

## 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

- Freudenberg SE, Weinheim, Germany – Member of the Supervisory Board

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

- Freudenberg & Co. Kommanditgesellschaft, Weinheim, Germany – Member of the Shareholders' Committee

#### Other Major Activities Beyond Supervisory Board Position

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