

Competence Profile

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
Environmental impact; Consumption of resources
International experience or personal background
Leadership in international enterprises; C-level experience

M. Grosse	R. Fáber	P. Boissel	S. Dexter	C. Dussart
		■		■
■	■	■	■	■
	■		■	■
■	■		■	
■	■	■	■	
		■	■	■
■	■	■	■	■
■				
			■	■
■	■	■	■	■
■	■	■	■	■

Competence Profile

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
Environmental impact; Consumption of resources
International experience or personal background
Leadership in international enterprises; C-level experience

R. Fernandes	A. Graffin	L. Kappich	C. Nowers
	■	■	
	■	■	■
■			■
			■
	■		
■	■	■	
■	■	■	■
■	■	■	
	■		■
	■	■	■