

## Curriculum Vitae | Ilke Panzer

### Member of the Supervisory Board of Sartorius AG

Year of birth: 1966

Nationality: German

Place of residence: Fredonia, Wisconsin, USA

Member since: April 2017

#### Professional Profile

- |            |   |
|------------|---|
| Since 2022 | Freelance consultant in the field of healthcare innovation  |
| 2018–2022  | Assurance Laboratories LLC, USA – CEO   |
| 2016–2018  | Versiti Inc., USA – Executive Vice President   Chief Innovation Officer  <br>General Manager, Diagnostic Laboratories   |
| 2006–2016  | BloodCenter of Wisconsin, USA<br>2009–2016: Senior Vice President, Diagnostic Laboratories<br>2006–2009: Vice President, Diagnostic Laboratories  |
| 2004–2006  | Johnson & Johnson, DePuy Orthopaedics, Inc., USA – Vice President, Worldwide<br>Research & Development  |
| 1992–2004  | General Electric Corporation, USA<br>2000–2004: General Manager, Global Ultrasound<br>1999–2000: Global Supply Chain Manager, Ultrasound<br>1998–1999: Quality Manager, Six Sigma Champion, Ultrasound<br>1996–1998: Engineering Manager, Computer Tomography<br>1992–1996: Edison Engineer, Electrical Distribution and Controls |

#### Educational Background

- Master of Science in Engineering, Computer and Systems Engineering National Technological University, Fort Collins, Colorado, USA
- Bachelor of Science in Engineering, Computer and Systems Engineering University of Connecticut

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

- None

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

- None

#### **Other Major Activities Beyond Supervisory Board Position**

- Vice Chair on the Board of the Wisconsin Technology Council
- Member of the Investment Advisory Committee of a healthcare venture fund