

Göttingen, February 16, 2024

## Sartorius releases Annual Report for 2023

The life science group Sartorius has released its Annual Report for 2023 today. The document is available under the following link: <https://ir-reports.sartorius.com/en/ag/fy-2023>.

### Financial calendar

March 28, 2024	Annual General Meeting
April 18, 2024	Publication of first-quarter figures (January to March 2024)
July 19, 2024	Publication of first-half figures (January to June 2024)
October 17, 2024	Publication of nine-month figures (January to September 2024)

### A profile of Sartorius

Sartorius is a leading international partner of life sciences research and biopharmaceutical manufacturing. With innovative laboratory instruments and consumables, the Group's Lab Products & Services division focuses on laboratories performing research and quality control at pharmaceutical and biopharmaceutical companies as well as academic research institutes. The Bioprocess Solutions division, with its broad product portfolio focusing on single-use solutions, helps customers manufacture biotech medications, vaccines, and cell and gene therapies safely, rapidly, and economically. The company, based in Göttingen, Germany, has a strong global reach with around 60 production and sales sites worldwide. Sartorius regularly expands its portfolio through the acquisition of complementary technologies. In 2023, the company generated sales revenue of around 3.4 billion euros. Currently, around 14,600 employees are working for customers around the globe.

### Contact

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

Follow Sartorius on [LinkedIn](#).