

# Curriculum Vitae | Prof. Dr. Klaus Trützschler

## Member of the Supervisory Board of Sartorius AG

Year of birth: 1948
Nationality: German

Place of residence: Essen, Germany Member since: November 2011

#### **Professional Profile**

Since 2012	Freelance business consultant
2008-2013	Westphalian Wilhelms University of Münster, Germany - Honorary Professor of the Chair for Business Administration, especially for Organization, Personnel and Innovation
2000-2012	Franz Haniel & Cie. GmbH, Duisburg, Germany – member of the Executive Board   CFO
1977-2000	Ruhrkohle AG   RAG Aktiengesellschaft, Essen, Germany
	1997–2000: Member of the Executive Board   CFO
	1997–2000: Member of the Executive Board   CFO 1995–1996: Chief Representative & Head of Company Development, RAG
	·

#### Educational Background

- Doctorate in Economics ("Dr. rer. pol.") Munich University of Applied Sciences (TU München),
   Germany
- Graduate Mathematician ("Dipl.-Mathematiker") Universities of Würzburg / Freiburg / Bonn,
   Germany
- Graduate Mathematician ("Dipl.-Wirtschaftsmathematiker") RWTH Aachen University / TU München

#### Membership in Other Supervisory Boards Mandated by German Law

None

As of January 2024 1 | 2

# Membership in Comparable Supervisory Committees of Business Enterprises Within Germany and Abroad

None

### Other Major Activities Beyond Supervisory Board Position

• Chairman of the Board of Trustees of Stiftung Universitätsmedizin Essen

As of January 2024 2 | 2