

Key Figures for the First Half of 2021

in millions of € unless otherwise specified	Group				Bioprocess Solutions				Lab Products & Services			
	6-mo. 2021	6-mo. 2020 ¹	Δ in %	Δ in % cc ²	6-mo. 2021	6-mo. 2020 ¹	Δ in %	Δ in % cc ²	6-mo. 2021	6-mo. 2020 ¹	Δ in %	Δ in % cc ²
Sales Revenue and Order Intake												
Order intake	2,178.8	1,244.8	75.0	82.4	1,803.4	984.9	83.1	91.0	375.5	259.9	44.5	49.9
Sales revenue	1,629.2	1,056.8	54.2	60.1	1,266.5	809.3	56.5	62.6	362.7	247.5	46.6	52.0
▪ EMEA ³	682.5	419.5	62.7	63.8	539.2	310.9	73.4	74.3	143.3	108.6	31.9	33.2
▪ Americas ³	519.0	373.1	39.1	52.0	405.2	304.8	32.9	45.3	113.9	68.3	66.6	82.3
▪ Asia Pacific ³	427.6	264.2	61.9	65.5	322.1	193.7	66.3	70.2	105.5	70.5	49.7	52.7
Results												
EBITDA ⁴	555.3	293.5	89.2		460.2	247.2	86.2		95.1	46.4	105.1	
EBITDA-Margin ⁴ in %	34.1	27.8			36.3	30.5			26.2	18.7		
Net result ⁵	259.4	124.3	108.7									
Financial data per share												
Earnings per ordinary share ⁵ in €	3.79	1.81	108.9									
Earnings per preference share ⁵ in €	3.80	1.82	108.3									
	June 30, 2021	Dec. 31, 2020 ¹										
Balance Sheet Financials												
Balance sheet total	5,213.5	4,693.5										
Equity	1,630.8	1,402.2										
Equity ratio in %	31.3	29.9										
Net debt	1,698.4	1,883.9										
Net debt to underlying EBITDA	1.8	2.6										

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

² In constant currencies

³ Acc. to customers' location

⁴ Adjusted for extraordinary items

⁵ After non-controlling interest, adjusted for extraordinary items and non-cash amortization, as well as based on the normalized financial result and the normalized tax rate

Key Figures for the Second Quarter of 2021

in millions of € unless otherwise specified	Group				Bioprocess Solutions				Lab Products & Services			
	Q2 2021 ¹	Q2 2020 ^{1,2}	Δ in %	Δ in % cc ³	Q2 2021 ¹	Q2 2020 ^{1,2}	Δ in %	Δ in % cc ³	Q2 2021 ¹	Q2 2020 ^{1,2}	Δ in %	Δ in % cc ³
Sales Revenue and Order Intake												
Order intake	1,040.2	615.4	69.0	75.6	850.3	478.8	77.6	84.7	189.9	136.6	39.0	43.6
Sales revenue	838.1	546.9	53.2	58.7	655.7	415.0	58.0	63.7	182.3	131.9	38.3	43.1
▪ EMEA ⁴	348.1	212.5	63.8		279.0	157.1	77.6		69.2	55.4	24.9	
▪ Americas ⁴	265.9	192.5	38.1		208.1	155.0	34.2		57.8	37.4	54.4	
▪ Asia Pacific ⁴	224.1	141.9	57.9		168.7	102.9	64.0		55.4	39.0	41.8	
Results												
EBITDA ⁵	291.8	155.7	87.4		243.1	127.3	90.9		48.7	28.4	71.6	
EBITDA-Margin ⁵ in %	34.8	28.5			37.1	30.7			26.7	21.5		
Net result ⁵	137.9	66.9	106.1									
Financial Data per Share												
Earnings per ordinary share ⁶ in €	2.02	0.98	106.0									
Earnings per preference share ⁶ in €	2.02	0.98	106.0									

1 Figures are not audited nor reviewed

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

3 In constant currencies

4 Acc. to customers' location

5 Adjusted for extraordinary items

6 After non-controlling interest, adjusted for extraordinary items and non-cash amortization, as well as based on the normalized financial result and the normalized tax rate