

Shareholder representatives

Member since

Independence (acc. DCGK C.6)

Overboarding (acc. DCGK C.4)

Gender

Year of birth

Nationality

Educational background

L. Kappich	D. Ebsworth	D. Favoccia	I. Panzer	F. Riemensperger	K. Trützschler
April 2007	January 2020	April 2017	April 2017	March 2022	Nov. 2011
no ¹	yes	yes	yes	yes	yes
no	no	no	no	no	no
male	male	female	female	male	male
1957	1954	1964	1966	1962	1948
German	British	German	German USA	German	German
Economics	Pharma	Law	Engineering	Computer Science	Mathematics

Corporate Governance; Compliance

Strategy development; Portfolio management; M&A

Industry know-how and customer-specific perspectives

Technology and product development

Capital Markets

Finance; Accounting and Financial Statement Audit;

ESG regulation

Human Resources Management;

Employee safety and engagement

Digital transformation; IT

Leadership in international enterprises; C-level experience

Environmental impact; Consumption of resources

International experience or personal background

■	■	■	■	■	■
■	■	■	■	■	■
	■		■		
	■	■	■	■	■
■	■	■			■
■	■		■	■	■
			■	■	
■			■		■
■			■		■
	■	■	■	■	

Employee representatives

Corporate Governance; Compliance

Strategy development; Portfolio management; M&A

Industry know-how and customer-specific perspectives

Technology and product development

Capital Markets

Finance; Accounting and Financial Statement Audit;

ESG regulation

Human Resources Management;

Employee safety and engagement

Digital transformation; IT

Leadership in international enterprises; C-level experience

Environmental impact; Consumption of resources

International experience or personal background

A. Becker	B. Henseler	D. Langosch	D. Müller	H. Ritzau	S. Wirth
		■			■
	■	■			■
			■	■	
		■	■		
■	■	■	■	■	■
■	■			■	
	■				
■					
	■				