

## 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: 2025 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