

## Sartorius employees with up to five decades of seniority celebrate company anniversaries

Göttingen, November 29, 2018 – On Wednesday evening at Sartorius Campus in Göttingen, Sartorius celebrated its 266 employees who each share up to five decades of seniority, representing a grand total of 6,590 years of experience and expertise. During this celebration, CEO Joachim Kreuzburg thanked long-term staff from the German sites in Göttingen, Guxhagen and Laupheim: "Your dedication and knowledge have contributed to making Sartorius so successful today."

Employees who have been with Sartorius for 10, 20, 25, 30, 35 or 40 years, or even longer, were recognized – ranging from a membrane specialist to a cell line expert. Fifty-one people with a seniority of at least 40 years received tribute, among them Dietmar Bode, Guenter Effler, Rudolf Kemmling and Herbert Engelhardt, who each have dedicated 47 years of service to the company, receiving special thanks for their commitment.

Afterwards, long-term staff put their Sartorius knowledge to the test in a company anniversary quiz. Both colleagues and managers alike congratulated them in a video message. Rounding off the evening, the participants exchanged anecdotes about the past and the present while enjoying dinner and live music.

### **A Profile Sartorius 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](mailto:timo.lindemann@sartorius.com) | [www.sartorius.com](http://www.sartorius.com)