

Any questions?

If you have questions concerning the scholarship or your application, our HR experts will be happy to help.

Sartorius Scholarship, Germany



Daniela Müller
Human Resources Management
Sartorius Corporate Administration GmbH
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49.551.308.2408
daniela.mueller@sartorius.com

Sartorius Scholarship, India



Bharathi Krishnan
Human Resources Management
Sartorius Stedim India Pvt. Ltd.
No. 69/2 & 69/3, NH 48
Jakkasandra, Nelamangala
Bangalore – 562 123
bharathi.krishnan@sartorius-stedim.com

Sartorius Corporate Administration GmbH
Corporate Communications
37079 Göttingen, Germany

To learn more about Sartorius as an employer,
visit us at our website www.sartorius.com/career.





Sartorius Scholarship

for Masters Students

From Campus to Company

Are you ready to move on to a master's program or have you started one while being eager to get a foothold in the working world at the same time? The Sartorius Scholarship is a unique opportunity to enhance your education and career opportunities with our financial, professional and personal support.

Our scholarship is designed to match your individual skills, interests and needs to unfold your full potential. Learn more about who we look for, what we offer and how to apply.



"When searching for scholarships, I was immediately impressed by Sartorius' offer. I appreciate the regular exchange with my mentor, experts and other scholars. I really enjoyed working on internationally oriented teams and projects. At the same time the family-like culture made me feel immediately integrated as part of the organization."

Johanna Wilzmann completed her bachelor's dual study program at a global healthcare supplier. After two years of professional experience, she enrolled in a master's program in business studies and became a Sartorius scholar. Her internship in the company helped her gain insights into the implementation, harmonization and optimization of international sales processes. She stayed in regular contact with her mentor at Sartorius and attended several trainings. After she finished her master's degree, Johanna Wilzmann has been hired within the Sales department.



Who We Look For

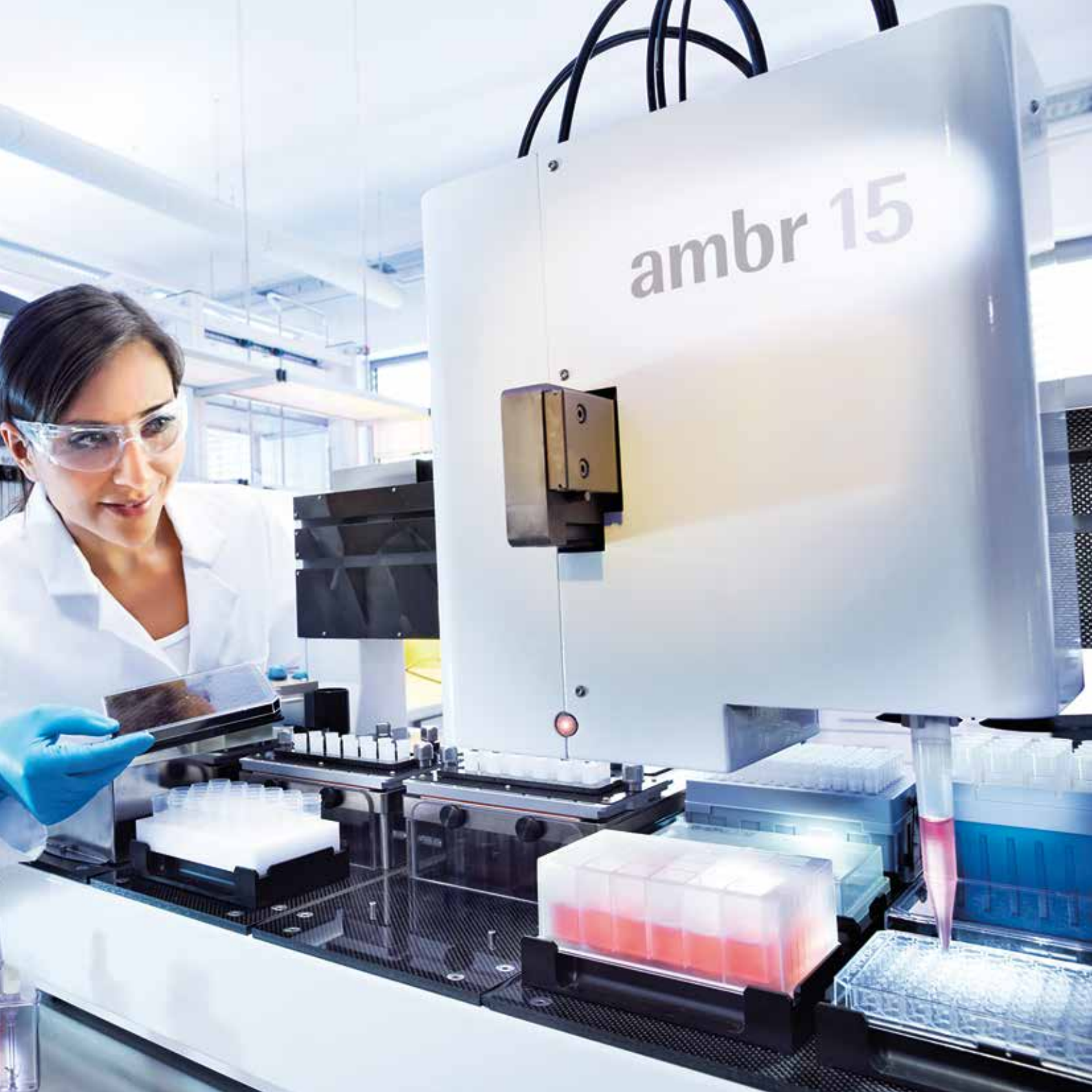
- » You have earned a bachelor's degree in natural sciences, engineering, business informatics or business studies with good or excellent grades
- » You are enrolled in a master's program in one of these fields at a German or Indian university, or intend to do so
- » You would like to gain professional experience at an internationally leading provider of pharmaceutical and laboratory equipment
- » Your English skills are strong; experience in working or studying abroad is a plus
- » You see challenges as opportunities, enjoy taking responsibility and act proactively
- » You identify with our corporate values of sustainability, openness and enjoyment

What We Offer

We build the scholarship around your scientific interests and career goals. For example, if you opt for an internship in R&D, you will take part in advancing new technologies and designing product prototypes. Working in Sales or Marketing departments will give you the chance to learn about product launches, marketing concepts and sales techniques in global markets. By working together with related units, you will gain insights in other departments as well. During the Sartorius Scholarship program you can write your thesis, and your results can even become part of a new Sartorius product. Get the best of both worlds: deepen your knowledge and put what you learned into practice.

- » A monthly financial grant during the master's program
- » An internship or master's thesis at Sartorius
- » A personal mentor for one-on-one consultation regarding professional and personal development
- » Training seminars, e.g. in project management and communication
- » Networking with other scholarship holders and experts
- » Good prospects for starting your career with us









How To Apply

Our application process will give you the opportunity to present your skills. Your online application needs to be written in English and should include the following:

- » A letter of motivation
- » A résumé (CV)
- » A letter of reference from your professor or university lecturer
- » A copy of your bachelor's degree certificate or of a comparable diploma
- » Job references, testimonials or letters of recommendation from your employer, if any, as well as any certificates of other training or completed education

Our group assessment center will give us both the opportunity to get to know each other personally. Here you will be asked to introduce yourself to our selection panel in an interview, make a presentation on a pre-announced topic, and take part in a role-playing situation together with other candidates. Once we have evaluated all results, we will make our decision and inform you promptly. We look forward to receiving your application online: <https://www.sartorius.com/sartorius/en/EUR/company/careers/students/sartorius-scholarship>

Who We Are

Sartorius is a leading international pharmaceutical and laboratory equipment supplier. With innovative products and high-quality services our two divisions, Bioprocess Solutions and Lab Products & Services, help customers all over the globe to implement complex and quality-critical processes in biopharmaceutical production and laboratory environments in a time- and cost-efficient way.





Enabling Medical Progress

A growing number of medications are biopharmaceuticals. These are produced using living cells in complex, lengthy and expensive procedures. To enable the still young biopharma sector to continue growing, it must find ways to manufacture such drugs more efficiently. The Bioprocess Solutions division provides the essential products and technologies to accomplish this.

From cell cultivation to harvesting, purification, final formulation and filling, Sartorius covers nearly all steps in the production process of a biopharmaceutical. Our comprehensive bioreactor portfolio makes us the worldwide leader in upstream processing. In the downstream sector, we are advancing technology by developing innovative and enhanced single-use technologies, such as membrane chromatography, body feed filtration for antibody harvesting, and freeze and thaw systems.



Host cell with modified DNA in a cell culture medium.





"I was surprised at how flexible Sartorius was in customizing its scholarship. As my master's program was very time-consuming, I couldn't afford to work several months as an intern. So my mentor at Sartorius organized one-day workshops every month, where I could talk to him and to experts from other R&D departments, present my ideas and collect useful information for my thesis."

Svenja Guillaume studied biotechnology and wrote her master's thesis at Sartorius. During her scholarship, she was based at the R&D Department for Membrane Technologies in Göttingen. Afterwards she finished off her doctoral thesis on the development of sterile filters for final filling applications in the biopharmaceutical industry and simultaneously assumed as an in-house application specialist for filtration technologies.

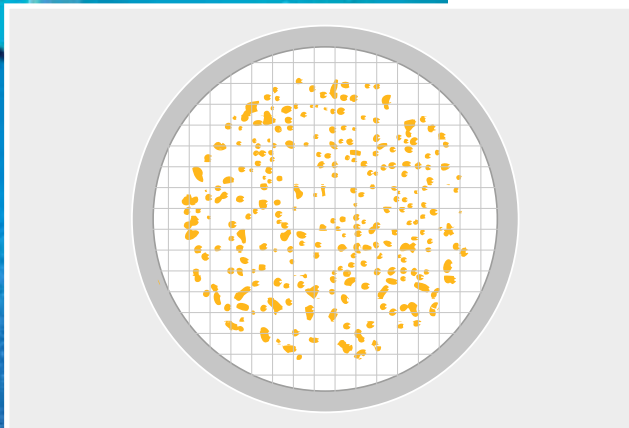




Making Lab Life Easier

Lab work is complex and demanding: Despite repetitive analytical routines, staff in quality assurance laboratories and public research institutes must perform each step in a highly concentrated and careful way for reliable and accurate results. Lab Products & Services, Sartorius' other division, helps them excel: with products that minimize human error, simplify workflows and reduce physical workloads.

Sartorius is well known for its premium lab instruments, such as balances, pipettes and lab water systems and among the leading suppliers of these products worldwide. It also offers a broad portfolio of consumables as well as services that cover the entire lifecycle of an instrument.



Nutrient pad for TTC assays



At Home Worldwide

With its own production facilities, sales subsidiaries and local commercial agencies Sartorius serves customers in more than 110 countries. They stand to benefit from this close proximity thanks to fast delivery, accelerated order processing and in-person service.

Sartorius focuses its business on three core markets: North America, Europe and Asia.

THE AMERICAS

North America, the world's largest pharmaceutical and laboratory market, where many key accounts are headquartered, is served from the main sales office in Bohemia, New York. The company's key manufacturing facility for the USA is its plant in Yauco, Puerto Rico, which supplies membrane filters and sterile single-use bags.



Bohemia

Yauco





Göttingen

EMEA

Based on its historical roots, Sartorius has the densest network of sites in Europe. The German university city of Göttingen has been home to the company's headquarters, its largest manufacturing and R&D site, up to this very day. This is where Sartorius develops and manufactures its two core technology products, membrane filters and laboratory balances.



Beijing

ASIA | PACIFIC

In the dynamically growing Asian market, Sartorius operates out of its sales hub in Shanghai. Its key production centers are in Beijing, China, and Bangalore, India, at which Sartorius manufactures stainless steel bioreactors, single-use bags and laboratory balances, among other products.

Bangalore



Working at Sartorius

Over 7,500 people worldwide contribute to the success of Sartorius with their ideas, expert knowledge and passion to excel. The company has been attracting employees who appreciate freedom, would like to grow personally and professionally and who put their hearts into their work.



"The interdisciplinary project work was lots of fun for me. Thanks to my internship in Business Development, I gained thorough insight into the company and learned a lot about successful project management. My participation in trainings, the research project in India and regular exchange with my mentor optimally rounded off my scholarship."

Markus Loidol earned his diploma in bioprocess and environmental process engineering, and completed two internships at Sartorius. He spent three months at the Research and Development Department in Bangalore, India, where he wrote his diploma thesis on the validation of bioreactors. In Göttingen, he supported the Business Development Department for four months. His work at this unit focused on the project involving the integration of an acquisition. During his time as an intern, Markus Loidol lived in the apartment shared by Sartorius Scholarship work-study students.



Make an Impact

Sartorius has all it takes to lead and shape markets internationally: cutting-edge technologies, a strong brand and the financial resources. We want to make a real difference in our industries and welcome people who share this ambition. Person-to-person communication, straightforward decision-making and the freedom to make your own footprint: this is the way to work and succeed at Sartorius.

Grow with Us

The biotech industry and laboratories around the world, our major markets, are fast-evolving sectors that we believe will continue to grow for decades. Supplying these markets calls for professionals who are at the same time highly quality-minded, creative and customer-focused. Depending on your personal goals and ambitions, you can grow your career in a variety of fields and geographies while staying with Sartorius.

Strong Individuals, Great Team

Our customers associate Sartorius with quality products and innovative solutions. This reputation is built on experienced professionals who are top-ranked in their fields, yet great team players at the same time. We believe in creating success together.

Strong Roots and Values

Founded in 1870 as a spin-off of the University of Göttingen in Germany, and having been a publicly-traded company for 25 years, Sartorius still offers a personal, family style culture. At our company, everyone is important. Sartorius' employees are united across the world by three common values: sustainability, openness and enjoyment.

Sartorius Scholarship

für Studierende in Deutschland

Das Sartorius Scholarship richtet sich an herausragende Master-studierende der Natur- und Ingenieurwissenschaften sowie der Wirtschaftsinformatik und der Betriebswirtschaftslehre, die ihre wissenschaftlichen Kenntnisse in der Praxis anwenden und

vertiefen möchten. Sartorius fördert seine Stipendiaten fachlich, finanziell und persönlich durch Weiterbildungsangebote und eine intensive Betreuung.

Alle Leistungen im Überblick

- » Monatliches Stipendium in Höhe von 800 Euro bis zum Studienabschluss
- » Praktikum oder Werkstudententätigkeit bei Sartorius in Göttingen oder Guxhagen
- » Die Möglichkeit, seine Abschlussarbeit bei Sartorius zu schreiben
- » Kostenloses Zimmer in einer der Sartorius Praktikanten-WGs oder Mietzuschuss und Unterstützung bei der Wohnungssuche
- » Weiterbildungsmöglichkeit durch Trainings und Seminare
- » Persönliche Betreuung durch Mentoren-Programm und Austausch innerhalb des Stipendiaten- und Expertennetzwerks
- » Perspektive einer Folgebeschäftigung

Mehr zum Sartorius Scholarship, Auswahlkriterien und Bewerbungsverfahren erfahren Sie in der Broschüre. Oder besuchen Sie unsere Homepage, um sich ein Bild von Sartorius als Arbeitgeber zu machen.

www.sartorius.de/karriere

