

Third Sartorius Research Xchange Forum 2019 Focuses on CAR-T Cell Therapy

- **Event for experts from academia and industry, including talks, live demo sessions, and panel discussion**
- **Keynote lecture to be held by Bruce Levine, PhD, Barbara and Edward Netter Professor in Cancer Gene Therapy, University of Pennsylvania**
- **Award ceremony of the Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy 2018**

Göttingen, Germany | Bohemia, USA, March 8, 2019 – Promoting interdisciplinary exchange and fostering potential collaborations between academia and industry on highly promising scientific fields relevant to both scientists and businesses: this is the objective of Sartorius in hosting its third Research Xchange Forum (RXF2019) on April 25, 2019, which will spotlight CAR-T cell therapy. The forum is being organized in collaboration with *Science*/AAAS. The American Association for the Advancement of Science (AAAS) is the world's largest general scientific society, and the forum will be held at the AAAS headquarters in Washington D.C. Experts from the academic, clinical and biotech sectors will offer insights on novel developments in CAR-T cell therapies.

Guest speakers will include Prof. Bruce Levine (University of Pennsylvania), Rick Morgan, Sr. VP Immunogenetics (Editas Medicine), and others from institutions such as UCLA, Boston University, MD Anderson Cancer Center, Horizon Discovery, Colorado State University, MaSTherCell and St. Jude Children's Research Hospital.

The Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy will be awarded in an invitation only post-event ceremony. Through sponsoring the award, Sartorius and *Science*/AAAS aim to recognize outstanding research contributions made in the regenerative medicine and cell therapy fields by the next generation of young scientists, the next thought leaders shaping our future as a society. To register free of charge for the RXF2019, please visit our website: <https://www.sartorius.com/us-en/company/conferences/rxf-2019>

A Profile of Sartorius

The Sartorius Group is a leading international partner of biopharmaceutical research and the industry. With innovative laboratory instruments and consumables, the Group's Lab Products & Services Division concentrates on serving the needs of laboratories performing research and quality control at pharma and biopharma companies and those of academic research institutes. The Bioprocess Solutions Division with its broad product portfolio focusing on single-use solutions helps customers to manufacture biotech medications and vaccines safely and efficiently. The Group has been annually growing by double digits on average and has been regularly expanding its portfolio by acquisitions of complementary technologies. In fiscal 2018, the company earned sales revenue of some 1.6 billion euros according to its preliminary figures. Currently, more than 8,100 people work at the Group's approximately 60 manufacturing and sales sites, serving customers around the globe.

About *Science*/AAAS

The American Association for the Advancement of Science (AAAS) is the world's largest general scientific society, and publisher of the journals, *Science*, *Science Translational Medicine*, *Science Signaling*, *Science Advances*, *Science Immunology*, and *Science Robotics*. AAAS was founded in 1848, and includes some 254 affiliated societies and academies of science, serving 10 million individuals. *Science*, founded by Thomas Edison, has the largest paid circulation of any peer-reviewed general science journal in the world, with an estimated total readership of more than 400,000. The non-profit AAAS—www.aaas.org—is open to all and fulfils its mission to "advance science and serve society" through initiatives in science policy, international programs, science education, and more. *Science's* daily online news is always free to the public, as are editorials, any paper with broad public health significance, and all research articles 12 months after publication. *Science* further participates in various efforts to provide free access for scientists in the world's poorest countries.

Contact

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