

Göttingen, December 17, 2020

## Sartorius refinances syndicated loan and increases financing flexibility

- Credit volume now amounts to 600 million euros
- International bank syndicate led by Commerzbank, BNP Paribas, and UniCredit

The life science group Sartorius refinanced its syndicated loan, increasing it to 600 million euros. The company signed a loan agreement with an international bank syndicate led by BNP Paribas, Commerzbank, and UniCredit. Sartorius has now adjusted its credit facility in place so far to accommodate its strong growth. As a result, the company has raised its credit line that will run until the end of 2021 by 200 million euros and extended the members of the bank syndicate by J.P. Morgan and UniCredit.

“The new loan agreement ensures our financing flexibility following the company’s three acquisitions in 2020, providing leeway for further strategic measures. We took out this loan under highly attractive terms in a market environment of tightening financial conditions. This reflects the confidence of banks in the financial strength and positive prospects of Sartorius,” said Rainer Lehmann, member of the Executive Board and Chief Financial Officer of Sartorius. The credit line is for a term of three years and can be extended twice for one year at a time.

### 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 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

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)

Follow Sartorius on [Twitter](#) @Sartorius\_Group and on [LinkedIn](#).