

## Curriculum Vitae | Abdelatif Zeroual

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

Director Representing Employees

Year of Birth: 1985

Nationality: French

First appointment to the Board of Directors: March 2026

End of term: 2029 Shareholders' Meeting

### Professional Profile

Since 2023	Sartorius Stedim Biotech, France - IT Engineering Expert
2015-2023	Sartorius Stedim Biotech, France - End User Support Specialist EMEA
2015	EDF CIPN, France - IT Technician
2015	Econocom-Osiastis, France - Bilingual English Support Technician
2014	Technicien Center de Service Techniques et Logiciels, France - General Council of the Jury
2014	Computacenter, France - IT Technician

### Educational Background

- Bachelor's degree in Sales, Marketing and Communication, University of Burgundy
- Advanced Technician Diploma in IT Systems and Network Administration, Afpa

### Other current directorships and positions outside the Group

- None