

## Sartorius enters into a Corporate Partnership Program with the European Molecular Biology Laboratory (EMBL)

- Sartorius and EMBL join forces to offer advanced training and science education
- Sartorius is now a member of the group of corporate EMBL partners aiming to connect industry to academia

Goettingen, February 17, 2017 – Sartorius, a leading international laboratory and pharmaceutical equipment supplier, yesterday announced a partnership with EMBL, the European Molecular Biology Laboratory. EMBL is one of the Europe's leading research institutions with global reach. It includes many research sites, such as Heidelberg and Hamburg in Germany, Hinxton in the UK, Grenoble in France and Monterotondo in Italy.

Sartorius has joined the Corporate Partnership Program, EMBL's platform for fostering advanced training collaboration and activities designed to connect industry to academia. The Corporate Partnership Program includes initiatives such as creating training courses, sponsoring conferences, organizing workshops for professionals and young scientists at the EMBL International Center for Advanced Training, EICAT, which offers numerous courses and conferences each year. In addition, Sartorius will also provide long-term support to EMBL's school education program by funding travel bursaries each year for teachers to attend courses run by EMBL's Learning Laboratory for the Life Sciences, ELLS.

Backing this well-known research institution, Sartorius is offering its expertise to enhance scientific education and practice. "By focusing on professional training at all levels, and ensuring that the next generation of young talent has access to the latest knowledge, we aim to support the scientific community to reach its maximum potential," stated Karen Storm, Vice President of Marketing of the Lab Products & Services Division at Sartorius.

In parallel, EMBL is actively expanding its footprint by strengthening activities beyond the academic sphere into industry. "On behalf of EMBL, we welcome Sartorius to our family of Corporate Partners – working together will enhance our program of internationally renowned events, and we look forward to integrate their technologies and expertise into our offering," added Prof. Matthias Hentze, Director, EMBL.

## **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 according to its preliminary figures. More than 6,900 people work at the Group's approximately 50 manufacturing and sales sites, serving customers around the globe.

---

## **Contact**

Dominic Grone; Senior Manager Corporate Communications  
Sartorius Corporate Administration GmbH, Goettingen, Germany  
Phone: +49.(0)551.308.3324; Fax +49.(0)551.308.3410  
Email: [dominic.grone@sartorius.com](mailto:dominic.grone@sartorius.com); [www.sartorius.com](http://www.sartorius.com)