

Key Figures for the First Half and Second Quarter of 2021

in millions of € unless otherwise specified	6 months 2021	6 months 2020 ¹	Δ in %	Δ in % cc ²
Order Intake and Sales Revenue				
Order intake	1,889.9	1,048.7	80.2	87.4
Sales revenue	1,351.9	869.1	55.6	61.1
▪ EMEA ³	580.1	340.3	70.5	71.3
▪ Americas ³	419.8	312.6	34.3	46.1
▪ Asia Pacific ³	352.0	216.2	62.8	66.1
Results				
EBITDA ⁴	487.4	262.3	85.8	
EBITDA margin ⁴ in %	36.1	30.2		
Net profit ⁵	321.4	167.4	91.9	
Earnings per share ⁵ in €	3.49	1.82	91.9	
	June 30, 2021	Dec. 31, 2020		
Balance Sheet Financials				
Balance sheet total	3,521.9	3,065.5		
Equity	1,693.1	1,482.9		
Equity ratio in %	48.1	48.4		
Net debt	355.2	527.3		
Ratio of net debt to underlying EBITDA	0.4	0.8		

in millions of € unless otherwise specified	Q2 2021 ⁶	Q2 2020 ^{1,6}	Δ in %	Δ in % cc ²
Order Intake and Sales Revenue				
Order intake	885.5	514.0	72.3	78.7
Sales revenue	696.7	447.0	55.9	61.1
▪ EMEA ³	296.0	172.7	71.4	
▪ Americas ³	215.8	159.9	35.0	
▪ Asia Pacific ³	184.9	114.4	61.6	
Results				
EBITDA ⁴	255.8	135.6	88.6	
EBITDA margin ⁴ in %	36.7	30.3		
Net profit ⁵	170.2	86.8	96.2	
Earnings per share ⁵ in €	1.85	0.94	96.2	

1 The previous year's figures have been restated due to finalization of the purchase price allocations for acquisitions of 2020

2 cc = in constant currencies

3 According to customer location

4 Underlying EBITDA = earnings before interest, taxes, depreciation and amortization, and adjusted for extraordinary items

5 Underlying net profit = net profit after non-controlling interest; adjusted for extraordinary items, amortization and based on a normalized financial result and normalized tax rate

6 Figures are not audited nor reviewed