

## Curriculum Vitae | Anne-Marie Graffin

Member of the Board of Directors of Sartorius Stedim Biotech S.A.  
Lead Independent Director

Year of Birth: 1961

Nationality: French

First appointment to the Board of Directors: April 2015

End of term: 2027 Shareholders' Meeting

### Professional Profile

Since 2011	Independent Non-Executive Board Member, Life Sciences Expert and Advisor
2011-2024	SMAG Consulting S.A.S., France - President
2009-2010	Sanofi Pasteur MSD, France - Executive Vice President
2006-2008	Sanofi Pasteur MSD, France - Executive Director Business Management
2001-2005	Sanofi Pasteur MSD, France - Head of range then Adult Vaccines Marketing Director Europe
1998-2000	Sanofi Pasteur MSD, France - Product Manager Adult Vaccines France
1991-1995	RoC S.A (Johnson & Johnson), France - Head of International Marketing Group
1988-1990	URGO Laboratories, France - Marketing Manager
1984-1987	International Distillers and Vinters, France - France Products Manager

### Educational Background

- Graduated from ESSEC Business School, France

### Other current directorships and positions outside the Group

- Chairwoman of the Board of Directors of Valneva SE<sup>1</sup>, France
- Member of the Supervisory Board of Nanobiotix S.A.<sup>1</sup>, France
- Member of the Board of Directors of Vetoquinol S.A.<sup>1</sup>, France

<sup>1</sup> Listed company