

Member of the Supervisory Board of Sartorius AG

Dr. Daniela Favoccia, resident of Frankfurt am Main, Germany
Born 1964

Professional Profile:

- Since 1993 Hengeler Mueller, Frankfurt am Main, Germany – Lawyer
Since 1999: Hengeler Mueller, Frankfurt am Main, Germany – Partner at Law
- 2010–2014: Hengeler Mueller, Frankfurt am Main, Germany – Co-Managing Partner at Law
- 1995–1996 Law Office in Los Angeles, California, USA – Lawyer

Educational Background:

- Doctor of Law (“Dr. jur.”) – University of Freiburg, Germany
- Studied law – University of Passau / Freiburg, 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:

- Lawyer and Partner at Law at Hengeler Mueller in Frankfurt am Main, Germany (Specialties: Mergers & acquisitions, corporate law, restructurings)