

Göttingen, February 17, 2023

Sartorius releases Annual Report for 2022

The life science group Sartorius has released its Annual Report for 2022 today. The document is available under the following link: <https://ir-reports.sartorius.com/en/ag/fy-2022>.

Financial calendar

| | |
|------------------|---|
| March 29, 2023 | Annual General Meeting |
| April 20, 2023 | Publication of first-quarter figures (January to March 2023) |
| July 21, 2023 | Publication of first-half figures (January to June 2023) |
| October 19, 2023 | Publication of nine-month figures (January to September 2023) |

A profile of Sartorius

The Sartorius Group is a leading international partner of life sciences 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 pharmaceutical and biopharmaceutical companies and those of academic research institutes. The Bioprocess Solutions division, with its broad product portfolio focusing on single-use solutions, helps customers manufacture biotech medications and vaccines safely and efficiently. The company is growing at double-digit rates on average per year and regularly expands its portfolio through the acquisition of complementary technologies. In fiscal 2022, the company generated sales revenues of around 4.2 billion euros. At the end of 2022, around 16,000 employees were working for customers around the globe at the group's 60 or so production and sales sites.

Contact

Philipp Grontzki
Head of External Communications
+49 (0)551.308.5581
philipp.grontzki@sartorius.com

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