

Sartorius Recognized as a Double Award Winner and Promoted to Platinum Supplier Status in Europe by Fisher Scientific

Göttingen, February 2, 2018 – Sartorius, a leading international pharma and laboratory equipment supplier, was a double award winner at the Fisher Scientific European Sales Conference in Malta, picking up the awards for Supplier Excellence and Best Booth. These awards coincided with the formal announcement of the promotion of Sartorius to Platinum supplier status in Europe.

John Collings, European Category Director at Fisher Scientific, announced the change in partnership and explained what it means for the relationship between the two companies. "This change is a significant long-term investment from both companies and will now help us move forward in strategic partnership. This promotion to Platinum will also drive ambitious revenue growth targets, through increased marketing, commercial, supply chain and strategic plans."

The Supplier Excellence Award recognized the excellent cooperation between Fisher Scientific and Sartorius. The year 2017 saw the Sartorius business with Fisher Scientific increase. In addition, the award also recognizes Sartorius for supply chain excellence, product quality and marketing support. The Best Booth award, which is voted for by the Fisher Scientific Commercial team, is awarded to the supplier whose booth best incorporates the four values of Fisher Scientific: involvement, intensity, innovation and integrity.

Daniela Winzker-Demes, Sartorius' Vice President of LPS Sales & Services EMEA, and John Bryson, Director Channel Management Europe, proudly accepted the awards. "The new status is an incentive for us to intensify our cooperation in order to continuously improve our customers' experience," said Winzker-Demes. John Bryson commented: "These awards, and the move to Platinum status, reflect our shared success and close partnership with Fisher Scientific. We look forward to another successful year in 2018 as we work together to drive further growth and deliver solutions to our customers."

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 according to preliminary figures. More than 7,500 people work at the Group's 50 manufacturing and sales sites, serving customers around the globe.

Contact: Timo Lindemann | Sartorius Corporate Communications

Phone: +49(0)551.308.4724 | Email: Timo.Lindemann@Sartorius.com | www.sartorius.com