

## Member of the Supervisory Board of Sartorius AG

Dr. Klaus Trützscher, Professor; resident of Essen, Germany  
Born 1948

### Professional Profile:

- 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  
1995–1996: Chief Representative & Head of Company Development (ZB Unternehmensentwicklung), RAG  
1992–1995: Rütgerswerke and double positions as CFO of Bakelite AG / CFO of Rütgerswerke AG  
1977–1992: Various positions (focal areas: corporate accounting, general accounting)

### 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

**Positions Held Within the Meaning of Sec. 125, Subsec. 1, Sentence 5, of the German Stock Corporation Law (AktG):**

- Wuppermann AG, Leverkusen, Germany – Supervisory Board (Chairman)
- Zwiesel Kristallglas AG, Zwiesel, Germany – Supervisory Board (Chairman)
- Odenwald Faserplattenwerk GmbH, Amorbach, Germany - Advisory Board

**Major Activities Beyond Supervisory Board Position:**

- Chairman of the University Council of the University of Duisburg-Essen
- Positions Held Within the Meaning of Sec. 125, Subsec. 1, Sentence 5, of the German Stock Corporation Law (AktG)