

Sartorius wins prize for financial market communication

- **First place in the German TecDax technology index at the Investors' Darling competition**

Goettingen, November 15, 2017 – Sartorius, a leading international pharma and laboratory equipment supplier, was awarded first place for its financial communication in the TecDax technology index at the Investors' Darling capital market competition. In the overall ranking of the 160 largest German joint stock corporations listed, Sartorius achieved seventh place and was the only TecDax corporate group to reach the Top Ten.

"We maintain an ongoing, open dialog with our investors and financial analysts, and our objective is to communicate transparent and understandable information. We are happy about this award as it shows we are moving forward on the right track," said Petra Kirchhoff, Vice President of Corporate Communications and Investor Relations at Sartorius.

The competition held by the German business journal "Manager Magazin" and the Leipzig Graduate School of Management (HHL) has been annually assessing the quality of financial communication of German joint stock corporations since 2014. "Pitted against established DAX30 companies, Sartorius leaves the competition behind regarding the quality of capital market communication. A very impressive achievement," commented Professor Henning Zülch, Chair of Accounting and Auditing at HHL.

The criteria assessed were financial reports, investor relations website and content provided, as well as how the company's reporting is perceived by the financial market. Sartorius moved up from third to first place among the TecDax companies by improving in all three areas compared with last year. Sartorius received the Investors' Darling award on November 14 at Metzler Bank in Frankfurt, Germany.

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. More than 6,900 people work at the Group's 50 manufacturing and sales sites, serving customers around the globe.

Contact: Timo Lindemann, Corporate Communications, Sartorius AG, Goettingen, Germany
Phone: +49.(0)551.308.4724; Email: Timo.Lindemann@Sartorius.com, www.sartorius.com