

Curriculum Vitae | Dr. Joachim Kreuzburg

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

Year of Birth: 1965 Nationality: German

First appointment to the Board of Directors: June 2007

End of term: 2025 Shareholders' Meeting

Professional Profile

June 29, 2007- March 26, 2023	CEO of Sartorius Stedim Biotech S.A. ¹ , Aubagne, France
Since Nov. 11, 2005	CEO and Executive Board Chairman of Sartorius AG ¹ , Germany
May 1, 2003- Nov. 10, 2005	Spokesman (Sprecher) of the Executive Board of Sartorius AG ¹ , Germany
Since Nov. 11, 2002	Member of the Executive Board of Sartorius AG ¹ , Germany
Since May 1, 1999	Sartorius AG ¹ , Germany. Most recent position before promotion to the Executive Board: Vice President, Finance and Investor Relations
1995-1999	University of Hannover, Germany - Research associate at the Faculty of Economics and Management
1992-1995	Institute for Solar Energy Research - Research associate in Hameln, Germany

Educational Background

- Diplom-Maschinenbau-Ingenieur, (university degree in mechanical engineering), University of Hanover, Germany
- Dr. rer. pol., Wirtschaftswissenschaften (doctorate in economics), University of Hanover, Germany

Other current directorships and positions within the Group

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

¹ Public listed

As of February 2024 1|2

- Managing Director of Sartorius Lab Holding GmbH, Germany
- Managing Director of SI Weende-Verwaltungs-GmbH, Germany
- Managing Director of SI Grone 1-Verwaltungs-GmbH, Germany
- Managing Director of SIV Grone 2 GmbH, Germany
- Managing Director of Sartorius Ventures GmbH, Germany
- Chairman of the Supervisory Board of Sartorius Corporate Administration GmbH, Germany
- Chairman of the Advisory Board of LabTwin GmbH, Germany
- Chairman of the Board of Directors of Sartorius North America Inc., United States

Other current directorships and positions outside the Group

- Member of the Supervisory Board (Aufsichtsrat) of Carl Zeiss AG¹, Germany
- Member of the Administrative Board (Verwaltungsrat) of Ottobock Management SE, Germany

As of February 2024 2 | 2