

# Sartorius AR Assistance Privacy Notice

This Privacy Notice provides information on the processing of your personal data in connection with your use of the Sartorius AR Assistance Application.

**Personal data** include any information related to an identified or identifiable natural person. An identifiable natural person is one who can be identified, directly or indirectly, in particular by reference to an identifier such as a name, an identification number, location data, an online identifier or to one or more factors specific to the physical, physiological, genetic, mental, economic, cultural or social identity of that natural person.

**Processing** means any operation or set of operations which is performed on personal data or on sets of personal data, whether or not by automated means, such as collection, recording, organization, structuring, storage, adaptation or alteration, retrieval, consultation, use, disclosure by transmission, dissemination or otherwise making available, alignment or combination, restriction, erasure or destruction.

## 1. Identity and contact details of the controller and the Data Protection Officer

Sartorius Stedim Biotech GmbH, August-Spindler-Str. 11, 39079 Göttingen, Germany ("**Sartorius**", "**we**", "**our**" or "**us**"), as controller of personal data, is responsible for the processing of your personal data in connection with your use of the Sartorius AR Assistance Application ("**App**").

Our Data Protection Officer and the Sartorius Group Data Protection Organization are available to you as a contact for all data protection-related matters and for exercising your rights. They may be contacted at the above address or at [dataprotection@sartorius.com](mailto:dataprotection@sartorius.com).

## 2. Information collected when downloading the App

Before you can download and install this App, you may need to enter into a usage agreement with an App Store operator (e.g. Google, Apple) to access its portal (e.g. Google Play, Apple App Store - "App Stores"). The App Store operator collects and processes data in connection with the use of the App Store, such as user name, email address and individual device identification number as the responsible Controller. We process the data only to the extent necessary to download the App to your device. We are

not party to the usage agreement with the App Store operator and have no influence on its data processing. In this respect, the privacy policy of the respective App store operator applies.

### **3. Processing of personal data**

The App runs offline. When using the App, Sartorius does not process any information, thus also no categories of personal data.

Sartorius may only process your personal data, if you contact us. Those processing activities are not subject to this privacy notice and are addressed, for example, in our [Sartorius Website Privacy Notice](#).

### **4. External links**

This Privacy Notice only applies to this App, not to websites or applications offered and operated by third parties. The App may provide links to other websites or applications of third parties which may be of interest to you. After you click the link, we no longer have any influence on the collection, processing and utilization of any personal data that is transferred to third parties, as our control of the conduct of third parties is then naturally withdrawn. We are not responsible for privacy practices or the content of external websites or applications.

### **5. Permissions**

Certain functions of this App can only be used if the corresponding permissions are granted. Permissions are interfaces to the operating system of your end device, through which the App can access data stored on your end device. The App works with the following permissions:

- Camera access: to place the hologram on the right place

You can manage and disable permissions within the settings of your operating system. Please note that after deactivating a permission, you may no longer be able to use all of the App's features.

### **6. Transfer and disclosure of personal data**

As we don't process any data, Sartorius does also not transfer or disclose (your personal) data to service providers or other third parties.

### **7. Children**

To use this App, you must be at least 16 years old.

### **8. Your rights**

Under applicable data protection law, you may have specific rights in relation to your personal data. In particular, and subject to the statutory requirements, you may have the following data protection rights:

- **Right of access:** You have the right to obtain information on the processing of your personal data and to receive a copy of these data.
- **Right to rectification:** You have the right to request that we correct or complete your inadequate, incomplete or inaccurate personal data.
- **Right to erasure:** Under certain circumstances, you have the right to request that we delete your personal data.
- **Right to restriction of processing:** Under certain requirements, you may request us to restrict the processing of your personal data.
- **Right to data portability:** You have the right to receive your personal data in a structured, common, and machine-readable format and request that these data are transferred to another data controller, if applicable under the specific circumstances.
- **Right to object:** You might have the right to object to the processing of your personal data by us, in particular, if the processing of your personal data is based on (i) the necessity of the performance of a task in the public interest, or (ii) legitimate interests. We will then stop the processing of your personal data unless we remain legally authorized to do so.
- **Right to lodge a complaint:** You also have the right to lodge a complaint with a supervisory authority or other applicable privacy regulator about our processing of your data. This can be for example the data protection authority in your country of residence. A list of all data protection authorities in the European Union can be found [here](#).
- **Right to withdrawal:** If data processing is based on your consent, you have the right to withdraw your consent at any time and free of charge, with effect for the future, i.e., your withdrawal does not affect the lawfulness of the processing based on consent before its withdrawal.

Status of this Privacy Notice: August 2024