

Curriculum Vitae | Annette Becker

Member of the Supervisory Board of Sartorius AG

Year of birth: 1964

Nationality: German

Place of residence: Gleichen, Germany

Member since: September 2002

Professional Profile

Since 1995 Sartorius Group, Göttingen, Germany

Since 2014: Member of the Group Employees' Council of Sartorius AG (2014–2022: Chairwoman, since 2022: Deputy Chairwoman)

Since 2008: Member and Chairwoman of the Employees' Council of Sartorius Corporate Administration GmbH; exempted from regular employment duties

1998–2008: Member of the Employees' Council of Sartorius AG (since 2002: Employees' Council member exempted from regular employment duties)

1996–2002: Assistant to the Employees' Council of Sartorius AG

1991–1995 Family phase (two children)

1983–1990 Self-employed (manufacture / sale of craft items, used furniture business, managing director of a moving and short-haul transportation company with car rental business)

Educational Background

- HR Specialist ("Personalfachkauffrau") [German Chamber of Industry and Commerce IHK degree]

Membership in Other Supervisory Boards Mandated by German Law

- None

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

- BKK Technoform, Göttingen, Germany – Member of the Board of Directors

Other Major Activities Beyond Supervisory Board Position

- Honorary judge at the Labor Court of the State of Lower Saxony in Hanover, Germany
- Development and sale of educational games on the subject of labor law and representation of labor interests
- Volunteer consultant at the German Metalworkers' Union Training Center in Berlin, Germany