

Member of the Supervisory Board of Sartorius AG

Dr. Thomas Scheper, Professor; resident of Hanover, Germany
Born 1956

Professional Profile:

- | | |
|------------|---|
| Since 1995 | Institute of Technical Chemistry at the Gottfried Wilhelm Leibniz University in Hanover, Germany - University Professor (currently head of the institute) |
| 1992-1995 | Westphalian Wilhelms University of Münster, Germany – Professor of Biotechnology |
| 1985-1992 | Gottfried Wilhelm Leibniz University in Hanover, Germany, and at Caltech, Pasadena, California, USA – Postdoctoral Position |

Educational Background:

- Doctorate in Natural Sciences (“Dr. rer. nat.”) – University of Hannover, Germany
- Graduate Chemical Engineer (“Dipl.-Chemiker”) – University of Hannover, Germany

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:

- University professor and head of the Institute of Technical Chemistry at the Gottfried Wilhelm Leibniz University in Hanover, Germany (research focus: bioprocess engineering, purification, bioanalytics, regenerative medicine / tissue engineering, enzyme technology)