

## Curriculum Vitae | Hermann Jens Ritzau

### Member of the Supervisory Board of Sartorius AG

Year of birth: 1966

Nationality: German

Place of residence: Katlenburg-Lindau, Germany

Member since: March 2021

#### Professional Profile

Since 1989     Sartorius Group, Göttingen, Germany

Since March 2020: Member and Chairman of the Employees' Council of Sartorius Lab Instruments GmbH & Co. KG as well as a member of the Group Employees' Council of Sartorius AG; exempted from regular employment duties as a full-time member

2002–2020: Member of Employees' Council of Sartorius AG, Sartorius Weighing Technology GmbH and of Sartorius Lab Instruments GmbH & Co. KG

1996–2020: Deputy Representative and Principal Representative of the Committee for Employees with Disabilities at Sartorius AG, Sartorius Weighing Technology GmbH and at Sartorius Lab Instruments GmbH & Co. KG

1989–2020: Employee in the departments of logistics, materials planning and services

#### Membership in Other Supervisory Boards Mandated by German Law

- None

#### Membership in Comparable Supervisory Committees of Business Enterprises Within Germany and Abroad

- None

#### Other Major Activities Beyond Supervisory Board Position

- Member of the pay scale bargaining committee for the German metal and electrical industries in Lower Saxony, Osnabrück-Emsland-Grafschaft Bentheim and Saxony-Anhalt, as well as for Volkswagen AG
- Honorary judge at the Labor Court of Göttingen, Germany