

Curriculum Vitae | Dr. Daniela Favoccia

Member of the Supervisory Board of Sartorius AG

Year of birth: 1964

Nationality: German

Place of residence: Frankfurt am Main, Germany

Member since: April 2017

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

Membership in Other Supervisory Boards Mandated by German Law

- None

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

- None

Other Major Activities Beyond Supervisory Board Position

- Member of the Government Commission “Regierungskommission” of the German Corporate Governance Code