

Sartorius Academia Innovation Award 2026

What would you do with an Ambr® 15 Cell Culture?



About the Sartorius Academia Innovation Award

The Sartorius Academia Innovation Award recognizes outstanding academic initiatives that advance bioprocess education and foster innovative research within academic institutions. The award includes a five year loan of an Ambr® 15 Cell Culture system from the Sartorius demo pool*, providing access to cutting edge technology that enables automated bioprocess development and enhances practical training opportunity. The participant ranked second shall receive a voucher equal to 60%, and the participant ranked third shall receive a voucher equal to 40%, each to be applied toward the purchase of an Ambr® 15 Cell Culture system.

Why Apply? Unlock New Possibilities with the Ambr® 15

The Ambr® 15 Cell Culture is more than a bioreactor system, it is a platform that can transform how academic institutions conduct research, train students, and accelerate innovation. With HT automated operation and DOE-MODDE capabilities, researchers can rapidly evaluate multiple process conditions, cell cultures, media formulations, and feeding strategies while using only a fraction of the resources required by shake flasks and traditional bioreactors.

Whether your vision is to develop the next generation of bioprocesses, explore novel therapeutic modalities, create innovative teaching programs, or establish a center of excellence in bioprocess development, the Ambr® 15 Cell Culture platform provides the technology foundation to turn ambitious ideas into reality.

*The EMEA Demo Pool is a centralized portfolio of demonstration equipment that enables customer to evaluate Sartorius solutions before making a purchasing decision

Sartorius Academia Innovation Award

Sartorius Stedim Biotech GmbH
August-Spindler-Straße 11
37079 Göttingen
Germany



For more information get in contact with us
or visit our website:

 EventsEMEA@sartorius.com

 [Sartoris Academia Innovation Award](#)

Who can participate?

Research driven academic institutions in the life sciences in the EMEA, particularly those focused on biopharma and proposing innovative ideas for an Ambr® 15 Cell Culture application, are eligible to participate. Entries may be submitted by individuals or teams, in English only, via the official registration form. Multiple submissions are not allowed.

Registration opens on **July 20, 2026**, and the submission deadline is **September 27, 2026**.

What do you need to apply?

You will need an innovative concept for how you would use the Ambr® 15 Cell Culture system, clearly highlighting the added value for your research. The registration form includes a mix of structured and open questions. Please take the time to complete all sections carefully and ensure you submit the most compelling and innovative application. Joint submissions with other teams are encouraged and will be viewed positively in the overall evaluation.

How are the entries evaluated?

An independent jury of experts from diverse fields will decide on the award and select the winner. Three criteria are key for their decision: degree of innovation, disruptive potential, and relevance to an existing market challenge.

The decisions of the jury are final and not subject to appeal.

Award Ceremony

The winners will be announced on **October 5, 2026** and awarded on October 29, 2026, as part of the Global Academia Summit at the Sartorius Campus in Göttingen.

Legal Notices

Participants grant Sartorius the non-exclusive right to store their contribution for one year after the submission. Submissions are valid without a signature. By submitting their entries, participants accept the terms of the competition and assure that their contribution is free from third-party rights. Those awarded with the loan of the Ambr® 15 Cell Culture system commit to attending the awards ceremony on October 29, 2026, in Göttingen.

Sartorius in Brief

Sartorius is a leading international partner of life sciences research and the biopharmaceutical industry. With innovative laboratory instruments and consumables, the Group's Lab Products & Services division focuses on laboratories performing research and quality control at pharmaceutical and biopharmaceutical companies as well as academic research institutes. The Bioprocess Solutions division, with its broad product portfolio focusing on single-use solutions, helps customers manufacture biotech medications, vaccines, and cell and gene therapies safely, rapidly, and economically. The company, based in Göttingen, Germany, has a strong global reach with around 60 production and sales sites worldwide. Sartorius regularly expands its portfolio through the acquisition of complementary technologies. In 2025, the company generated sales revenue of around 3.5 billion euros. Currently, more than 14,000 employees are working for customers around the globe.

Sartorius Academia Innovation Award

Sartorius Stedim Biotech GmbH
August-Spindler-Straße 11
37079 Göttingen
Germany



For more information get in contact with us
or visit our website:

 EventsEMEA@sartorius.com

 [Sartoris Academia Innovation Award](#)