

Mitglied des Aufsichtsrats der Sartorius AG

Ilke Panzer, Fredonia, WI (USA)
geboren 1966

Beruflicher Werdegang:

- seit 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

Ausbildung:

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

Mandate i.S.d. § 125 Abs. 1 Satz 5 AktG:

- keine

Wesentliche Tätigkeiten neben dem Aufsichtsratsmandat:

- CEO von Assurance Laboratories LLC
- Vice Chair im Board des Wisconsin Technology Council
- Mitglied im Strategic Planning Committee of the National Blood Foundation
- Mitglied im Investment Advisory Committee eines Healthcare Venture Funds