

Apply now

# LifeScienceXplained 2026

## Sartorius Award for New Communication

### About LifeScienceXplained

The life sciences have the potential to change our lives. Medical knowledge is expanding rapidly worldwide. New types of therapy offer hope to millions of people; with innovations such as gene and cell therapies, diseases that have not been treatable until now can become curable.

The LifeScienceXplained | Sartorius Award for New Communication aims to promote the exchange between science and society on current topics in the life sciences and to foster an understanding of the complex concepts. It is endowed with 15,000 euros in prize money.

### Who can participate?

Communicators who use creative formats to make complex life science content easy and understandable can apply now.

Entries can be submitted by individuals or teams in various formats, including videos, podcasts, science slams, texts, and comics, in either German or English. Submissions must be specific thematic contributions, as entire social media channels are not eligible for evaluation. Multiple entries are permitted. The date of first publication must be between August 1, 2025 and August 1, 2026, and submitted at [www.sartorius.com/lsx](http://www.sartorius.com/lsx). The deadline for submission is August 12, 2026.

### How are the entries evaluated?

An independent jury of experts from diverse fields will decide on the award. They will nominate three finalists and select the winner. Three criteria are key for their decision: verifiability, comprehensibility, and creativity.

### Award Ceremony

The winner will be announced at the award ceremony on November 5, 2026 as part of the Göttinger Literaturherbst together with the NDR Sachbuchpreis at the Sartorius Campus in Göttingen.

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



### LifeScienceXplained 2026

Sartorius Award for New Communication

Sartorius Corporate Administration GmbH  
Corporate Communications  
Otto-Brenner-Straße 20  
37079 Göttingen

Lisa Kottas  
[LifeScienceXplained@sartorius.com](mailto:LifeScienceXplained@sartorius.com)  
[www.sartorius.com/lsx](http://www.sartorius.com/lsx)

# SARTORIUS

## Legal Notices

Participants grant Sartorius the non-exclusive right to use their contribution free of charge for one year after the publication of the shortlist in the name of the LifeScienceXplained Award and to make it publicly available worldwide. Entries may be edited as needed and excerpts may be used. 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 selected for the shortlist commit to being available for a video production starting at the end of September and to attending the awards ceremony on November 5, 2026, in Göttingen. The winner agrees to be a member of the jury the following year and to attend the jury meeting in person or virtually.

## 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.



## LifeScienceXplained 2026

Sartorius Award for New Communication

Sartorius Corporate Administration GmbH  
Corporate Communications  
Otto-Brenner-Straße 20  
37079 Göttingen

Lisa Kottas  
LifeScienceXplained@sartorius.com  
www.sartorius.com/lxs

# SARTORIUS