

Member of the Supervisory Board of Sartorius AG

Petra Kirchhoff, resident of Göttingen, Germany
Born 1969

Professional Profile:

- Since 2001 Sartorius Group, Göttingen, Germany
Since 2011: Head of Corporate Communications and Investor Relations
- 2003–2011: Head of Corporate Communications
- Since 2001: Head of Sartorius College
- 1999–2001 Ministry of Science and Culture of Lower Saxony, Hanover, Germany –
Head of Press and PR Activities / Spokesperson and Head of the Minister’s Office
- 1998–1999 Aries M Buchwald e. K., Kassel, Germany – Project Consultant
- 1996–1998 Gesellschaft für Wirtschaftsförderung und Stadtentwicklung Göttingen mbH (subsidiary
of the City of Göttingen for Business Development and Urban Planning), Göttingen,
Germany – Assistant to Executive Management

Educational Background:

- Graduate Political Economist (“Dipl.-Volkswirtin”) - University of Göttingen, Germany

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

- AWO Göttingen gGmbH, Göttingen, Germany – Supervisory Board
- SüdniedersachsenStiftung, Göttingen, Germany – Foundation’s Board of Directors
- Niedersächsische Börse zu Hannover, Hanover, Germany – Stock Exchange Council of Lower Saxony

Major Activities Beyond Supervisory Board Position:

- Head of Corporate Communications and Investor Relations
- Positions held within the meaning of Sec. 125, Subsec. 1, Sentence 5, of the German Stock Corporation Law (AktG)