

Curriculum Vitae | Dr. Alexandra Gatzemeyer

Member of the Executive Board of Sartorius AG

Head of Lab Products & Services Division

Alexandra Gatzemeyer, born in 1979, is Head of the Lab Products & Services division and has been Member of the Executive Board of Sartorius AG since May 1, 2023.

During her more than 20 years with the company, she has held various senior leadership roles, including Product Management, Human Resources Management, and as Managing Director of a country organization.

Before her appointment to the Executive Board, Alexandra Gatzemeyer led the Sales & Services organization of the Lab Products & Services division at Sartorius. She holds a Ph.D. in chemistry and pharmacology from the St. Petersburg State Chemical and Pharmaceutical University.

Personal Information

Year of birth: 1979

Nationality: German

End of current ordering period: April 2031

Professional Profile

Since 2023	Head of Lab Products & Services Division Member of the Executive Board of Sartorius AG ¹ , Germany
2020-2023	Head of Sales and Service LPS Division, Sartorius Group, Germany
2018-2020	Head of Product Management Lab Essentials, Sartorius Group, Germany
2017-2018	Global HR Business Partner LPS, Sartorius Group, Germany
2012-2017	Managing Director, Sartorius Group, Russia
2011-2017	Managing Director RU / CIS, Sartorius Group, Germany
2009-2011	Market Manager RU CIS, Sartorius Group, Germany
2004-2009	Sales Representative Sartorius Group, Russia

¹ Public listed

Educational Background

- Ph.D. in chemistry and pharmacology from the St. Petersburg State Chemical and Pharmaceutical University, Russia

Other current directorships and positions within the Group

- Vice Chairwoman of the Supervisory Board of Sartorius Corporate Administration GmbH, Germany
- Member of the Board of Directors of Sartorius BioAnalytical Instruments, Inc., USA
- Member of the Board of Directors of Sartorius Corporation, USA
- Member of the Board of Directors of Sartorius Liquid Handling Oy, Finland
- Vice Chairwoman of the Board of Directors of Sartorius Scientific Instruments (Beijing) Co., Ltd., China
- Member of the Board of Directors of Sartorius ForteBio (Shanghai) Co., Ltd., China
- Member of the Board of Directors of Sartorius Lab (Shanghai) Trading Co., Ltd., China
- Member of the Board of Directors of Sartorius (Shanghai) Trading Co., Ltd., China
- Member of the Board of Directors of Sartorius Korea LLC, South Korea
- Chairwoman of the Executive Board Sartorius France S.A.S., France

Other current directorships and positions outside the Group

- Member of the Board of Directors of BICO Group AB¹, Sweden
- Member of the Advisory Board of labforward GmbH, Germany

¹ Public listed