

## Member of the Supervisory Board of Sartorius AG

Manfred Zaffke, resident of Osterode am Harz, Germany

Born 1957

### Professional Profile:

Since 2008	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

### Positions Held Within the Meaning of Sec. 125, Subsec. 1, Sentence 5, of the German Stock Corporation Law (AktG):

- Demag Cranes & Components GmbH, Düsseldorf, Germany – Supervisory Board
- GMH GUSS GmbH, Schwerte, Germany – Supervisory Board (Vice Chairman)
- Konecranes Holding GmbH, Mannheim – Supervisory Board

**Major Activities Beyond Supervisory Board Position:**

- Managing Director of the German Metalworkers' Union IG Metall in southern Lower Saxony-Harz
- Positions held within the meaning of Sec. 125, Subsec. 1, Sentence 5, of the German Stock Corporation Law (AktG)