

Individual Meeting Attendance of the Supervisory Board Members 2021

Supervisory Board Member	Meetings	Executive Task Committee	Audit Committee	Nomination Committee
Lothar Kappich	8 / 8	4 / 4	4 / 4	5/5
Manfred Zaffke	8 / 8	3 / 4	4 / 4	
Annette Becker (on the Executive Committee since March 1, 2021)	8 / 8	3/4		
Uwe Bretthauer (until February 28, 2021)	1 / 8	1 / 4	1 / 4	
David Raymond Ebsworth	8 / 8			
Daniela Favoccia	8 / 8			4 / 5
Petra Kirchhoff	8 / 8			
Karoline Kleinschmidt	8 / 8			
Dietmar Müller (on the Audit Committee since March 1, 2021)	8 / 8		3/4	
Ilke Hildegard Panzer	8 / 8			
Hermann Jens Ritzau (since March 1, 2021)	7 / 8			
Thomas Scheper	7 / 8			
Klaus Rüdiger Trützscher	8 / 8	4 / 4	4 / 4	5/5

The Conciliation Committee did not meet in 2021.