

Member of the Supervisory Board of Sartorius AG

Ilke Panzer, resident of Fredonia, Wisconsin, USA
Born 1966

Professional Profile:

- Since 2018 Assurance Laboratories LLC, USA – CEO
- 2016–2018 Versiti Inc., USA – Executive Vice President / Chief Innovation Officer /
General Manager, Diagnostic Laboratories
- 2006–2016 BloodCenter of Wisconsin, USA
2009–2016: Senior Vice President, Diagnostic Laboratories

2006–2009: Vice President, Diagnostic Laboratories
- 2004–2006 Johnson & Johnson, DePuy Orthopaedics, Inc., USA – Vice President, Worldwide
Research & Development
- 1992–2004 General Electric Corporation, USA
2000–2004: General Manager, Global Ultrasound

1999–2000: Global Supply Chain Manager, Ultrasound

1998–1999: Quality Manager, Six Sigma Champion, Ultrasound

1996–1998: Engineering Manager, Computer Tomography

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

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

- None

Major Activities Beyond Supervisory Board Position:

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