

Curriculum Vitae | Beatrix Henseler

Member of the Supervisory Board of Sartorius AG

Year of birth: 1965

Nationality: German

Place of residence: Mettmann, Germany

Member since: September 2025

Professional Profile

Since 2019	Sartorius Group, Göttingen, Germany – Head of Human Resources
2015-2019	CLIP TRAINING, Mettmann, Germany – Managing Director and Founder
2013-2019	Consultant and Interim Manager, Mettmann, Germany
2010-2013	TK Maxx GmbH & Co KG, Düsseldorf, Germany – Vice President Human Resources
2002-2010	Douglas Holding AG, Hagen, Germany
	2007-2010: Director Human Resources & Organization
	2004-2007: Director Human Resources
	2002-2004: Head of Human Resources
2000-2002	ESCADA AG, Munich, Germany – International HR Director
1997-2000	Philip Morris GmbH, Munich, Germany – EVP Policies & Systems, Head of Central HR
1991-1997	Mars GmbH, Viersen, Germany – Management Positions in Sales and HR
1990-1991	H. Bahlsen Keksfabrik KG, Hannover, Germany – Marketing Trainee
1989-1990	FAA company for vocational training, Gelsenkirchen, Germany – Lecturer

Educational Background

- Postgraduate Studies in Organizational Psychology, Universität Dortmund, Germany
- Economics / Business Administration („Diplom-Kauffrau“), Universität Essen, Germany
- Business Administration, Universität Bielefeld, Germany

Membership in Other Supervisory Boards Mandated by German Law

- Sartorius Stedim Biotech GmbH, Göttingen, Germany – Member of the Supervisory Board

Membership in Comparable Supervisory Committees of Business Enterprises Within Germany & Abroad

- None

Other Major Activities Beyond Supervisory Board Position

- None