

Boston Consulting Group again ranks Sartorius shares as among the best in the world

- Second place among German stocks
- Average total shareholder return (TSR) of 36.6% over the past five years

Göttingen, July 13, 2018 – Shares of the Sartorius Group, a leading international partner of the biopharmaceutical industry and the research sector, again won a top slot in the global Value Creators rankings of the Boston Consulting Group (BCG). In this annual analysis, an eight-member team of experts determines which of the more than 2,400 shares worldwide deliver the best TSR. After Sartorius had won the ranking spot a year ago as the German share with the highest return, the company was pleased to be ranked second this year. Sartorius also performed well in the rankings by industry: Only four German stocks number among the best of their sector and Sartorius was one of these, ranking ninth in medical technology.

The key metric for determining rankings is how much value a share generates by price increases and dividends. In assessing top value creators, BCG analysts tracked TSR over a five-year period from 2013 through 2017, in which the Sartorius share delivered a total return averaging 36.6% per year for its shareholders.

A Profile

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

Contact

Petra Kirchhoff | Corporate Communications
+49 (0)551.308.1686 | petra.kirchhoff@sartorius.com | www.sartorius.com