

## Curriculum Vitae | Pascale Boissel

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

Year of birth: 1966

Nationality: French

First appointment to the Board of Directors: March 2019

End of term: 2025 Shareholders' Meeting

### Professional Profile

- |           |   |
|-----------|---|
| 2017-2021 | NOVADISCOVERY – Part time Chief Financial Officer, France                             |
| 2017-2018 | ENYO PHARMA – Part time Chief Financial Officer, France                               |
| 2012-2016 | BIOASTER Institute – Chief Financial Officer & Deputy Chief Executive Officer, France |
| 2009-2012 | IPSOGEN – Chief Financial Officer, France   |

### Educational Background

- Graduated from HEC (Ecole des hautes Etudes de Commerciales): MBA in Finance & Audit  
Graduated with a CPA diploma (diplôme d'expertise comptable & commissariat aux comptes)

### Other current directorships and positions outside the Group

- Member of the Board of Directors of Poxel S.A.<sup>1</sup>, France
- Member of the Supervisory Board of Innate Pharma S.A.<sup>1</sup>, France

<sup>1</sup> Public listed