

Data protection information for the LifeScienceXplained | Sartorius Prize for New Communication competition

In the following data protection information, Sartorius Corporate Administration GmbH ("Sartorius") informs you about the processing of your personal data when you participate in our competition LifeScienceXplained | Sartorius Prize for New Communication.

A. General data protection information

1. Name and contact details of the controller

The controller is:

Sartorius Corporate Administration GmbH
Otto-Brenner-Str. 20,
37079 Göttingen, Germany info@sartorius.com

2. Contact details of the data protection officer

The data protection officer can be contacted as follows:

Sartorius Corporate Administration GmbH
Data Protection Officer
Otto-Brenner-Str. 20,
37079 Göttingen, Germany dataprotection@sartorius.com

3. Scope of the processing of personal data

When you take part in the competition, we process the data and information you provide in the registration form. This includes the data required for participation, such as

- Personal identification data (e.g. surname, first name, title, date of birth);
- Contact details (postal address, telephone number, e-mail address);
- Photos, film and video recordings (e.g. the video recording sent for participation in the competition), blog, slam, podcast, article or book.
- Curriculum vitae

as well as any data and information voluntarily provided by you during participation.

4. Purpose of the processing of personal data

Your personal data will be processed for the purpose of running the competition, in particular to determine and notify the winners.

We may subsequently collect and process additional data, e.g. your bank details, for the purpose of sending and delivering prizes.

5. Legal basis for the processing

The legal basis for the processing is the fulfillment of the contractual obligation arising from participation in the competition (Art. 6 para. 1 sentence 1 lit. b GDPR)

6. Data erasure and storage duration

The participant's data will be stored for the duration of the competition and as long as the processing of your data is necessary for the establishment of the legal relationship with the participant, the subsequent implementation and processing of the competition (in particular selection and evaluation of the entries, as well as information of the finalists selected on the shortlist, invitation to the award ceremony or provision of the prize). In principle, the data of participants not taking part in the final will be deleted no later than 6 months after the end of the competition. The data of participants in the final will be stored until consent is withdrawn, but will be deleted after 2 years at the latest. The prize winner's data may be retained for longer, e.g. to answer queries about the prize or to fulfill the prize performance. In addition, Sartorius stores the personal data collected in the context of competitions to the extent and for as long as this is necessary to fulfill contractual or legal obligations. Sartorius processes personal data to fulfill commercial or tax law obligations to provide evidence and retain data. The periods for storage and retention stipulated therein documentation is six years in accordance with commercial law requirements pursuant to Section 257 HGB and up to ten years due to tax requirements pursuant to Section 147 AO. The periods begin at the end of the calendar year in which the document was created.

7. Categories of recipients

7.1 Disclosure to third parties.

In addition, we will pass on your data to third parties if this is permitted by legal or contractual provisions and/or if you have given your consent. Under this condition, the data may be passed on to the following categories of recipients: Public bodies and institutions (e.g. tax authority, data protection supervisory authority) to process official inquiries, insofar as this is in your interest or we are legally obliged to do so. The legal basis is Art. 6 para. 1 sentence 1 lit. f or c GDPR.

7.2 Processor (Art. 28 GDPR).

Service providers commissioned by Sartorius to support Sartorius in the execution of the business relationship will have access to the data. If other categories of processors are also used, please refer to the respective special data protection notices.

8. Transfer to third countries

Data is not transferred to third countries (countries outside the EU or the European Economic Area - EEA).

9. Obligation to provide personal data

Unless otherwise stipulated in the Special Data Protection Notice, you are neither legally nor contractually obliged to provide your data.

10. Rights of data subjects

In accordance with Art. 15 GDPR, you have the right to obtain information about the personal data stored about you. If incorrect personal data has been processed, you have the right to rectification in accordance with Art. 16 GDPR. If the legal requirements are met, you can request the erasure or restriction of processing and object to data processing (Art. 17, 18 and 21 GDPR). According to Art. 20 GDPR, you can assert the right to data portability for data that is processed automatically on the basis of your consent or a contract with you.

Information about your right to object in accordance with Art. 21 GDPR

You have the right to object, on grounds relating to your particular situation, at any time to processing of personal data concerning you which is based on point (f) of Article 6(1) GDPR (data processing on the basis of a balancing of interests).

If you object, Sartorius will no longer process your personal data unless Sartorius can demonstrate compelling legitimate grounds for the processing which override your interests, rights and freedoms or for the establishment, exercise or defense of legal claims.

Right to withdraw your consent in accordance with Art. 7 (3) GDPR

If you have consented to the processing of your personal data, you have the right to withdraw your consent at any time. The withdrawal of consent does not affect the lawfulness of processing based on your consent before its withdrawal. The revocation can be made informally. You can contact us using the contact details provided in section A point 2.

You can assert your rights against Sartorius using the contact details given under point 1.

If you are of the opinion that data processing violates data protection law, you have the right to lodge a complaint with a data protection supervisory authority of your choice. (Art. 77 GDPR).

Of course, you can also lodge a complaint with the Sartorius data protection officer at any time (dataprotection@sartorius.com).

B. Special data protection information: Participation in the LifeScienceXplained | Sartorius Prize for New Communication competition

You have the opportunity to take part in the LifeScienceXplained | Sartorius Prize for New Communication competition by making complex life science content simple and easy to understand in your blog, video, slam, podcast, article or book.

Sartorius processes the relevant personal data that you provide to us by participating in our LifeScienceXplained competition in order to carry out the competition (e.g. to determine whether you are eligible to participate in the competition). For this purpose, we process the following personal data from our participants Personal identification data (e.g. surname, first name, title, date of birth); contact details (postal address, telephone number, e-mail address); photos, film and video recordings (e.g. the video recording sent to participate in the competition), blog, slam, podcast, article or book.

The legal basis is Art. 6 para. 1 sentence 1 lit. b GDPR. Data processing is necessary for the performance of the contract for participation in the LifeScienceXplained | Sartorius Prize for New Communication competition.

You can only take part in the competition if you provide us with the necessary data.

C. Special data protection information: Determination and notification of the winner and payment of the prize

We will contact the prize winner and the participants in the LifeScienceXplained competition using the details provided when registering as a participant. So that we can pay the prize to the winner, we ask them to provide us with their bank details.

Sartorius processes the relevant personal data that you provide to us for participation in a competition in order to carry out the competition (e.g. to determine and notify the prize winner and the participants in the award ceremony, as well as to pay out or deliver the prize). We process the following personal data: Personal identification data (e.g. surname, first name, title); contact data (telephone number, e-mail address); financial identification data (e.g. bank details).

The legal basis for data processing is Art. 6 para. 1 sentence 1 lit. b GDPR. Data processing is necessary for the performance of the competition contract.

Without your data, we will not be able to contact you as a prize winner or as a participant in the award ceremony and pay out the prize to the winner.

D. Special data protection information: Publication of the finalists' entries

We would like to publish the finalists' entries as a token of our appreciation, as part of the event, in print and online media and on our website www.sartorius.com.

For this purpose, we contact the prize winner and the participants in the award ceremony to ask them separately for their consent to the publication of the submitted entry. For this purpose, we process the following personal data from our participants Personal identification data (e.g. surname, first name, title) contact details (telephone number, e-mail address); photos, film and video recordings (e.g.

SARTORIUS

the video recording sent), blog, slam, podcast, article, infographic, comic or book. The legal basis is your express consent in accordance with Art. 6 para. 1 sentence 1 lit. a GDPR.

You can revoke your consent at any time with effect for the future, see Section A point 10.