

Curriculum Vitae | Sabrina Wirth

Member of the Supervisory Board of Sartorius AG

Year of birth: 1979

Nationality: German

Place of residence: Nienburg/Weser, Germany

Member since: March 2022

Professional Profile

- | | |
|------------|--|
| Since 2020 | German Metalworkers' Union (IG Metall) Lower Saxony and Saxony-Anhalt – Political secretary for organizational policy in the district management |
| 2016-2020 | IG Metall Nienburg-Stadthagen – Chief authorized representative and treasurer |
| 2008-2016 | IG Metall Nienburg-Stadthagen – Political secretary in the branch office |
| 2007-2008 | IG Metall – Trainee |
| 2005-2007 | University information center (HIB) in Osnabrück – Employee |
| 2003-2007 | IG Metall Leer-Papenburg – Clerk for youth affairs |
| 2001 | KS Gleitlager GmbH, Papenburg – Industrial clerk |

Educational Background

- Social sciences (B.A.) – University of Osnabrück
- Business administration (pre-degree certificate) – University of Osnabrück
- Industrial clerk – KS Gleitlager GmbH, Papenburg

Membership in Other Supervisory Boards Mandated by German Law

- None

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

- None

Other Major Activities Beyond Supervisory Board Position

- None