

Individual Meeting Attendance of the Supervisory Board Members 2024

Supervisory Board	Meetings	Executive Committee	Audit Committee	Nomination Committee
Lothar Kappich	8/8	10/10	4/4	2/2
Manfred Zaffke	8/8	10/10	4/4	
Annette Becker	8/8	10/10		
David Raymond Ebsworth	8/8			
Daniela Favoccia	8/8			1/2
Petra Kirchhoff	8/8			
Dietmar Müller	8/8		4/4	
Ilke Hildegard Panzer	8/8			
Frank Riemensperger	8/8			
Hermann-Jens Ritzau	8/8			
Klaus Rüdiger Trützschler	8/8	10/10	4/4	2/2
Sabrina Wirth	8/8			

The Conciliation Committee did not meet in 2024.