

Göttingen, Germany | February 6, 2026

## The Supervisory Board of Sartorius AG proposes a dividend of 0.74 euros per preference share

At today's meeting, the Supervisory Board of Sartorius AG approved the Executive Board's recommendation to submit a proposal to the Annual General Meeting on March 26, 2026, to pay a dividend of 0.74 euros per preference share and 0.73 euros per ordinary share for fiscal 2025, the same amount as in 2024. The total distributed profit would accordingly be 50.8 million euros. The dividend will be paid on March 31, 2026.

This media 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.

### Financial calendar

February 16, 2026	Publication of the 2025 Annual Report
March 26, 2026	Annual General Meeting
April 23, 2026	Publication of the first quarter results for January to March 2026
July 23, 2026	Publication of half-year results for January to June 2026
October 22, 2026	Publication of nine-month results for January to September 2026

### A profile of Sartorius

Sartorius is a leading international partner to the biopharmaceutical research and manufacturing industries. The Lab Products & Services Division focuses on innovative laboratory instruments and consumables for research and quality assurance laboratories in pharmaceutical and biopharmaceutical companies as well as academic research institutions. The Bioprocess Solutions Division supports customers with a broad product portfolio focused on single-use solutions for the safer, faster, and more sustainable production of biotech drugs, vaccines, and cell and gene therapies. With around 60 production and sales locations worldwide, the Göttingen-based company has a strong global presence. Sartorius regularly supplements its portfolio with acquisitions of complementary technologies. In 2025, according to preliminary figures, the company generated sales revenue of around 3.5 billion euros. More than 14,000 employees serve customers around the globe.

Visit our [newsroom](#) and follow us on [LinkedIn](#).

### Contact

Verena Sattel  
External Communications  
+49 551 308 9261  
[verena.sattel@sartorius.com](mailto:verena.sattel@sartorius.com)