

	L. Kappich	M. Zaffke	A. Becker	D. Ebsworth	D. Favoccia	P. Kirchhoff
Corporate governance and strategy development	●	●		●	●	
Customer-specific perspectives				●		
Technology and product development				●		
International markets				●		●
Financial economy	●	●		●		●
Corporate and capital market law					●	●
Human resources	●	●	●	●		
Employee-specific perspectives		●	●			●
Digitalization			●			
Competence sustainability and regulatory affairs	●				●	●
International experience or life backgrounds				●	●	
	D. Müller	I. Panzer	H. Ritzau	K. Trützscher	F. Riemensperger	S. Wirth
Corporate governance and strategy development		●		●	●	●
Customer-specific perspectives	●	●	●			
Technology and product development		●			●	
International markets		●			●	
Financial economy	●			●		
Corporate and capital market law				●		
Human resources		●		●	●	●
Employee-specific perspectives	●		●			●
Digitalization		●	●		●	
Competence sustainability and regulatory affairs				●		
International experience or life backgrounds		●			●	