

Competence Profile 2024

Corporate Governance
Strategy development
Customer-specific perspectives
Technology and product development
International markets
Finance, audit and risk management
Corporate and capital market law
Employee-specific perspectives
Digitalization
Sustainability and ESG

Corporate Governance
Strategy development
Customer-specific perspectives
Technology and product development
International markets
Finance, audit and risk management
Corporate and capital market law
Employee-specific perspectives
Digitalization
Sustainability and ESG

Joachim Kreuzburg	Pascale Boissel	Susan Dexter	René Fáber
■	■		
■	■	■	■
		■	■
■		■	■
■		■	■
■	■	■	
	■	■	
		■	■
■		■	

Romaine Fernandes	Anne-Marie Graffin	Lothar Kappich	Henri Riey
	■	■	■
	■	■	■
■			
■	■	■	■
	■		■
■	■	■	
			■
			■