

Call for tenders

Apply now

LifeScienceXplained 2024

Sartorius Award for New Communication

- **LifeScienceXplained | Sartorius Award for New Communication honors outstanding contributions about the life sciences**
- **Comprehensible, explanatory formats such as videos, podcasts, texts, or illustrations are eligible**
- **In its fourth year, it is now also open for entries in English language and endowed with €15,000 in prize money**
- **The deadline for entries is August 4, 2024**

Life sciences and related technologies have the potential to transform lives and already today bring hope of recovery to millions of people. With innovations such as gene and cell therapies, previously incurable diseases like cancer or dementia might now be treated or even cured. The LifeScienceXplained | Sartorius Award for New Communication aims to advance the dialogue between science and society on current topics in the life sciences and to promote understanding.

Communicators who use creative formats to make complex life science content understandable can now apply for this year's LifeScienceXplained | Sartorius Award for New Communication. In its fourth year, it is now also open for entries in English language and endowed with €15,000 in prize money.

Entries can be submitted by individuals or teams; all formats, from videos, to podcasts, to texts, to comics, in German or English, are welcome. Entries must be published between August 1, 2023, and August 1, 2024, and submitted at www.sartorius.com/lSX. The deadline is August 4, 2024.

An independent jury of experts from various fields will decide on the award. They will nominate a shortlist and select the winning entry. Three criteria are key for their decision: verifiability, comprehensibility, and creativity. The winner will be announced at the award ceremony on November 7, 2024, along with the jury's reasoning.

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. If selected to the shortlist of the top five contributions, they commit to be available for a video production in calendar week 39 and to attend the award ceremony on November 7, 2024, in Göttingen.



LifeScienceXplained 2024

Sartorius Award for New Communication

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

Mihaela Mitrovic
LifeScienceXplained@sartorius.com
www.sartorius.com

SARTORIUS