

## Curriculum Vitae | Lothar Kappich

Member of the Board of Directors of Sartorius Stedim Biotech S.A.

Year of Birth: 1957

Nationality: German

First appointment to the Board of Directors: September 2017

End of term: 2029 Shareholders' Meeting

### Professional Profile

Since 2017	Chairman of the Supervisory Board of Sartorius AG <sup>1</sup> , Germany
2007-2017	Member of the Supervisory Board of Sartorius AG <sup>1</sup> , Germany
1990-2017	ECE Projektmanagement G.m.b.H. & Co. KG, Germany, latest position Managing Director of ECE's HR & Corporate Services as well as Managing Director of numerous subsidiaries at the ECE group
1988-1990	Schering AG, Germany, Controller in the Central Controlling Department

### Educational Background

- Doctorate (Dr. rer. pol.) in economics (subject of the doctoral dissertation: Theory of International Business Activity), University of Hannover, Germany

### Other current directorships and positions within the Group

- Chairman of the Supervisory Board of Sartorius AG<sup>1</sup>, Germany

### Other current directorships and positions outside the Group

- None

<sup>1</sup> Public listed