

## Curriculum Vitae | Dominik Langosch

### Member of the Supervisory Board of Sartorius AG

Year of birth: 1984

Nationality: German

Place of residence: Duderstadt, Germany

Member since: April 2025

#### Professional Profile

- |            |   |
|------------|---|
| Since 2021 | German Metalworkers' Union (IG Metall) in southern Lower Saxony-Harz, Northeim, Germany – Managing Director               |
| Since 2016 | IG Metall Coburg/Kassel, Germany – Political secretary  |
| 2014–2015  | Mining, Chemical and Energy Industrial Union (IG BCE), Alsdorf, Germany – Political secretary                             |
| 2008–2011  | Construction-Agriculture-Environment Industrial Union (IG BAU), Dortmund/Frankfurt, Germany – Trainee and union secretary |

#### Educational Background

- Psychotherapist – further education, Göttingen, Germany
- Economic and Social Policy/ Law – Europäische Akademie der Arbeit Frankfurt, Germany
- Psychology/Business Administration – SRH Riedlingen/Hannover, Germany
- Housekeeper – University Medical Center Göttingen, Germany

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

- Demag Cranes & Components GmbH, Wetter, Germany
- Novelis Deutschland GmbH, Göttingen, Germany

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

- None

#### Other Major Activities Beyond Supervisory Board Position

- IG Metall, Lower Saxony and Saxony Anhalt, Germany – Member of the regional commission
- Federal Employment Agency, Göttingen, Germany – Member of the administrative committee