

Sartorius and *Science* magazine to award prize for regenerative medicine and cell therapy

- Award entails publication in *Science* magazine and prize money of U.S. \$25,000
- Exceptional scientists are invited to apply

Goettingen, Germany, June 27, 2017 – Sartorius, a leading international pharmaceutical and laboratory equipment provider, together with the renowned *Science* magazine will award a Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy. This prestigious annual award is geared towards exceptional scientists who focus on achieving advancements in basic or translational research in these areas. All scientists who have earned a Ph.D. or an M.D. within the past ten years may submit an essay on regenerative medicine, cell therapy, gene therapy or immunotherapy, as well as on materials or tissue engineering.

The winner of the Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy will be awarded U.S. \$25,000 and have his or her essay printed in *Science* magazine. In addition, the winner will receive an attractive Sartorius product package gratis. Moreover, the grand prize essay and those of up to three runners-up will be published in *Science* Online. Runners-up will receive prize money of U.S. \$5,000.

Regenerative medicine and cell therapy address stimulation of the body's repair mechanisms. Furthermore, these therapeutic approaches can lead to technologies allowing tissue and transplantable organs to be grown in the lab. "This area will play a major role in the future. Sartorius' innovative bioanalytical tools help scientists answer fundamental and complex biological questions. We are excited to enable medical progress in these fields, both with our technologies and this award. The prize will be conferred as a token of appreciation to scientists," said Gerry MacKay, member of the Sartorius Group Executive Committee and Executive Vice President of Marketing, Sales and Services for Sartorius' Lab Products & Services Division.

Founded by Thomas Edison, *Science* is published by the American Association for the Advancement of Science (AAAS), the world's largest general scientific society. "*Science* is delighted to join Sartorius in awarding a prize in the exciting area of regenerative medicine and cell therapy. We look forward to promoting basic or translational research in fields ranging from gene therapy to materials engineering, which has the potential to improve human health," said Valda J. Vinson, Deputy Editor Research of *Science*.

Applicants need to submit a 1,000-word essay describing their research and implications for regenerative medicine and cell immunotherapy, as well as supporting documents. The application deadline is October 1, 2017. The awards ceremony will be held in 2018 in Goettingen, Germany.

For more information about the Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy, visit <https://www.passionforscience.com/win-it>

Image files

Logo of the Sartorius & *Science* Prize for Regenerative Medicine & Cell

[Please click here to download](#)

A Profile of Sartorius

The Sartorius Group is a leading international pharmaceutical and laboratory equipment provider with two divisions: Bioprocess Solutions and Lab Products & Services. Bioprocess Solutions with its broad product portfolio focusing on single-use solutions helps customers produce biotech medications and vaccines safely and efficiently. Lab Products & Services, with its premium laboratory instruments, consumables and services, concentrates on serving the needs of laboratories performing research and quality assurance at pharma and biopharma companies and on those of academic research institutes. Founded in 1870, the company earned sales revenue of more than 1.3 billion euros in 2016. Over 6,900 people work at the Group's 50 manufacturing and sales sites, serving customers around the globe.

Contact

Timo Lindemann | Sartorius | Senior Manager Corporate Communications
+49 (0)551.308.4724 | timo.lindemann@sartorius.com
