

Sartorius Eureka Grant:

Supporting the Next Step in Scientific Discovery



About the Sartorius Eureka Grant

One experiment. One dataset. One breakthrough. The Sartorius Eureka Grant will award EUR 5,000 to help one researcher turn promising science into meaningful impact.

Who should apply?

Passionate researchers who are:

- Affiliated with academic institutions, research institutes, biotech companies or startups
- Principal investigators, postdoctoral researchers, scientists

Eligible regions:

Japan, India, Southeast Asia, Australia & New Zealand

How to apply?

Simply scan the QR code on the right to get started.

How applications will be evaluated

Applications will be assessed based on the scientific merit, innovation, and potential impact of the proposed research.

Eligible projects should be:

- Be at the data-generation stage
- Include a clear experimental plan utilizing at least one Sartorius BioAnalytical instrument
- Demonstrate a clear project completion plan by June 2027

Mark your calendar

- Application window: 1 July – 31 October 2026
- Recipient announcement: By 31 December 2026

Fuel your experiment.,
Apply for the Sartorius
Eureka Grant today:



Please review the **Terms and Conditions** on the application form before submission.

*For further inquiries, please
contact Jane Tan at:
Email: Jane.tan@Sartorius.com*

Let's spark your next breakthrough!