

Individual Meeting Attendance of the Supervisory Board Members 2020

Supervisory Board Member	Meetings	Executive Task Committee	Audit Committee
Lothar Kappich	7 / 7	6 / 6	4 / 4
Manfred Zaffke	7 / 7	6 / 6	4 / 4
Annette Becker	6 / 7		
Uwe Bretthauer	7 / 7	6 / 6	4 / 4
Michael Dohrmann (until May 15, 2020)	4 / 4		
David Raymond Ebsworth	7 / 7		
Daniela Favoccia	7 / 7		
Petra Kirchhoff	7 / 7		
Karoline Kleinschmidt	7 / 7		
Dietmar Müller (since May 16, 2020)	2 / 3		
Ilke Hildegard Panzer	7 / 7		
Thomas Scheper	7 / 7		
Klaus Rüdiger Trützschler	7 / 7	5 / 6	4 / 4

The Nomination Committee and Conciliation Committee did not meet in 2020.