

Member of the Supervisory Board of Sartorius AG

Karoline Kleinschmidt, resident of Hanover, Germany
Born 1965

Professional Profile:

- Since 2006 German Metalworkers' Union IG Metall, District Management of Lower Saxony and Saxony-Anhalt – Union Secretary
- 2001–2005 IG Metall in Berlin – Union Secretary
- 1993–2001 City of Norderstedt, Germany – Employee
- 1990–1993 Stiftung Berufliche Bildung, Foundation for Vocational Training Hamburg, Germany – Educator
- 1988–1989 Flohr Otis, Berlin, Germany – Employee

Educational Background:

- Graduate Social Economist (“Dipl.-Sozialwirtin”) – Hamburg University of Applied Sciences and Politics, Germany
- Energy Electronics Engineer – Flohr Otis, Berlin, Germany

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:

- Union Secretary of the German Metalworkers' Union IG Metall, District Management of Lower Saxony and Saxony-Anhalt