

Nobel Prize Laureate in Chemistry, Dr. Stefan Hell, Opens the Second Research Xchange Forum at Sartorius

- Experts from science and industry discuss trends and challenges in regenerative medicine and cell therapy
- Microbiologist Kole Roybal awarded the Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy

Göttingen, March 20, 2018 – How to win a Nobel Prize: Dr. Stefan W. Hell gave an inspirational speech on this subject today at the opening of the second Research Xchange Forums hosted by Sartorius. A professor and scientist at the Max Planck Institute for Biophysical Chemistry in Germany, he addressed an audience of more than 140 guests at Sartorius College in Göttingen, describing the path he took that led him to winning a Nobel Prize in 2014 – and what the secrets of such success are. This year's two-day Forum at Sartorius has been focusing on the interdisciplinary exchange between science and industry on the topics of regenerative medicine and cell therapy.

At this expert conference that will continue on Wednesday, a U.S. \$40,000 Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy was awarded for the first time. Microbiologist Kole Roybal of the University of California at San Francisco won first prize for his development of a new class of T-cell immunotherapies that help the immune system to better identify cancer cells. As a result, precision cell therapies can be initiated to combat cancer. The three finalists, Shruti Naik of Rockefeller University of New York, Fotios Sampaziotis of the University of Cambridge, UK, and Will Mclean, formerly at Harvard Medical School, Boston, USA, also presented their research results.

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 1.4 billion euros in 2017. More than 7,500 people work at the Group's approximately 50 manufacturing and sales sites, serving customers around the globe.

Contact

Timo Lindemann | Sartorius Corporate Communications

+49 (0)551.308.4724 | timo.lindemann@sartorius.com