

Curriculum Vitae | Manfred Zaffke

Member of the Supervisory Board of Sartorius AG

Year of birth: 1957

Nationality: German

Place of residence: Osterode am Harz, Germany

Member since: March 2014

Professional Profile

- | | |
|------------|--|
| Since 2022 | German Metalworkers' Union IG Metall in southern Lower Saxony-Harz – Project Secretary with special duties |
| 2008–2022 | German Metalworkers' Union IG Metall in southern Lower Saxony-Harz – Managing Director |
| 1992–2008 | IG Metall in Goslar-Osterode, Germany – Managing Director |
| 1988–1992 | IG Metall in Osterode, Germany – Union Secretary |
| 1984–1987 | IG Metall in Siegburg, Germany – Union Secretary |
| 1976–1980 | Kamax GmbH & Co. KG – Toolmaker |

Educational Background

- Graduate Political Economist (“Dipl.-Volkswirt”) – Hamburg University of Applied Sciences and Politics, Germany
- Toolmaker – Kamax GmbH & Co. KG

Membership in Other Supervisory Boards Mandated by German Law

- Demag Cranes & Components GmbH, Düsseldorf, Germany – member of the Supervisory Board
- Konecranes Holding GmbH, Mannheim, Germany – member of the Supervisory Board

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

- None

Other Major Activities Beyond Supervisory Board Position

- Honorary member of the Executive Board of the German Metalworkers' Union IG Metall
- Shareholder of IGETMET