

Göttingen, September 9, 2020

## Sartorius successfully places note loan with a volume of 750 million euros

- Transaction oversubscribed more than three times
- Volume increased from 300 million euros to 750 million euros

Sartorius, a leading international partner of life science research and the biopharmaceutical industry, successfully placed a note loan with a volume of 750 million euros. The note loan was issued with maturity terms of 3, 5, 7 and 10 years and was significantly oversubscribed. “We see such great interest as clear proof of our attractiveness for investors,” said Rainer Lehmann, Chief Financial Officer of Sartorius.

In view of high demand, the targeted minimum volume of 300 million euros was significantly exceeded. Most of the funds will be used for early repayment of the bridge loan that Sartorius had taken out at the end of April in connection with the acquisition of selected life science businesses from Danaher.

This press release contains forward-looking statements about the future development of the Sartorius Group. Forward-looking statements are subject to known and unknown risks, uncertainties and other factors that could cause actual results to differ materially from those expressed or implied by such statements. Sartorius assumes no liability for updating such statements in light of new information or future events. This is a translation of the original German-language press release. Sartorius shall not assume any liability for the correctness of this translation. The original German press release is the legally binding version.

### 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

Timo Lindemann  
Corporate Communications  
+49 (0)551.308.4724  
[timo.lindemann@sartorius.com](mailto:timo.lindemann@sartorius.com)  
sartorius.com

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