

## Curriculum Vitae | Christopher Nowers

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

Year of Birth: 1963

Nationality: British

First appointment to the Board of Directors: March 2025

End of term: 2028 Shareholders' Meeting

### Professional Profile

|           |   |
|-----------|---|
| 2020-2025 | ONK Therapeutics, USA – CEO and Board Member  |
| 2018-2020 | Cell Medica, UK – CEO and Board Member  |
| 2016-2018 | Kite Pharma Europe, UK – Head of Europe   |
| 2011-2016 | Bristol Myers Squibb, USA – VP Early Oncology, VP Global Commercialisation Yervoy, Head of Immunology and Hematology France |
| 2009-2011 | Genomic Health, USA – VP Sales and Marketing  |
| 2007-2009 | Independent Biopharma Consultant, USA   |
| 2006-2007 | Avantogen Oncology, USA – CEO and Board Member  |
| 2003-2006 | Amgen, USA – Global Commercial Leader Neurology, Brand Director Nephrology  |
| 1994-2003 | Amgen Europe, Switzerland – Positions in Oncology, Hematology, and Rheumatology   |
| 1987-1994 | Zeneca Pharma, UK – Positions in Sales, Marketing, and Product Management   |
| 1986-1987 | The Wellcome Foundation, UK – Sales Representative  |

### Educational Background

- Bachelor of Science Hons degree in Biochemistry, University of Kent, UK

### Other current directorships and positions outside the Group

- None