

## Curriculum Vitae | Dr. Michael Grosse

### Chairman of the Executive Board of Sartorius AG

Michael Grosse holds a doctorate in mechanical engineering and has held various management and board positions in the packaging industry for the pharmaceutical and food sector over the past 20 years.

From 2020 to 2023, he was Chairman of the Management Board and Chief Executive Officer of Syntegon Technology GmbH, an international market leader in process and packaging solutions based in Germany, which was spun off from the Bosch Group in 2019. Prior to that, Grosse worked for the Swiss company Tetra Pak for many years, where he became a member of the Executive Board in 2006, and where he was responsible for the expansion of the global service business and held various management positions in product development. Before joining Tetra Pak, he worked in the automotive industry at BMW and Ford.

#### Personal Information

Year of birth: 1967

Nationality: German

End of current ordering period: June 30, 2028

#### Professional Profile

Since July 2025	Chief Executive Officer at Sartorius AG <sup>1</sup> , Germany
2024–2025	Independent Senior Advisor, Germany
2020–2023	Chief Executive Officer at Syntegon Technology GmbH, Germany
2003–2019	Several positions at Tetra Pak International S.A.
	2018–2019   Executive Vice President Services, Switzerland
	2011–2018   Executive Vice President Development & Service Operations, Switzerland
	2006–2011   Executive Vice President Development & Engineering, Sweden
	2003–2006   Vice President Development, Italy
1999–2003	Several positions at Ford Motor Company <sup>1</sup> , Germany
	2001–2003   Chief Carline Engineer Large & Luxury Cars
	1999–2001   Product & Business Manager Ford Focus
1993–1999	Several positions at BMW AG <sup>1</sup> , Germany
	1997–1999   Manager Production Technologies
	1995–1997   Manager of Corporate Product Strategy

1991–1993      Managing Director at BSB GmbH, Germany

### **Educational Background**

1991–1994      Dr. Ing./PhD, Mechanical Engineering, Technical University Braunschweig, Germany

1992–1994      Business Administration, University of Regensburg, Germany

1987–1991      Dipl. Ing./ Master, Mechanical Engineering, Technical University of Munich, Germany

### **Other current directorships and positions**

- Member of the Board of Advisors at Mustad Hoofcare Group, Switzerland

<sup>1</sup> Public listed