

Individual Meeting Attendance of the Supervisory Board Members 2022

Supervisory Board Members	Supervisory Board Meetings	Executive Committee	Audit Committee
Lothar Kappich	6/6	3/3	4/4
Manfred Zaffke	6/6	3/3	4/4
Annette Becker	6/6	3/3	
David Raymond Ebsworth	6/6		
Daniela Favoccia	6/6		
Petra Kirchhoff	6/6		
Karoline Kleinschmidt (until March 25, 2022)	1/6		
Dietmar Müller	6/6		4/4
Ilke Hildegard Panzer	6/6		
Frank Riemensperger (until March 25, 2022)	5/6		
Hermann-Jens Ritzau	6/6		
Thomas Scheper (until March 25, 2022)	1/6		
Klaus Rüdiger Trützschler	6/6	3/3	4/4
Sabrina Wirth (until March 25, 2022)	5/6		

The Nominations and Conciliation Committees did not meet in 2022.