

## Curriculum Vitae | Dietmar Müller

### Member of the Supervisory Board of Sartorius AG

Year of birth: 1976

Nationality: German

Place of residence: Gleichen, Germany

Member since: May 2020

#### Professional Profile

- Since 1997     Sartorius Group, Göttingen, Germany
- Since 2022: Chairman of the Group Employees' Council of Sartorius AG
- Since 2018: Member and Chairman of the Employees' Council of Sartorius Stedim Biotech GmbH as well as a member of the Group Employees' Council of Sartorius AG; exempted from regular employment duties as a full-time member
- 2016–2018: In-house Sales Representative, BPS Central Europe
- 2012–2016: Team Lead, Commercial Operations Central Europe
- 1999–2012: Customer Service and Support

#### 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

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

- None

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

- Gesellschaft für Gemeindeentwicklung und Wirtschaftsförderung Gleichen mbH, Gleichen, Germany – deputy member of the General Assembly for Community and Business Development of Gleichen

#### Other Major Activities Beyond Supervisory Board Position

- Honorary local political offices: local mayor of Etzenborn; chairman of the parliamentary group in the municipal council of Gleichen