

Curriculum Vitae | Dr. René Fáber

Member of the Executive Board of Sartorius AG

Head of Bioprocess Solutions Division

René Fáber, born in 1975, has been Head of Bioprocess Solutions Division and Member of the Executive Board of Sartorius AG since 2019. In the same year, he has also became Member of the Board of Directors of Sartorius Stedim Biotech S.A. and was appointed as CEO of the company in 2023.

René Fáber studied chemistry in Bratislava, Slovakia, and earned his Ph.D. in polymer chemistry from the Technical University of Munich in Germany.

He joined Sartorius in 2002, where he initially worked as a scientist on membrane modification. As Vice President, he held various management positions in R&D, Marketing for Filtration and Fermentation Technologies, as well as in Key Account Management. In his most recent position, he was responsible for the Product Development unit of the Bioprocess Solutions Division.

Personal Information

Year of birth: 1975

Nationality: Slovak

End of current ordering period: December 31, 2026

Professional Profile

Since March 27, 2023	CEO of Sartorius Stedim Biotech S.A., France
Since 2019	Head of Bioprocess Solutions Division of Sartorius Group, Member of the Executive Board of Sartorius AG, Germany
2004-2018	Several positions at Sartorius Group (esp. Sartorius Stedim Biotech GmbH, Germany)
2018	Head of Product Development, Bioprocess Solutions Division
2017-2018	Vice President Marketing and Product Management Fermentation Technologies
2016-2018	Key Account Manager Roche/Genentech
2014-2017	Vice President Marketing and Product Management Filtration Technologies
2012-2014	Value Creation Agent in Supplier Relationship Center of Roche and Genentech, San Francisco, USA
2010-2013	Vice President R&D Process Technologies

2006-2010 | Director development and production of surface modified membranes

2004-2006 | Scientist R&D Membrane Modification

2002-2004 Vivascience - Post-doctoral researcher

2001-2002 Rhodia - Scientist at French specialty chemical group, Slovakia

Educational Background

- Master's degree in chemistry at the Technical University in Bratislava, Slovakia
- Doctoral degree in polymer chemistry at the Technical University of Munich, Germany

Other current directorships and positions within the Group

- Chairman of the Executive Board of Sartorius Stedim Biotech S.A.¹, France
- Chairman of the Supervisory Board of Sartorius Stedim Biotech GmbH, Germany
- Chairman of the Advisory Board of Sartorius CellGenix GmbH, Germany
- Member of the Board of Directors of Sartorius Stedim (Shanghai) Trading Co., Ltd., China
- Vice Chairman of the Board of Directors of Sartorius Stedim Biotech (Beijing) Co., Ltd., China
- Member of the Board of Directors of Albumedix Ltd., United Kingdom
- Member of the Board of Directors of Sartorius Stedim BioOutsource Ltd., Scotland, UK
- Member of the Board of Directors of Sartorius Korea Biotech LLC, South Korea
- Member of the Board of Directors of Sartorius Korea Operations LLC, South Korea
- Member of the Board of Directors of Sartorius Stedim North America Inc., USA
- Chairman of the Executive Committee of Sartorius Stedim FMT S.A.S., France
- Chairman of the Advisory Board of Sartorius BIA Separations d.o.o., Slovenia

Other current directorships and positions outside the Group

- Member of the Advisory Board of Curexsys GmbH, Germany

¹ Public listed