

## Curriculum Vitae | Dr. Lothar Kappich

### Member of the Supervisory Board of Sartorius AG

Year of birth: 1957

Nationality: German

Place of residence: Hamburg, Germany

Member since: April 2007

#### Professional Profile

- Since 2017 Independent Business Consultant
- 1998-2017 ECE Projektmanagement GmbH & Co. KG, Hamburg, Germany – Managing Director
- 1995-2017 Managing Director at diverse subsidiaries of ECE Group
- 1993-1998 ECE Consulting GmbH, Hamburg, Germany – Managing Director
- 1990-1992 ECE Projektmanagement GmbH & Co. KG, Hamburg, Germany – Assistant to the 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

#### Membership in Other Supervisory Boards Mandated by German Law

- None

#### Membership in Comparable Supervisory Committees of Business Enterprises Within Germany and Abroad:

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

#### Other Major Activities Beyond Supervisory Board Position

- Executor of Horst Sartorius’ will