

Göttingen, Germany | April 24, 2026

Sartorius AG Supervisory Board extends appointment of board member Florian Funck

The Supervisory Board of Sartorius AG resolved to extend the appointment of Dr. Florian Funck as Chief Financial Officer (CFO) and member of the Executive Board ahead of schedule to March 31, 2032.

Funck (55) studied business administration at the University of Münster, where he also earned his doctorate. He subsequently spent more than two decades with the Haniel Group, most recently serving as CFO from 2011 to 2023. Funck was appointed to the Executive Board of Sartorius in April 2024; in this role, he is responsible for Finance & Controlling, IT & Process, Corporate Compliance, Internal Audit and Corporate Sourcing. In addition, Funck is a member of the Supervisory Board of Vonovia SE and an honorary professor at the University of Münster.

The Executive Board of Sartorius AG also comprises Chairman and Chief Executive Officer Dr. Michael Grosse, who oversees Strategy, Human Resources, Corporate Research, Legal Affairs, and Communications & IR, as well as the two Heads of Division Dr. René Fáber (Bioprocess Solutions) and Dr. Alexandra Gatzemeyer (Lab Products & Services).

Further Information

[CV of Florian Funck](#)

Financial calendar

July 23, 2026	Publication of half-year results for January to June 2026
October 22, 2026	Publication of nine-month results 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, the company generated sales revenue of around 3.5 billion euros. More than 14,000 employees serve customers around the globe.

Visit our [newsroom](#) or follow Sartorius on [LinkedIn](#).

Contact

Leona Malorny

Head of External Communications

+49 551 308 4067

leona.malorny@sartorius.com