



Sartorius Group
First-Quarter Report
January to March 2009

Business Development and Outlook

- **Growth trend in the Biotechnology Division continues**
- **EBITA margin for the Biotechnology Division significantly increases from 10.0% to 13.0%**
- **Global recession leads to sharp slump in demand and negative earnings contribution for the Mechatronics Division**
- **Solid Group operating earnings**
- **Comprehensive package of cost-cutting measures planned for the Mechatronics Division**

Order Intake

For the Biotechnology Division, order intake from January to March 2009 grew 6.9% (currency-adjusted: +3.9%) from €97.6 million a year ago to €104.4 million. In particular, business with single-use products for the biopharmaceutical industry fueled this gain at clearly double-digit growth rates. The dynamic growth trend of the last months of 2008 thus continued on into the first quarter of 2009. In North America, where measures for inventory reduction on the part of our customers were largely completed by the end of 2008, demand even increased by more than a third compared with the year-earlier quarter. In Europe and Asia | Pacific, business with single-use products also grew; however, overall order intake fell in this region. This can be attributed to the exceptionally high order intake for equipment business in the previous year.

In the Mechatronics Division, we reported a sharp slump in demand due to the global recession. For this division, first-quarter order intake slid to €52.6 million from €64.9 million a year ago. This is a minus of 18.9% (currency-adjusted: -21.8%). During the past months, our customers primarily in the chemical industry drastically cut back their orders. The global decline in demand affected all business areas, except for service business, which saw a moderate gain.

Overall, the Sartorius Group received orders valued at €157.0 million (previous year: €162.5 million). Compared with a year earlier, this signifies a decrease of 3.4% (currency-adjusted: -6.3%).

Sales Revenue

Sales revenue of the Biotechnology Division rose 6.6% from €90.6 million to €96.6 million. On the basis of constant currencies, sales revenue grew 4.1%. This growth was driven by products in the segment of single-use technology. We reported the highest rates of increase in North America and Asia | Pacific.

The cyclical slump in demand in the Mechatronics Division is reflected accordingly in the development of its sales revenue. During the first quarter, the division's sales revenue fell 15.6% (currency-adjusted: -18.6%) from €58.6 million to €49.5 million.

At the Group level, first-quarter sales revenue was at €146.0 million, up from €149.2 million a year ago. Thus, consolidated sales revenue dipped 2.1% relative to the previous year (currency adjusted: -4.8%).



Key Figures

€ in millions (unless otherwise specified)	Group				Biotechnology				Mechatronics			
	1st quarter 2009	1st quarter 2008	Change in %	Change in % (const. currencies)	1st quarter 2009	1st quarter 2008	Change in %	Change in % (const. currencies)	1st quarter 2009	1st quarter 2008	Change in %	Change in % (const. currencies)
Sales revenue	146.0	149.2	-2.1	-4.8	96.6	90.6	6.6	4.1	49.5	58.6	-15.6	-18.6
- Europe ¹⁾	85.0	92.3	-8.0	-6.3	56.7	57.6	-1.6	0.3	28.3	34.7	-18.6	-17.4
- North America ¹⁾	30.7	27.6	11.1	-3.2	23.7	19.2	23.5	7.7	7.0	8.4	-17.3	-27.9
- Asia Pacific ¹⁾	24.9	24.3	2.8	-3.8	12.1	11.1	9.7	6.6	12.8	13.2	-2.9	-12.5
- Other Markets ¹⁾	5.4	5.0	9.2	8.7	4.0	2.6	50.7	50.7	1.4	2.3	-37.8	-38.8
EBITA	10.0	12.2	-18.2		12.6	9.1	38.3		-2.6	3.1	-181.7	
EBITA margin	6.9%	8.2%			13.0%	10.0%			-5.2%	5.4%		
Net profit ²⁾	1.8	4.2	-57.1									
Earnings per share in € ²⁾	0.11	0.25	-57.1									

¹⁾ Acc. to customers' location

²⁾ Excluding the non-cash items of amortization and interest for share price warrants

Earnings Development

The Biotechnology Division boosted its first-quarter operating earnings (EBITA = earnings before interest, taxes and amortization) from €9.1 million a year ago to €12.6 million in 2009. As a result, its EBITA margin significantly improved from 10.0% to 13.0%. In addition to sales revenue growth and the economies of scale resulting from this gain as well as favorable exchange rates, the enhanced product mix favoring single-use products especially contributed to the division's significant boost in profitability.

In the Mechatronics Division, the decline in sales revenue resulted in negative operating earnings of €2.6 million (previous year: +€3.1 million).

Accordingly, operating earnings at the Group level were at €10.0 million. EBITA of the comparable year-earlier period was €12.2 million. The EBITA margin fell from 8.2% a year ago to 6.9%.

Excluding the two non-cash items of amortization of €1.8 million (previous year: €1.6 million) and interest of €0.3 million (previous year: €0.5 million) for share price warrants granted to Sartorius Stedim Biotech shareholders, consolidated net profit after minority interest totaled €1.8 million (previous year: €4.2 million). The corresponding earnings per share were at €0.11 (previous year: €0.25).

Balance Sheet and Cash Flow

In the first three months of the current fiscal year, we reduced gross debt from €239.5 million to €232.7 million. Thus, the balance sheet total of the Sartorius Group decreased from €865.0 million as of the reporting date on December 31, 2008, to €858.4 million as of March 31, 2009.

As of March 31, 2009, equity was at €337.1 million (December 31, 2008: €333.4 million). Accordingly, the equity ratio rose from 38.5% as of December 31, 2008, to 39.3% as of March 31, 2009, and continues to remain at a comfortable level. We disclosed the share price warrants, which were granted to shareholders of Sartorius Stedim Biotech S.A. as part of the Stedim transaction, at a fair value of €19.60 per warrant in the current liabilities. This put the total value of all such warrants at €104.8 million (December 31, 2008: €103.7 million).

Net debt for the quarter ended March 31 amounts to €216.2 million (December 31, 2008: €217.6 million). The ratio of net debt to EBITDA is 2.8 compared with 2.7 as of December 31, 2008. The interest coverage ratio – EBITDA divided by cash interest expense of the past 12 months – is 5.4 (December 31, 2008: 5.5). Gearing, which is calculated as the ratio of net debt to equity, is 0.6 (December 31, 2008: 0.7). These key balance sheet ratios for assessing the company's financial condition are thus robust.

Net cash flow from operating activities at €14.9 million (previous year: –€3.0 million) was significantly positive. The increase over the year-earlier figure can be essentially explained by the low cash outflow for building up inventories and a positive quarter-on-quarter effect of income tax. At €7.8 million, cash outflows from investing activities were insignificantly higher than in the year-earlier period (€7.5 million). Due to repayment of bank liabilities, cash flows from financing activities amounted to –€12.3 million (previous year: €6.9 million).

Assessment of Business Development

We are satisfied with the dynamic development of business for the Biotechnology Division. In particular, we take the clearly double-digit growth rates in our business with single-use products and the overproportionate increase in earnings as confirmation of our business model and our excellent prospects for the future.

Amid the unsatisfactory development of earnings and the market conditions that are expected to remain exceptionally difficult for the Mechatronics Division, the Executive Board announced a package of measures for cutting costs on a lasting basis as outlined in the following.



Cost-Cutting Program

The package of cost-cutting measures entails a sustainable reduction of the cost base in the Mechatronics Division by more than €25 million. The goal of these measures is to steer the Mechatronics Division clearly back into the profit zone in 2010. Cutting staff costs by more than €20 million will play a central role in this effort. The company plans to intensively use various options, such as partial retirement, variable pay for employees under the collective bargaining agreement and termination of fixed-term employment contracts. Still, it is expected that the planned cost cuts will not be achievable without layoffs. The corresponding negotiations will begin shortly with employee representative committees. One-time expenses that will be incurred in this context are estimated at about €10–15 million.

Outlook

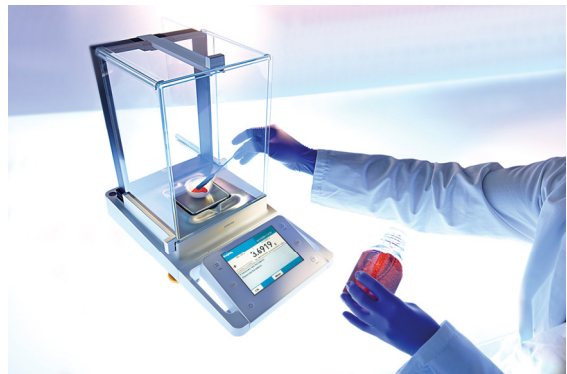
In view of the positive development of first-quarter business and the number of promising projects on which we are currently working together with our customers, we project that sales revenue will grow and earnings will increase overproportionately for the Biotechnology Division in fiscal 2009.

Amid the exceptionally tough market conditions, we anticipate a double-digit decline in sales revenue and a negative earnings contribution for the Mechatronics Division for the current year. By implementing the package of measures planned, we are striving to report positive earnings for this division in 2010.

Given the continued high uncertainty regarding the global economy, Sartorius does not consider it possible to make a precise quantitative forecast for 2009.

Dividends for 2008

The Annual Shareholders' Meeting resolved on April 23, 2009, to pay dividends of €0.42 per preference share and €0.40 per ordinary share to its shareholders for fiscal 2008. These dividends were disbursed on April 24, 2009. The total profit distributed amounted to €7.0 million (previous year: €11.4 million), which equates to a dividend payout ratio of 38.4% (previous year: 37.8%) relative to consolidated net profit excluding non-cash amortization and interest on share price warrants.



Balance Sheet

Assets	March 2009		December 2008	
	€ in mn	%	€ in mn	%
Non-current Assets				
Goodwill	293.2	34.2	292.5	33.8
Intangible assets	121.5	14.2	122.4	14.2
Property, plant and equipment	157.1	18.3	155.7	18.0
Financial assets	5.4	0.6	5.1	0.6
	577.3	67.3	575.8	66.6
Non-current trade and other receivables	3.4	0.4	3.5	0.4
Deferred tax assets	16.8	2.0	14.1	1.6
	597.4	69.6	593.3	68.6
Current Assets				
Inventories	97.4	11.3	95.6	11.1
Trade receivables	124.0	14.4	125.1	14.5
Current tax assets	6.2	0.7	13.5	1.6
Other assets	16.8	2.0	15.5	1.8
Cash and cash equivalents	16.5	1.9	21.9	2.5
	260.9	30.4	271.7	31.4
Total Assets	858.4	100.0	865.0	100.0
Equity and Liabilities				
Equity and Liabilities	March 2009		December 2008	
	€ in mn	%	€ in mn	%
Equity				
Issued capital	17.0	2.0	17.0	2.0
Capital reserves	87.0	10.1	87.0	10.1
Retained earnings (including net profit)	187.7	21.9	185.7	21.5
Minority interest	45.4	5.3	43.7	5.0
	337.1	39.3	333.4	38.5
Non-current Liabilities				
Pension provisions	35.6	4.1	35.1	4.1
Deferred tax liabilities	40.2	4.7	40.0	4.6
Other provisions	8.6	1.0	8.5	1.0
Loans and borrowings	181.9	21.2	192.7	22.3
Other liabilities	0.2	0.0	0.3	0.0
	266.4	31.0	276.6	32.0
Current Liabilities				
Provisions	7.7	0.9	8.9	1.0
Trade payables	39.5	4.6	44.4	5.1
Loans and borrowings	50.9	5.9	46.8	5.4
Current tax liabilities	9.5	1.1	9.8	1.1
Other liabilities	147.3	17.2	145.1	16.8
	254.8	29.7	255.0	29.5
Total Equity and Liabilities	858.4	100.0	865.0	100.0

Income Statement

	1st quarter 2009		1st quarter 2008	
	€ in mn	%	€ in mn	%
Sales revenue	146.0	100.0	149.2	100.0
Cost of sales	-80.0	-54.8	-79.1	-53.0
Gross profit on sales	66.0	45.2	70.1	47.0
Selling and distribution costs	-36.0	-24.7	-37.4	-25.1
Research and development costs	-10.0	-6.9	-10.3	-6.9
General administrative expenses	-10.3	-7.1	-10.6	-7.1
Other operating income and expenses	0.4	0.3	0.5	0.3
	-56.0	-38.4	-57.8	-38.8
Earnings before interest, taxes and amortization (EBITA)	10.0	6.9	12.2	8.2
Amortization ¹⁾	-1.8	-1.2	-1.6	-1.0
Earnings before interest and taxes (EBIT)	8.2	5.6	10.7	7.2
Interest and similar income	0.1	0.1	0.5	0.3
Interest and similar expenses	-4.3	-3.0	-4.1	-2.8
Financial result	-4.2	-2.9	-3.7	-2.5
Profit before tax	4.0	2.7	7.0	4.7
Income tax expense	-3.2	-2.2	-2.3	-1.6
Deferred tax income expenses	2.1	1.4	0.3	0.2
Other taxes	-0.6	-0.4	-0.7	-0.5
	-1.7	-1.2	-2.7	-1.8
Net profit	2.3	1.6	4.3	2.9
Attributable to:				
Equity holders of the parent	0.8	0.5	3.2	2.1
Minority interest	1.5	1.0	1.1	0.7
Earnings per ordinary share (€)	0.05		0.19	
Earnings per preference share (€)	0.05		0.19	

¹⁾ Amortization refers to goodwill and to the purchase price allocation (PPA) to intangible assets according to IFRS 3.

Cash Flow Statement

	1st quarter 2009	1st quarter 2008
	€ in mn	€ in mn
Cash flows from operating activities		
Net profit	0.8	3.2
Minority interest	1.5	1.1
Tax expenses	1.7	2.7
Financial expenses	4.2	3.7
Depreciation amortization of fixed assets	7.8	7.2
Increase decrease in provisions	-0.9	-1.2
Increase decrease in receivables and other assets	-0.4	2.4
Increase decrease in inventories	-0.6	-12.8
Increase decrease in liabilities (without loans and borrowings)	-3.0	1.0
Income taxes received paid	3.8	-10.3
Net cash flow from operating activities	14.9	-3.0
Cash flows from investing activities		
Payments for financial assets	-0.3	-0.6
Payments for property, plant and equipment	-5.5	-4.8
Income from the disposal of fixed assets	0.5	0.2
Payments for intangible assets	-2.5	-2.4
Net cash flow from investing activities	-7.8	-7.5
Cash flows from financing activities		
Changes in capital	0.0	0.0
Interest received	0.1	0.5
Interest paid and other financial charges	-5.3	-3.8
Payments for derivative financial instruments	-0.4	-0.1
Dividends paid to:		
Shareholders of the parent company	0.0	0.0
Minority shareholders	0.0	0.0
Changes in minority interest	0.2	0.0
Loans and borrowings repaid raised	-6.9	10.4
Net cash flow from financing activities	-12.3	6.9
Net increase decrease in cash and cash equivalents	-5.2	-3.7
Cash and cash equivalents at the beginning of the period	21.9	17.7
Net effect of currency translation on cash and cash equivalents	-0.3	-0.2
Cash and cash equivalents at the end of the period	16.5	13.8
Gross debt owed to banks	232.7	217.6
Net debt owed to banks	216.2	203.7

Segment Reports

Segment Report by Division

€ in millions	Biotechnology			Mechatronics			Group		
	1st quarter 2009	1st quarter 2008	Change	1st quarter 2009	1st quarter 2008	Change	1st quarter 2009	1st quarter 2008	Change
Order intake	104.4	97.6	7%	52.6	64.9	-19%	157.0	162.5	-3%
Sales revenue	96.6	90.6	7%	49.5	58.6	-16%	146.0	149.2	-2%
- as a total %	66.1%	60.7%		33.9%	39.3%		100.0%	100.0%	
EBITDA	16.3	12.6	30%	-0.2	5.3	-105%	16.0	17.8	-10%
- as a % of sales revenue	16.9%	13.9%		-0.5%	9.0%		11.0%	12.0%	
Depreciation and amortization	5.5	5.0	9%	2.3	2.1	9%	7.8	7.2	9%
EBITA	12.6	9.1	38%	-2.6	3.1	-182%	10.0	12.2	-18%
- as a % of sales revenue	13.0%	10.0%		-5.2%	5.4%		6.9%	8.2%	
EBIT	10.8	7.5	44%	-2.6	3.1	-182%	8.2	10.7	-23%
- as a % of sales revenue	11.2%	8.3%		-5.2%	5.4%		5.6%	7.2%	
R&D costs	5.9	6.1	-3%	4.1	4.2	-3%	10.0	10.3	-3%
No. of employees at March 31	2,382	2,297	4%	2,253	2,249	0%	4,635	4,546	2%

Segment Report by Region

€ in millions	Europe			North America		
	1st quarter 2009	1st quarter 2008	Change	1st quarter 2009	1st quarter 2008	Change
Sales revenue						
- acc. to customers' location	85.0	92.3	-8%	30.7	27.6	11%
- as a total %	58.2%	61.9%		21.0%	18.5%	
- acc. to company location	97.9	105.2	-7%	30.8	28.0	10%
EBITA	6.6	8.9	-26%	1.8	0.9	93%
- as a % of sales revenue	6.7%	8.4%		5.8%	3.3%	
No. of employees at March 31	3,149	3,030	4%	518	573	-10%

€ in millions	Asia Pacific			Other Markets			Group		
	1st quarter 2009	1st quarter 2008	Change	1st quarter 2009	1st quarter 2008	Change	1st quarter 2009	1st quarter 2008	Change
Sales revenue									
- acc. to customers' location	24.9	24.3	3%	5.4	5.0	9%	146.0	149.2	-2%
- as a total %	17.1%	16.3%		3.7%	3.3%		100.0%	100.0%	
- acc. to company location	17.3	15.9	9%	0.0	0.0		146.0	149.2	-2%
EBITA	1.5	2.3	-36%	0.1	0.1	89%	10.0	12.2	-18%
- as a % of sales revenue	8.7%	14.7%					6.9%	8.2%	
No. of employees at March 31	890	849	5%	78	94	-17%	4,635	4,546	2%

Statement of Changes in Equity

€ in millions									
	Issued capital	Capital reserves	Hedging reserves	Pension reserves	Earnings reserves and retained profits	Difference resulting from currency translation	Total	Minority interest	Total equity
Balance at Jan. 1, 2008	17.0	87.0	2.3	-1.7	201.2	-14.5	291.3	42.8	334.1
Cash flow hedges	0.0	0.0	2.2	0.0	0.0	0.0	2.2	0.5	2.8
Actuarial gains losses from pension provisions	0.0	0.0	0.0	-0.2	0.0	0.0	-0.2	0.2	0.0
Currency translation differences	0.0	0.0	0.0	0.0	0.0	-3.2	-3.2	-1.2	-4.4
Net investment in a foreign operation	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0
Tax effects due to net income recognized directly in equity	0.0	0.0	-0.7	0.0	0.0	0.0	-0.6	-0.2	-0.8
Net income recognized directly in equity	0.0	0.0	1.6	-0.2	0.0	-3.2	-1.8	-0.6	-2.5
Net profit for the period	0.0	0.0	0.0	0.0	3.2	0.0	3.2	1.1	4.3
Total recognized income and expense	0.0	0.0	1.6	-0.2	3.2	-3.2	1.3	0.5	1.8
Dividends	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0
Acquisition of additional shares in subsidiaries	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0
Change in minority interest	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0
Other changes in equity	0.0	0.0	0.0	0.0	0.1	0.0	0.1	-0.1	0.0
Balance at March 31, 2008	17.0	87.0	3.8	-1.9	204.5	-17.8	292.7	43.2	335.9

€ in millions									
	Issued capital	Capital reserves	Hedging reserves	Pension reserves	Earnings reserves and retained profits	Difference resulting from currency translation	Total	Minority interest	Total equity
Balance at Jan. 1, 2009	17.0	87.0	-0.6	-1.3	200.2	-12.6	289.7	43.7	333.4
Cash flow hedges	0.0	0.0	-0.8	0.0	0.0	0.0	-0.8	-0.2	-1.0
Actuarial gains losses from pension provisions	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0
Currency translation differences	0.0	0.0	0.0	0.0	0.0	2.0	2.0	0.4	2.3
Net investment in a foreign operation	0.0	0.0	0.0	0.0	-0.1	0.0	-0.1	0.0	-0.1
Tax effects due to net income recognized directly in equity	0.0	0.0	0.2	0.0	0.0	0.0	0.3	0.1	0.3
Net income recognized directly in equity	0.0	0.0	-0.6	0.0	-0.1	2.0	1.3	0.2	1.6
Net profit for the period	0.0	0.0	0.0	0.0	0.8	0.0	0.8	1.5	2.3
Total recognized income and expense	0.0	0.0	-0.6	0.0	0.7	2.0	2.1	1.7	3.9
Dividends	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0
Acquisition of additional shares in subsidiaries	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0
Change in minority interest	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0
Other changes in equity	0.0	0.0	0.0	0.0	-0.1	0.0	-0.1	0.0	-0.1
Balance at March 31, 2009	17.0	87.0	-1.2	-1.3	200.8	-10.6	291.8	45.4	337.1

Key Figures

€ in millions (unless otherwise specified)	1st quarter 2009	1st quarter 2008	Change in %
Results			
Order intake	157.0	162.5	-3.4
Sales revenue	146.0	149.2	-2.1
EBITDA	16.0	17.8	-10.1
EBITA	10.0	12.2	-18.2
EBIT	8.2	10.7	-22.7
Profit before tax	4.0	7.0	-42.6
Net profit after minority interest	0.8	3.2	-75.1
Earnings per share (in €)	0.05	0.19	-75.1
As a % of sales revenue			
EBITDA	11.0	12.0	
EBITA	6.9	8.2	
EBIT	5.6	7.2	
Profit before tax	2.7	4.7	
Net profit after minority interest	0.5	2.1	
Finances			
Capital expenditures	8.0	7.1	12.4
As a % of sales revenue	5.5	4.8	
Depreciation and amortization	7.8	7.2	8.8
Net cash flow from operating activities	14.9	-3.0	
Net debt	216.2	203.7	6.1
Employees as of March 31	4,635	4,546	2.0
R&D			
R&D costs	10.0	10.3	-3.1
As a % of sales revenue	6.9	6.9	

Notes to the Interim Financial Statements

Recognition and Measurement Principles

The annual consolidated financial statements of Sartorius AG for the period ended December 31, 2008, were prepared in accordance with the accounting standards of the International Accounting Standards Board (IASB) – the International Financial Reporting Standards (IFRS). In the present interim financial statements that were prepared in conformance with the requirements of IAS 34 "Interim financial reporting," basically the same accounting and measurement principles were applied on which the past consolidated financial statements of fiscal 2008 were based.

Furthermore, all interpretations of the International Financial Reporting Interpretations Committee (IFRIC) to be applied effective March 31, 2009, were observed. An explanation of the individual accounting and measurement principles applied is given in the Notes to the Consolidated Financial Statements for the year ended December 31, 2008.

Other Disclosures

This first-quarter report was not reviewed by the group auditor.

Contacts

Sartorius AG
Weender Landstrasse 94–108
37075 Goettingen, Germany

Phone: +49.551.308 0
Fax: +49.551.308 3289
www.sartorius.com

Investor Relations

Jörg Pfirrmann
Senior Vice President Finance
Phone +49.551.308 4076
joerg.pfirrmann@sartorius.com

Andreas Wiederhold
Phone +49.551.308 1668
andreas.wiederhold@sartorius.com

Henriette Meyer
Phone +49.551.308 3232
henriette.meyer@sartorius.com

Group Corporate Communications

Petra Kirchhoff
Vice President
Phone +49.551.308 1686
petra.kirchhoff@sartorius.com

Dominic Grone
Phone +49.551.308 3324
dominic.grone@sartorius.com

This is a translation of the original German-language first-quarter financial report "Sartorius Konzern 3-Monatsbericht." Sartorius AG shall not assume any liability for the correctness of this translation. The original German interim report is the legally binding version. Furthermore, Sartorius AG reserves the right not to be responsible for the topicality, correctness, completeness or quality of the information provided. Liability claims regarding damage caused by the use of any information provided, including any kind of information which is incomplete or incorrect, will therefore be rejected.

Forward-looking Statements Contain Risks

This interim report contains various statements concerning the Sartorius Group's future performance. These statements are based on assumptions and estimates. Although we are convinced that these forward-looking statements are realistic, we cannot guarantee that they will actually apply. This is because our assumptions harbor risks and uncertainties that could lead to actual results diverging substantially from the expected ones. It is not planned to update our forward-looking statements.

Throughout the entire report, differences may be apparent as a result of rounding during addition.

Financial Schedule

May 14, 2009

Investors' day "Biotech Suppliers Finance Day" at the ACHEMA 2009 in Frankfurt am Main, Germany

July 2009*

Publication of the first-half figures for 2009

October 2009*

Publication of the nine-month figures for 2009

November 9–11, 2009

Company presentation at the German Equity Forum, Fall 2009, in Frankfurt am Main, Germany

* Tentative date scheduled