

## 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  
1995–1996: Chief Representative & Head of Company Development, 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

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

- Lennertz & Co. Capital GmbH, Hamburg – Chairman of the Supervisory Board

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

- None

## Other Major Activities Beyond Supervisory Board Position

- Chairman of the Board of Trustees of Stiftung Universitätsmedizin Essen