

Second Sartorius Research Xchange Forum focuses on regenerative medicine and cell therapy

- Event for experts from academia and industry on March 20 and 21, 2018
- Lectures by academia and industry experts
- Guest lecture to be held by Nobel Prize laureate Dr. Stefan W. Hell
- Presentation of "The Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy"

Goettingen, Germany, November 29, 2017 – Promoting interdisciplinary exchange between science and industry on the topics of regenerative medicine and cell therapy: this is the objective Sartorius will be pursuing by hosting its second Research Xchange Forum on March 20 and 21, 2018, at Sartorius College in Goettingen. Experts from the biotechnology sector will report on the new developments in cell therapy. Sartorius has gained guest speakers Dr. Priscilla N. Kelly, editor of the journal *Science*, and Dr. Stefan W. Hell, the winner of the Nobel Prize in Chemistry 2014 and a professor at the Max Planck Institute for Biophysical Chemistry. As part of the Forum, the Sartorius & *Science* Prize will be conferred, which Sartorius together with the scientific journal offers to recognize research contributions made in regenerative medicine and cell therapy. The prize winner and the three finalists will present on their research results.

To attend the Forum, register at the website <https://promotions.sartorius.com/rxf2018>; registrations received up to January 15 will receive an early-bird discount.

Further information:

Research Xchange Forum 2018: <https://promotions.sartorius.com/rxf2018>

The Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy:
<https://www.passionforscience.com/win-it#>

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. More than 6,900 people work at the Group's 50 manufacturing and sales sites, serving customers around the globe.