

Member of the Supervisory Board of Sartorius AG

Uwe Bretthauer, resident of Göttingen, Germany
Born 1958

Professional Profile:

- Since 1988 Sartorius Group, Göttingen, Germany
Since 2014: Member and Chairman of the Employees' Council of Sartorius Lab Instruments GmbH & Co. KG; exempted from regular employment duties
- 2011-2014: Member and Chairman of the Employees' Council of Sartorius Weighing Technology GmbH and of the Group Employees' Council of Sartorius AG; exempted from regular employment duties
- 2002-2011: Member and Chairman of the Employees' Council of Sartorius AG and of the Group Employees' Council of Sartorius AG; exempted from regular employment duties
- 1988-2002: Scientific R&D Engineer
- 1978-1982 Thies Clima, Göttingen, Germany – Electronics Technician

Educational Background:

- Graduate Electrical Engineer ("Dipl.-Ingenieur Elektrotechnik") – University of Applied Technology Gesamthochschule Kassel, Germany
- Energy Electronics Engineer ("Energiegeräteelektroniker") – Robert Bosch GmbH, Göttingen, Germany
- Electronics Engineer ("Elektrogeräteelektroniker") – Robert Bosch GmbH, Göttingen, 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:

- Chairman of the Employees' Council of Sartorius Lab Instruments GmbH & Co. KG; exempted from regular employment duties