

## Chairman of the Executive Board | CEO

**Dr. Joachim Kreuzburg**

### **Career**

Joachim Kreuzburg, born in 1965, has been at the helm of the Executive Board of Sartorius AG since May 2003. He is responsible for Strategy, Human Resources and Corporate Research, as well as Legal, Compliance and Communications. Since 2007, Mr. Kreuzburg has also been Chairman of the Board of Directors and CEO of Sartorius Stedim Biotech S. A.

After earning a university degree in mechanical engineering in Hanover, Germany, Mr. Kreuzburg worked at the Lower Saxony Institute for Solar Energy Research in Hamlin, Germany, and at the Faculty of Economics at the University of Hanover, where he received his doctorate in 1999. In the same year, Mr. Kreuzburg began his career at Sartorius. Initially, he held various positions in Controlling, and then became Commercial Managing Director of the subsidiary GWT based in Hamburg, Germany. Afterwards, he was in charge of Financing and Investor Relations before being appointed to the Executive Board of Sartorius AG in November 2002.

### **Other Activities**

In addition to his work for Sartorius, Mr. Kreuzburg is also a member of the Supervisory Board of Carl Zeiss AG, a non-executive director on the Management Board of Ottobock and a member of the Economic Advisory Board of Norddeutsche Landesbank.