

Göttingen, March 17, 2020

## Sartorius postpones Annual Shareholders' Meeting

Sartorius will postpone its Annual General Shareholders' Meeting formerly scheduled for March 26, 2020, to a later date due to the coronavirus pandemic and the associated restrictive rules imposed on meetings. This was resolved today by the Executive Board of the company.

As a result of the cancellation of the Annual General Shareholders' Meeting on March 26, 2020, the resolution on the appropriation of the retained profit for 2019 and the payment of dividends, among other things, will also be deferred.

The Annual General Shareholders' Meeting is to be scheduled for a new date within the period of the first eight months of the current fiscal year as provided for under the German Stock Corporation Law. Sartorius will inform its shareholders and the public in a timely manner about its further planning.

### A profile of Sartorius

The Sartorius Group is a leading international partner of life science research and the biopharmaceutical 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 2019, the company earned sales revenue of some 1.83 billion euros. At the end of 2019, more than 9,000 people were employed at the Group's approximately 60 manufacturing and sales sites, serving customers around the globe.

### Contact

Petra Kirchhoff  
Head of Corporate Communications and Investor Relations  
+49 (0)551.308.1686  
[petra.kirchhoff@sartorius.com](mailto:petra.kirchhoff@sartorius.com)  
sartorius.com

Follow Sartorius on Twitter [@Sartorius\\_Group](https://twitter.com/Sartorius_Group) and on LinkedIn.