goals and changing priorities in life, you can grow your career in diverse ways – from technical to geographical.

### Add the Missing View

Sartorius offers the perfect size and culture for ambitious doers and diverse minds: every voice is heard. Your perspective could be the missing piece to the next game-changing discovery. Yet we believe that challenges are best mastered together: When experienced specialists fuse their expertise and views. That is why we foster exchange across fields, borders, cultures, and social backgrounds.



**Valérie Morillon**

Leads the virtual sales team for lab products and services in the EMEA region

“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 the bioprocess solutions division in Göttingen, Germany

“I started as a clerk at Sartorius in 2013 – today, I am responsible for bio-process operations at the Göttingen site with almost 1,400 colleagues. This journey 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 the manufacturing of products that are used in the fight against diseases worldwide.”



**Shelly Ren**

Responsible for the bioprocess marketing team in China

“Sartorius is a great company for life science talents to bring innovative technologies to the biopharmaceutical industry. I deeply appreciate the open, flexible, encouraging working environment the company offers, helping me to make my career dream a reality.”



Read more career stories on the Sartorius blog.  
[www.sartorius.com/blog](http://www.sartorius.com/blog)

# Benefit From Attractive Working Conditions

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.

# Advance Your Career

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.



Join Our  
Global Team  
and Grow  
With Us.

Visit [www.sartorius.com/careers](http://www.sartorius.com/careers)  
or find us on LinkedIn.



Simplifying Progress