

Curriculum Vitae | Romaine Fernandes

Member of the Board of Directors of Sartorius Stedim Biotech S.A.

Director Representing Employees

Year of Birth: 1969

Nationality: French

First appointment to the Board of Directors: October 2023

End of term: 2026 Shareholders' Meeting

Professional Profile

2019-2023	Sartorius Stedim Biotech, France - Purchasing Platform and Insurance Expert
2016-2019	Sartorius Stedim Biotech, France - Central Purchasing Manager
2014-2016	Sartorius Stedim Biotech, France - Assistant Manager in General Services
2009-2010	Biopharm Services, United Kingdom - Assistant Manager
2003-2014	Sartorius Stedim Biotech, France - Administration and Reception
1990-1995	Cathay Pacific Airways, Hong Kong - Stewardess
1990	Hotel Oberoi Sheraton, India - Cashier & Accounting

Educational Background

- Bachelor of Commerce at the University of Mumbai, India
- Diploma in Tourism, Bombay, India

Other current directorships and positions outside the Group

- None