

Key Figures

All figures are stated in millions of € according to IFRS, unless otherwise specified

	2024	Δ in %	2023	2022	2021	2020
Order intake, sales revenue and earnings						
Order intake ¹	2,781.6	12.3	2,476.1	3,314.8	3,664.4	2,381.0
Sales revenue	2,780.0	0.2	2,775.5	3,492.7	2,887.0	1,910.1
Underlying EBITDA ^{2,3}	779.0	-0.8	785.4	1,221.4	1,033.4	604.7
Underlying EBITDA ^{2,3} as % of sales revenue	28.0	-0.3pp	28.3	35.0	35.8	31.7
Net profit after non-controlling interest	175.1	-43.6	310.3	876.1	414.4	335.9
Underlying net profit after non-controlling interest ⁴	337.5	-12.5	385.9	796.6	687.8	383.8
Research and development costs	144.1	11.3	129.5	132.4	110.5	84.5
Financial data per share						
Earnings per share (in €)	1.81	-46.3	3.37	9.51	4.50	3.64
Underlying Earnings per share (in €) ⁴	3.49	-16.7	4.19	8.64	7.46	4.16
Dividend per share (in €)	0.69 ⁵	0.00	0.69	1.44	1.26	0.68
Balance sheet						
Balance sheet total	8,256.4	6.8	7,730.1	5,065.4	3,951.1	2,856.7
Equity	4,023.8	50.5	2,673.8	2,514.2	1,733.2	1,461.0
Equity ratio (in %) ⁶	48.7	14.1pp	34.6	49.6	43.9	51.1
Financials						
Capital expenditures as % of sales revenue	12.2	-4.9pp	17.1	12.3	11.2	8.3
Depreciation and amortization	301.7	27.4	236.8	179.9	141.5	100.3
Cash flow from operating activities ⁷	815.1	9.2	746.4	612.3	701.9	416.9
Net debt ⁸	2,190.6	-38.6	3,565.2	1,028.6	401.9	527.3
Ratio of net debt to underlying EBITDA ^{2,3,9}	2.8		4.5	0.8	0.4	0.8
Total number of employees as of December 31						
	9,901	-7.1	10,662	11,934	10,409	7,566

1 All customer orders contractually concluded and booked during the respective reporting period.

2 Earnings before interest, taxes, depreciation, and amortization and adjusted for extraordinary items.

3 For more information on EBITDA, net profit, and the underlying presentation, please refer to the Group Business Development chapter and to the Glossary.

4 Profit for the period after non-controlling interest, adjusted for extraordinary items and amortization, as well as based on the normalized financial result and the normalized tax rate.

5 Amount suggested by the Board of Directors (Conseil d'administration) and subject to approval by the Annual General Shareholders' Meeting.

6 Equity in relation to the balance sheet total.

7 Interest received are reported under cash flows from operating activities since fiscal 2022. The prior year figure was restated accordingly.

8 The net debt excludes the liability for the remaining purchase price for acquisitions; 2024: €79.6 million, 2023: €80.6 million, 2022: €245.1 million, 2021: €518.7 million, 2020: €127.8 million.

9 Quotient of net debt and underlying EBITDA over the past 12 months, including the pro forma amount contributed by acquisitions for this period.

The figures for the financial year 2023 have been revised due to finalization of the purchase price allocation for Polyplus.