



## Standard Terms and Conditions of the Sartorius Eureka Grant

that all information provided is true, accurate, complete, and not misleading.

### GENERAL TERMS AND CONDITIONS

#### 4) Selection Criteria

- a) Applications will be evaluated on the basis of the scientific merit, innovation, and potential impact of the proposed research, as well as the critical role of Sartorius BioAnalytical instruments in enabling or advancing the project outcomes.
- b) Additional preference may be given to proposals with a strong scientific rationale, clear objectives, a defined experimental plan, and a feasible timeline to project completion.
- c) Sartorius may take into account additional factors reasonably related to the purpose and integrity of the Grant Program, including compliance, feasibility, substantiation of statements made, and alignment with the eligibility requirements.

#### 5) Selection Decision; No Entitlement; Limitation of Challenge

- a) Sartorius shall have sole and reasonable discretion in evaluating applications and selecting the recipient.
- b) There is no legal entitlement to the grant, even if an applicant meets all eligibility criteria or receives a favorable evaluation in one or more categories.
- c) Sartorius is not obliged to disclose detailed scoring, internal deliberations, comparative assessments, or reasons for selecting or not selecting an applicant, except where disclosure is required by mandatory law.
- d) The selection decision of Sartorius is final. To the maximum extent permitted by applicable mandatory law, any legal recourse against the selection decision is excluded.

- e) For the avoidance of doubt, participation in the application process does not give rise to any claim for award, continuation of the Grant Program, reimbursement of costs, damages, or conclusion of any further agreement, except as expressly provided in these Terms and Conditions.

#### 6) Grant Amount and Permitted Use

- a) One selected applicant will receive a grant in the amount of EUR 5,000.

#### 1) Organizer and Purpose of the Program

- a) The Sartorius Eureka Grant 2026 (the "Grant Program") is organized by Sartorius Singapore Pte. Ltd. ("Sartorius").
- b) The purpose of the Grant Program is to support promising scientific research projects by awarding one grant in the amount of **EUR 5,000** to an eligible applicant for the funding of critical experiments and/or publication-related costs, in accordance with these Terms and Conditions.
- c) The Grant Program is a voluntary support initiative of Sartorius. It does not constitute a public funding scheme, prize competition governed by mandatory local consumer promotion rules, or a contractual offer to conclude a funding agreement with any applicant.

#### 2) Eligibility

- a) Applicants must be researchers affiliated with an academic institution, research institute, biotech company, or startup and must be a principal investigator, post-doctoral researcher, or scientist with institutional approval to conduct the proposed work.
- b) The Grant Program is open only to applicants located in Japan, India, South East Asia, Australia and New Zealand.
- c) Employees of Sartorius, its affiliates, and any persons directly involved in the administration or evaluation of the Grant Program may not apply.

#### 3) Application Process

- a) Applications must be submitted through the official Sartorius Eureka Grant application platform by **31 October 2026**.
- b) Incomplete, inaccurate, misleading, or late submissions will not be considered.
- c) Each applicant may submit only one application per research project. Multiple submissions for the same project may result in disqualification.
- d) By submitting an application, the applicant confirms

- b) The grant funds are intended solely to support research and publication-related expenses as specified by Sartorius, including, where applicable:
  - I) purchase of at least one reagent or consumable to be used with Sartorius BioAnalytical instruments (including the IncuCyte®, Octet® BLI, iQue® and CellCelector); and/or
  - II) Consumables to support the proposed study; and/or
  - III) article processing charges for a publication, if intended publication is shown to utilize such instruments in the relevant research work.
- c) The applicant must have access to the Sartorius BioAnalytical instrument used in the proposed research work.
- d) The grant must be used by **30 June, 2027**, unless Sartorius expressly agrees otherwise in writing.
- e) The recipient shall maintain reasonable records of the use of funds and provide supporting documentation upon request by Sartorius.

## 7) Recipient Obligations

- a) The recipient shall use the grant exclusively for the approved purpose and in compliance with all applicable laws, regulations, internal institutional rules, and ethical requirements.
- b) The recipient shall promptly notify Sartorius of any material change affecting eligibility, project feasibility, use of funds, institutional approvals, or compliance status.
- c) Upon reasonable request, the recipient shall provide a short update regarding the status of the funded project and the use of funds.

## 8) Publicity and Communications

- a) If the applicant has expressly agreed in the application form to future feature use by Sartorius, Sartorius and the recipient may separately agree on appropriate use of the recipient's name, institution, project title, and non-confidential project description in scientific communications, marketing materials, case studies, or event presentations.

- b) No obligation exists for the recipient to provide testimonials, endorsements, or promotional statements unless separately agreed in writing.
- c) Sartorius shall not publish confidential information of the applicant without appropriate permission, except where required by law.

## 9) Intellectual Property and Confidentiality

- a) All intellectual property rights in the application materials and the project shall remain with the applicant or the relevant rights holder.
- b) By submitting an application, the applicant grants Sartorius a non-exclusive, royalty-free right to review, internally circulate, and evaluate the submitted materials for the purposes of administering the Grant Program.
- c) Applicants shall not include third-party confidential information or personal data beyond what is reasonably necessary for the application unless lawfully disclosed and properly authorized.
- d) Unless expressly agreed otherwise in writing, the application shall not create any confidentiality obligations beyond those mandatorily imposed by applicable law.

## 10) Reservation of Rights; Modification, Suspension, Termination

- a) Sartorius reserves the right, at its reasonable discretion, to amend, suspend, postpone, or terminate the Grant Program, in whole or in part, at any time, in particular where necessary for legal, regulatory, operational, compliance, or integrity reasons.
- b) Sartorius further reserves the right to reject, withdraw, revoke, or reallocate the grant award if:
  - I) the applicant or recipient does not meet or no longer meets the eligibility requirements;
  - II) the application contains false, inaccurate, incomplete, or misleading information;
  - III) the applicant or recipient breaches these Terms and Conditions or applicable law; or
  - IV) the grant cannot be validly or lawfully

administered as intended.

- c) Sartorius may request repayment of all or part of the grant if the recipient obtained the grant through misrepresentation or used the funds in material breach of these Terms and Conditions.

## 11) Liability

- a) Sartorius shall not be liable for any costs, expenses, or losses incurred in connection with participation in the application process, preparation of application materials, or non-selection.
- b) To the maximum extent permitted by applicable law, Sartorius shall not be liable for indirect, incidental, special, consequential, or punitive damages arising out of or in connection with the Grant Program.
- c) Nothing in these Terms and Conditions shall exclude or limit liability to the extent such exclusion or limitation is prohibited by mandatory law, including liability for willful misconduct, fraud, or, where mandatorily applicable, death or personal injury caused by negligence.

## 12) Data Protection

- a) Personal data submitted in connection with the Grant Program shall be processed by Sartorius for the purposes of administering the Grant Program, evaluating applications, contacting applicants, documenting the award process, and meeting legal or compliance obligations.
- b) Further information regarding the processing of personal data, including the identity of the controller, categories of personal data, retention periods, and data subject rights, shall be provided in the applicable privacy notice referenced in the application platform.

## 13) Compliance with Laws and Institutional Rules

- a) The applicant is solely responsible for ensuring that participation in the Grant Program and acceptance and use of the grant are permitted under the applicant's institutional policies, funding rules, internal approval requirements, and applicable anti-corruption, sanctions, export control, and ethics rules.
- b) Sartorius may request confirmation of relevant institutional approvals where appropriate.

## 14) Governing Law and Jurisdiction

- a) These Terms and Conditions and any non-contractual obligations arising out of or in connection with them shall be governed by the laws of Singapore, excluding its conflict-of-laws rules and excluding the United Nations Convention on Contracts for the International Sale of Goods (CISG).
- b) To the extent legally permissible, the exclusive place of jurisdiction for all disputes arising out of or in connection with the Grant Program shall be Singapore.

## 15) Severability

- a) If any provision of these Terms and Conditions is or becomes invalid, illegal, or unenforceable, the validity and enforceability of the remaining provisions shall not be affected.
- b) Any invalid, illegal, or unenforceable provision shall, to the extent permitted by law, be deemed replaced by a valid and enforceable provision that most closely reflects the commercial and legal purpose of the original provision.

## 16) Acceptance of Terms

- a) By submitting an application, the applicant acknowledges and agrees to these Terms and Conditions.