

# Information on Data Protection for Event Participants of the Capital Markets Day of Sartorius Aktiengesellschaft

Data Protection is a top priority at Sartorius (hereinafter referred to as “we”), which includes transparent information about the use of your personal data. Therefore, we would like to provide you with an overview of how your personal data is processed by Sartorius, and your rights under Data Protection law. It goes without saying that we only process personal data in accordance with the applicable legal provisions on the protection of personal data and data security.

## Who is responsible for data processing and who is the data protection officer?

The responsible controller for data processing is:

### **Sartorius AG**

Otto-Brenner-Strasse 20

37079 Goettingen

Germany

Phone: +49 (0)551 308-0

E-mail: [info@sartorius.com](mailto:info@sartorius.com)

You can reach our Data Protection Officer at:

### **Data Protection Officer**

Otto-Brenner-Strasse 20

37079 Goettingen

Phone: +49 (0)551 308-0

E-mail: [dataprotection@sartorius.com](mailto:dataprotection@sartorius.com)

### **What categories of data do we use and where do they come from?**

We collect personal data directly from event participants as part of the online registration process via the event website.

The categories of personal data processed from event participants include:

- Contact information (first and last name, address, telephone number, mobile phone number, fax number, e-mail address);
- Professional data (position/role);
- Interests;
- Video recording;
- Spoken contributions to the floor during the event. Can also be made anonymously.

The provision of certain data is necessary in order to be able to properly prepare, conduct, and follow up on the event. Without the provision of certain data, participation in the event is not possible or could be impaired if the processing of certain data is not consented to or data is not given. This expressly does not apply to the processing of video recordings and spoken contributions.

### **For what purposes and on what legal basis is the data processed?**

We process your personal data in compliance with the provisions of the EU General Data Protection Regulation (EU-GDPR) and all other applicable laws.

The data processing serves the following purposes:

- Communication with event participants regarding participation;
- Responding to inquiries regarding the event such as venue and times;
- Preparation, implementation, and follow-up of the event;
- Safeguarding our legitimate interests, namely reporting and providing information about Sartorius events, company-related events, and for contemporary historical documentation;
- Public relations.

The processing of the personal data categories above is necessary to achieve the following purposes. The legal basis for data processing is—unless expressly stated otherwise – our legitimate interest (Art. 6 (1) (f) EU-GDPR). The legitimate interest is to ensure the smooth running of the event. In individual

cases, we obtain explicit consent pursuant to Art. (1) (a) EU-GDPR (e.g., in the case of video and image recordings and verbal contributions). If we wish to process the personal data of event participants for a purpose not mentioned above, they will be informed in advance.

#### **To whom is the personal data disclosed?**

For the purposes mentioned above Sartorius may transfer or disclose your personal data to:

- Other companies of the Sartorius Group or third parties in the context of your event participation;
- Third parties which provide IT services to Sartorius, and which process such data only for the purpose of such services (e.g., hosting or IT maintenance and support services); and/or
- Third parties in connection with complying with legal obligations or establishing, exercising or defending rights or claims (e.g., for court and arbitration proceedings, to law enforcement authorities and regulators, to attorneys and consultants).

Sometimes the recipients to whom Sartorius transfers or discloses your personal data are located in countries in which applicable laws do not offer the same level of data protection as the laws of your home country.

In such cases, if required by applicable law, Sartorius takes measures to implement appropriate and suitable safeguards for the protection of your personal data. Thus, Sartorius transfers personal data to external recipients in third countries only if the transfer is subject to appropriate safeguards, e.g., EU Standard Contractual Clauses or Binding Corporate Rules.

#### **How long will the personal data be stored?**

If no explicit storage period is specified at the time of collection (e.g., as part of a declaration of consent), the personal data of event participants will be deleted insofar as it is no longer required to fulfill the purpose for which it was stored, unless legal retention obligations (e.g. retention obligations under commercial and tax law) prevent deletion. In this case, the data will only be processed to a limited extent. In the context of the publication of the event, non-anonymized questions and spoken contributions will be edited out.

### What Data Protection rights do I have?

Under applicable data protection law, you may have the right to:

- Obtain confirmation from Sartorius as to whether or not personal data concerning you is being processed, and where that is the case, access to the personal data;
- Obtain from Sartorius the rectification of inaccurate personal data concerning you;
- Obtain from Sartorius the erasure of your personal data;
- Obtain from Sartorius restriction of processing regarding your personal data;
- Data portability concerning personal data, which you actively provided;
- Object, on grounds relating to your particular situation, to processing of personal data concerning you; and
- Complain to a supervisory authority (e.g., to the State Commissioner for Data Protection of Lower Saxony, Prinzenstrasse 5, 30159 Hanover, Germany).

In case you declared your consent for the processing of certain personal data by Sartorius, you have the right to withdraw the consent at any time with effect for the future, i.e., the withdrawal of the consent does not affect the lawfulness of processing based on the consent before its withdrawal. If the consent is withdrawn, Sartorius may only further process the personal data where there is another legal basis for the processing.