

## Member of the Supervisory Board of Sartorius AG

Dr. Lothar Kappich, resident of Hamburg, Germany  
Born 1957

### Professional Profile:

Since 2017	Independent Business Consultant
1998-2017	ECE Projektmanagement GmbH & Co. KG, Hamburg, Germany – Managing Director
1993-1998	ECE Consulting GmbH, Hamburg, Germany – Managing Director
1990-1992	ECE Projektmanagement GmbH & Co. KG, Hamburg, Germany – Assistant CEO
1988-1990	Schering AG, Berlin, Germany – Controller
1984-1988	Institute of Quantitative Economic Research “Institut für Quantitative Wirtschaftsforschung” (Department of International Business Relations “Internationale Wirtschaftsbeziehungen”), University of Hannover, Germany – Scientific Associate

### Educational Background:

- Doctorate in Economics (“Dr. rer. pol.”) - University of Hannover, Germany
- Graduate Economist (“Dipl.-Ökonom”) - University of Hannover, Germany
- Industrial Business Manager (“Industriekaufmann”) - XOX Nabisco GmbH, Celle, Germany

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

- Sartorius Stedim Biotech S.A., France (stock exchange listed / a subgroup of Sartorius AG) – Board of Directors

### Major Activities Beyond Supervisory Board Position:

- Independent Business Consultant
- Executor of Horst Sartorius’ will
- Position held within the meaning of Sec. 125, Subsec. 1, Sentence 5, of the German Stock Corporation Law (AktG)