

Curriculum Vitae | Dr. Michael Grosse

Chairman of the Board of Directors of Sartorius Stedim Biotech S.A.

Year of birth: 1967

Nationality: German

First appointment to the Board of Directors: July 2025

End of term: 2027 Shareholders' Meeting

Professional Profile

Since July 2025	Chief Executive Officer at Sartorius AG ¹ , 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 ¹ , Germany 2001-2003 Chief Carline Engineer Large & Luxury Cars 1999-2001 Product & Business Manager Ford Focus
1993-1999	Several positions at BMW AG ¹ , 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 within the Group

- Chairman of the Executive Board (Vorstandsvorsitzender) of Sartorius AG¹, Germany

Other current directorships and positions outside the Group

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

¹ Public listed