

## Individual Meeting Attendance of the Supervisory Board Members in 2025

Supervisory Board	Meetings	Executive Committee	Audit Committee	Nomination Committee	Virtual Meetings
Lothar Kappich	6/6	4/4	4/4	8/8	5/22
Manfred Zaffke until March 31, 2025	2/6	2/4	1/4	--	2/5
Dominik Langosch from April 2, 2025	4/6	--	3/4	--	4/7
Annette Becker	6/6	4/4	--	--	3/10
David Raymond Ebsworth	6/6	--	--	--	2/6
Daniela Favoccia	6/6	--	--	8/8	1/14
Beatrix Henseler from September 2, 2025	2/6	--	--	--	0/2
Petra Kirchhoff until July 31, 2025	4/6	--	--	--	1/4
Dietmar Müller	6/6	2/4	4/4	--	4/12
Ilke Hildegard Panzer	6/6	--	--	--	1/6
Frank Riemensperger	6/6	--	--	--	1/6
Hermann-Jens Ritzau	6/6	--	--	--	1/6
Klaus Rüdiger Trützschler	6/6	4/4	4/4	8/8	7/22
Sabrina Wirth	3/6	--	--	--	1/3

The Conciliation Committee did not meet in 2025.