

Göttingen, December 7, 2020

Sartorius pays employees a coronavirus bonus

- Thanks and appreciation for commitment and flexibility of employees during the coronavirus pandemic as well as recognition of additional burdens entailed
- Bonus payments to total around ten million euros

The life science company Sartorius will pay a total of around ten million euros in coronavirus bonuses to its employees in appreciation of their extraordinary commitment and special achievements.

“The year 2020 was highly challenging for all of us, both privately and in our jobs. Particularly our employees in production, logistics, labs and other areas who did not have any option of working from home faced especially difficult conditions. As Sartorius with its products is directly involved in the development and manufacture of coronavirus vaccines and Covid-19 therapeutics, it was crucial to maintain our delivery capability. We were able to master this challenge successfully thanks to the great commitment and flexibility of all our employees. For this reason, we have decided to pay out a total of around ten million euros as one-time special bonuses in recognition of what has been achieved,” said Sartorius Executive Board Chairman and CEO Joachim Kreuzburg.

The special bonuses approved by the Executive Board apply to the more than 10,000 employees of Sartorius worldwide, in addition to the company’s temporary agency workers, and will be paid out in December 2020.

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

Petra Kirchhoff

Head of Corporate Communications

+49 (0)551.308.1686

petra.kirchhoff@sartorius.com

www.sartorius.com

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