

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: 2029 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.¹, France
- Member of the Supervisory Board of Innate Pharma S.A.¹, France

¹ Public listed