

## Chairman of the Executive Board | CEO

**Dr. Joachim Kreuzburg**

### **Career**

Joachim Kreuzburg, born in 1965, has been Chairman of the Executive Board of Sartorius AG since 2005 and is responsible for Strategy, Human Resources and Corporate Research, as well as Legal, Compliance and Communications. Since 2007, he 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, he 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, he began his career at Sartorius. Initially, Kreuzburg held various positions in Controlling, then became of 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, which he has been heading since May 2003.

### **Other Activities**

In addition to his commitment to Sartorius AG, Kreuzburg is, inter alia, a member of the Supervisory Board of Carl Zeiss AG, Vice Chairman of the Supervisory Board of Ottobock SE & Co. KGaA and a member of the Economic Advisory Board of the Nord LB bank.