

Curriculum Vitae | Dr. René Fáber

Chief Executive Officer of Sartorius Stedim Biotech S.A.

Year of Birth: 1975

Nationality: Slovak

First appointment to the Board of Directors: March 2019

End of term: 2029 Shareholders' Meeting

Professional Profile

Since 2023	CEO of Sartorius Stedim Biotech S.A. ¹ , France
2022 – 2023	Deputy CEO of Sartorius Stedim Biotech S.A. ¹ , France
Since 2019	Head of Bioprocess Solutions Division of Sartorius Group, Member of the Executive Board of Sartorius AG ¹ , Germany
2004-2018	Various 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

¹ Public listed

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

- Member of the Executive Board of Sartorius AG¹, Germany
- 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 Albumedix Ltd., United Kingdom
- Member of the Board of Directors of Sartorius Stedim BioOutsource Ltd., United Kingdom
- Member of the Board of Directors of Sartorius Stedim North America Inc.
- 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 (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 Polyplus-Transfection S.A., France
- President and 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

- None

¹ Public listed