

## Curriculum Vitae | Henri Riey

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

Year of Birth: 1961

Nationality: Monegasque

First appointment to the Board of Directors: June 2007

End of term: 2025 Shareholders' Meeting

### Professional Profile

2004-2013	Hendyplan S.A. <sup>1</sup> , Beglium, CFO
1999-2004	Barclay Bank, France - Executive Vice President of Barclays Asset Management; in charge of all fund management businesses
1996-1999	Barclay Bank, France - Head of Research of Barclays Asset Management Europe
1988-1996	Barclays Bank, France - Fund Manager, responsible for the European Equity Fund Management Team
1985-1988	BNP Paribas, France - Fund Manager

### Educational Background

- Diplôme Institut Supérieur de Gestion, France (degree earned at the French Higher Institute of Business Management "Institut supérieur de gestion")

### Other current directorships and positions outside the Group

- President of Aidea
- President of Groupe HR S.A.S., France
- President of Association Monegasque de Cindynique
- Director and secretary-treasurer of The Princess Grace Foundation, Monaco

<sup>1</sup> Public listed