

Data protection information

on the processing of personal data in the context of the Sartorius Eureka Grant Program

Who is responsible for data processing and who can I contact?	Sartorius Singapore Pte. Ltd. 30 Pasir Panjang Road, #06-32, Mapletree Business City, Singapore 117440 You can reach our data protection officer at: lyekiat@asiacloud.com.sg +65 6471 6180
What do we process your data for (purpose of processing) and on what legal basis?	We process your personal data to administer the Sartorius Eureka Grant Program and to conduct the prize drawing. We process your data for administering and evaluating the Eureka Grant Program.
Categories of personal data	Full name, Job title, Organization or Institution, Department, Country, Work email address, Contact number, Project title, Signature, research proposal, project abstract and supporting information
Who receives my data?	Within the Sartorius Group the organisational units provided with your data are those that require them for the purpose of processing them. Your personal data may be provided to external processors contracted by the company for these same express purposes. These are firms providing services in the fields of IT services, telecommunications, printing services, etc.
How long will my data be stored?	We store your data for the duration of the Sartorius Eureka Grant Program and delete it as soon as the purpose for which it was collected no longer applies and there is no legal obligation to retain it.
Cross-border transfers	Some of the recipients of your data listed above may be located in countries outside of Singapore, where data protection laws may be less comprehensive. In such cases, we ensure that appropriate safeguards are in place to protect your personal data in accordance with our legal obligations.
What data protection rights do I have?	Every data subject has the right to information about the processing of their personal data, the right to rectification, the right to erasure, the right to restriction of processing, the right to object to processing and the right to data portability within the scope of the statutory provisions. There is also a right to lodge a complaint with a data protection supervisory authority. To assert your rights, please contact our data protection officer at lyekiat@asiacloud.com.sg. In addition to contacting our data protection organization, you also have the option of contacting the responsible Personal Data Protection Commission at any time.