

Curriculum Vitae | 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 18 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: Russian

End of current ordering period: May 2026

Professional Profile

Since May 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
2017-2020	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

Educational Background

¹ Public listed

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

Other current directorships and positions outside the Group

- None

Past directorships (held during the past five years) outside the Group

- None