

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 Investment Advisory Committee eines Healthcare Venture Funds