

Member of the Supervisory Board of Sartorius AG

Annette Becker, resident of Göttingen, Germany
Born 1964

Professional Profile:

- Since 1995 Sartorius Group, Göttingen, Germany
Since 2014: Chairwoman of the Group Employees' Council of Sartorius AG
- 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]

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

- None

Major Activities Beyond Supervisory Board Position:

- Chairwoman of the Employees' Council of Sartorius Corporate Administration GmbH / Group Employees' Council of Sartorius AG; exempted from regular employment duties
- 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 IG Metall-Bildungsstätte in Berlin, Germany