

Curriculum Vitae | Frank Riemensperger

Member of the Supervisory Board of Sartorius AG

Year of birth: 1962

Nationality: German

Place of residence: Dietzenbach, Germany

Member since: March 2022

Professional Profile

- | | |
|------------|---|
| Since 2022 | Founder and Managing Director, 440.digital GmbH, consulting and investments in digital start-ups |
| 1989–2022 | Accenture

2009–2021: CEO of Accenture Germany. Member of the Global Leadership Council, then of the Global Management Committee of the Accenture Group.

2005–2009: Head of the Technology Business Unit for Accenture’s global sector group Manufacturing Industry (Products)

2002–2007: Deputy Spokesman for management in Germany

1998: Appointed as a partner to the company

Project head for complex, digital transformations in large corporations. |
| 1987–1988 | Integrata Unternehmensberatung AG, Tübingen – Executive Assistant |

Educational Background

- Graduate Degree in Computer Science (“Diplom-Informatiker”) in 1987 from the University of Furtwangen (HFU), University of Applied Sciences

Membership in Other Supervisory Boards Mandated by German Law

- Drägerwerk Verwaltungs AG, Lübeck – member of the Supervisory Board
- Dräger Safety Verwaltungs AG, Lübeck – member of the Supervisory Board

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

- DRM Datenraum Mobilität GmbH, München – member of the Supervisory Board

Other Major Activities Beyond Supervisory Board Position

- Member of the Executive Committee of Acatech (German Academy of Technical Sciences)
- Member of steering committee of the platform Learning Systems (AI), "Plattform Lernende Systeme (KI)"
- Member of the Board of Trustees of Senckenberg Gesellschaft
- Member of the Feldafing district group (discussion group focusing on CTO and digitalization)
- Member of the Baden-Baden Entrepreneur Talks (BBUG)