

## Member of the Supervisory Board of Sartorius AG

Dietmar Müller, Resident of Gleichen, Germany  
Born 1976

### Professional Profile

Since 1997 Sartorius Group, Göttingen, Germany  
Since 2018: Member and Chairman of the Employees' Council of Sartorius Stedim Biotech GmbH; exempted from regular employment duties  
2016–2018: In-house Sales Representative, BPS Central Europe  
2012–2016: Team Lead, Commercial Operations Central Europe  
2004–2012: Customer Support for Austria  
1999–2003: Junior Customer Service for Germany

### Educational Background

- Business Economist ("Betriebswirt") – degree earned at the VWA vocational academy in Göttingen, Germany
- Industrial Business Manager ("Industriekaufmann") – vocational degree completed at Sartorius AG, Göttingen, Germany

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

- Sartorius Stedim Biotech GmbH, Göttingen, Germany - Supervisory Board
- Gesellschaft für Gemeindeentwicklung und Wirtschaftsförderung Gleichen mbH, Gleichen, Germany - General assembly for community and business development

### Major Activities Beyond Supervisory Board Position

- Chairman of the Employees' Council of Sartorius Stedim Biotech GmbH
- Honorary local political offices: local mayor of Etzenborn; chairman of the parliamentary group in the municipal council of Gleichen
- Member of the committee of the Rhume entertainment association

- Positions held within the meaning of Sec. 125, Subsec. 1, Sentence 5, of the German Stock Corporation Law (AktG)