

Curriculum Vitae | Dr. Joachim Kreuzburg

Chairman of the Executive Board of Sartorius AG

Joachim Kreuzburg, born in 1965, has been at the helm of the Executive Board of Sartorius AG since 2003. On the Executive Board, he is responsible for Strategy, Human Resources, Corporate Research, and Legal and Compliance, as well as Communications & Investor Relations. Since 2007, he has also been Chairman of the Board of Directors of the publicly listed subgroup Sartorius Stedim Biotech S.A.

After earning a university degree in mechanical engineering in Hannover, Joachim Kreuzburg first worked at the Lower Saxony Institute for Solar Energy Research in Hameln and at the Faculty of Economics at the University of Hannover, where he received his doctorate in 1999.

In his subsequent career at Sartorius, he held various leadership positions in Finance and Controlling before being appointed to the Executive Board of Sartorius AG in 2002.

In addition to his activities at Sartorius, Joachim Kreuzburg is, inter alia, a member of the Supervisory Board of Carl Zeiss AG, Non-Executive Director of the Administrative Board of Ottobock Management SE, Member of the Senate of acatech – National Academy of Science and Engineering, Member of the Executive Board of Stifterverband für die Deutsche Wissenschaft, and Member of the Foundation Council of the University of Göttingen.

Personal Information

Year of birth: 1965

Nationality: German

End of current ordering period: November 10, 2025

Professional Profile

June 29, 2007-
March 26, 2023

CEO of Sartorius Stedim Biotech S.A.¹, France

Since Nov. 11,
2005

CEO and Executive Board Chairman of Sartorius AG¹, Germany

May 1, 2003-
Nov. 10, 2005

Spokesman (Sprecher) of the Executive Board of Sartorius AG¹, Germany

Since Nov. 11,
2002

Member of the Executive Board of Sartorius AG¹, Germany

Since May 1,
1999

Sartorius AG¹, Göttingen, Germany. Most recent position before promotion to the Executive Board: Vice President, Finance, and Investor Relations

1995-1999 University of Hanover, 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, Dr. rer. pol. (University degree in mechanical engineering, doctorate in economics), University of Hanover, Germany

Other current directorships and positions within the Group

- Chairman of the Board of Directors of Sartorius Stedim Biotech S.A.¹, France
- 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
- 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

Other current directorships and positions outside the Group

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

¹ Public listed