

## Curriculum Vitae | Cécile Dussart

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

Year of Birth: 1964

Nationality: French

First appointment to the Board of Directors: March 2025

End of term: 2028 Shareholders' Meeting

### Professional Profile

Since 2022	EUROAPI S.A. <sup>1</sup> , France – Board Member and Chair of the ESG Committee NEOBRAIN S.A.S., France – Member of the Advisory Committee HEC PARIS and ESCP Business School – Keynote Speaker
2020-2023	Galderma, France – Global Head of Operations and ESG
2013-2019	Galderma, France – Global Head of Operations
2009-2012	Galderma, France – Plant Director
2006-2008	Galderma, France – Human Resources Director
2001-2006	Galderma, France – Human Resources Manager
1997-2000	Galderma, France – Senior Global Brand Manager
1990-1996	Sanofi, France – Global Brand Manager

### Educational Background

- Doctor of Pharmacy (Pharm.D.) from Paris-Sud University
- Master's Degree in Pharmaceutical Marketing

### Other current directorships and positions outside the Group

- Independent Director and Chair, ESG Committee at EUROAPI<sup>1</sup>
- Member of the Advisory Committee at NEOBRAIN

<sup>1</sup> Public listed