

Sartorius receives award for outstanding performance in its production chain

Göttingen, November 7, 2018 – Manufacturing operations at Sartorius received a prize for outstanding performance in its value creation chain. At the competition for the Factory of the Year 2018 Award hosted annually by the German production trade journal Produktion and the global management consulting firm A.T. Kearney, the Sartorius manufacturing building for laboratory instruments in Göttingen won the GEO Award – Global Excellence in Operations. This prize will be conferred to Sartorius in March 2019 at the GEO Conference in Ludwigsburg, Germany.

Each of the competing production chains was assessed by performance indicators that were benchmarked, such as customer satisfaction, quality or innovation. These, in turn, involved several assessment criteria, including cost management, resource efficiency and customer focus.

"Our thanks go out to the entire team that faced up to the challenges posed by the competition and to all staff members who have shaped the transformative activities over the past years and helped to implement the changes and improvements. With our drive to provide customers with the best possible service, I see the Global Excellence in Operation Award as confirmation that we are on the right track, and as an incentive to go forward," said Matthias Wessel, Head of Operations of the Sartorius Lab Products & Services Division.

Around 450 employees manufacture laboratory instruments, such as balances, ultrapure water systems and moisture analyzers in the building covering some 25,000 square meters. As high-precision weighing instruments are also calibrated inside, the building is protected from temperature fluctuations and by a vibration-decoupled foundation.

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 2017, Sartorius earned sales revenue of more than 1.4 billion euros. Currently, more than 8,000 people work at the Group's approximately 60 manufacturing and sales sites, serving customers around the globe.

Contact

Timo Lindemann | Corporate Communications

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