

# Curriculum Vitae | Dr. Florian Funck

Member of the Executive Board of Sartorius AG  
Chief Financial Officer

Florian Funck, born 1971, has been a Member of the Executive Board of Sartorius AG and CFO since April 1, 2024.

Funck studied business administration at the University of Münster where he also earned his doctorate. Before joining Sartorius, he was with Haniel Group for over two decades, where he held several leadership roles: In 2004, he was appointed as CFO to the Management Board of TAKKT AG, a publicly listed company where Haniel is the majority shareholder. Between 2011 and 2023, he was CFO of Haniel Group.

He is a member of the supervisory board of Vonovia SE, and honorary professor at the University of Münster.

**Personal Information**

Year of birth: 1971  
Nationality: German  
End of current ordering period: March 31, 2027

**Professional Profile**

Since April 2024	Member of the Executive Board of Sartorius AG <sup>1</sup> , Germany
2011- 2024	Franz Haniel & Cie. GmbH - CFO, Germany
2004-2011	TAKKT AG <sup>1</sup> – Management Board Member (CFO), Germany
1999-2004	Franz Haniel & Cie. GmbH - Head of Corporate Controlling, Strategic Planning, Group Accounting and Risk Management, Germany
1994-1998	Research Assistant at the Institute for Industrial Management, University of Münster, Germany

**Educational Background**

- Dr. rer. pol., Business Administration, University of Münster, Germany

<sup>1</sup> Publicly listed

#### Other current directorships and positions within the Group

- Chairman of the Supervisory Board of Sartorius Corporate Administration GmbH, Germany

#### Other current directorships and positions outside the Group

- Member of the Supervisory Board of Vonovia SE<sup>1</sup>