



Sartorius AG
2008 Financial Statements



Sartorius AG
2008 Financial
Statements

Forward-looking Statements Contain Risks

This annual report contains 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.

This is a translation of the original German-language financial statements. Sartorius shall not assume any liability for the correctness of this translation. The original German financial statements are the legally binding version. Furthermore, Sartorius 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.

Throughout these financial statements, differences may be apparent as a result of rounding during addition.

Contents

Financial Statements and Notes

- 4 Balance Sheet
- 5 Income Statement
- 6 Notes to the Financial Statements

Supplementary Information

- 26 Executive Board and
Supervisory Board | Positions Held
- 28 About This Publication

As of fiscal 2001, Sartorius AG's financial statements and the consolidated financial statements are no longer described in combination. Because Sartorius AG prepared the consolidated financial statements according to the "International Financial Reporting Standards" (IFRS), the Group financial statements are explained separately.

Balance Sheet

Assets	Notes	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
A. Fixed assets			
I. Intangible assets	(4)	989	1,248
II. Property, plant and equipment	(5)	17,956	17,179
III. Financial assets	(6)	368,158	311,320
		387,103	329,747
B. Current assets			
I. Inventories	(7)	17,249	16,618
II. Trade and other receivables	(8)	40,677	37,371
III. Securities	(9)	15,069	16,082
IV. Cash on hand, deposits in banks		2,044	4,887
		75,039	74,958
C. Prepaid expenses	(10)	1,429	44
		463,571	404,749
Equity and Liabilities			
	Notes	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
A. Equity			
I. Issued capital	(11)	18,720	18,720
II. Capital reserves		101,397	101,397
III. Earnings reserves	(12)	25,276	25,276
IV. Retained profit incl. net profit	(12)	27,810	31,610
		173,203	177,003
B. Provisions	(13)	128,328	77,141
C. Liabilities	(14)	162,040	150,605
		463,571	404,749

Income Statement

	Notes	2008 € in K	€ in K	2007 € in K
1. Sales revenue	(18)	129,687		175,888
2. Increase in inventories of finished goods and work in progress	(19)	447		715
			130,134	176,603
3. Work performed by the enterprise and capitalized	(20)	113		169
4. Other operating income	(21)	14,535		38,914
			14,648	39,083
			144,782	215,686
5. Raw materials and supplies	(22)	48,404		65,296
6. Employee benefits expense	(23)	50,176		65,706
7. Depreciation and amortization expense	(24)	3,459		6,065
8. Other operating expenses	(25)	41,689		54,245
9. Income from investments	(26)	18,699		5,879
10. Impairment losses on financial assets and current marketable securities	(27)	1,527		0
			126,556	185,433
11. Earnings before interest and taxes			18,226	30,253
12. Interest and similar income	(28)	914		1,441
13. Interest and similar expenses	(28)	11,142		8,248
			-10,228	-6,807
14. Profit before tax			7,998	23,446
15. Income tax expense	(29)	233		1,089
16. Other taxes		143		161
			376	1,250
17. Net profit			7,622	22,196
18. Profit brought forward			20,188	9,414
19. Retained profit			27,810	31,610

Notes to the Financial Statements

Cash Flow Statement	2008 € in K	2007 € in K
Net profit for the period	7,622	22,196
Depreciation and amortization of fixed assets	3,459	6,065
Increase in long-term provisions	554	1,124
Other material non-cash income (-) expenses (+)	-10,779	-5,061
Cash earnings acc. to the DVFA SG*	856	24,324
Decrease (-) Increase in short-term provisions	-2,281	-9,245
Gains resulting from the disposal of fixed assets	-1	-17,526
Increase (-) in inventories	-631	-2,464
Increase (-) in short- and long-term receivables and other assets	-4,692	-7,053
Decrease (-) Increase in liabilities (excluding liabilities to banks)	-3,871	13,463
Cash flows from operating activities	-10,620	1,499
Proceeds from fixed asset disposals	117	18,585
Payments (-) for intangible assets	-388	-793
Payments (-) for property, plant and equipment	-3,598	-7,364
Payments (-) for financial assets	-7,239	-107,157
Cash flows from investing activities	-11,108	-96,729
Dividend payments (-)	-11,421	-10,740
Additions to and repayment (-) of financial liabilities	30,306	110,464
Cash flows from financing activities	18,885	99,724
Change in cash and cash equivalents	-2,843	4,494
Cash and cash equivalents at beginning of period	4,887	393
Cash and cash equivalents at end of period	2,044	4,887
Composition of cash and cash equivalents	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
Cash on hand, deposits in banks	2,044	4,887

The cash flow statement shows the flow of funds and payments according to operating activities, investing activities and financing activities in a table.

Cash flows from operating activities are calculated according to the indirect method, which means that expenses not affecting cash payments are added to the net profit for the year, whereas income not affecting cash payments is deducted.

* DVFA | SG = The Methods Committee of the Society of Investment Professionals in Germany, also commonly referred to as the German association for financial analysis and asset management

Notes to the Individual Balance Sheet Items

1. Accounting and Valuation Methods

The financial statements for Sartorius AG were prepared according to the rules of the German Commercial Code (HGB) and the German Stock Corporation Law (AktG).

With economic effect as of April 1, 2007, the Biotechnology Division of Sartorius AG became an independent entity. As a result of this carve-out, it was incorporated into the subsidiary Sartorius Stedim Biotech GmbH, Goettingen. Furthermore, effective May 1, 2007, various administrative functions were outsourced at book value to the subsidiary Sartorius Corporate Administration GmbH, Goettingen. This restructuring within the Group resulted in significant changes to Sartorius AG's revenue and cost structures. For this reason, the income statement for the reporting year of 2008 is comparable only to a limited extent with that of the previous year.

The only change made from the previous year in the accounting and valuation methods is that impairment losses of capital assets having minor importance were adjusted to the amended tax regulations.

2. Structure of the Balance Sheet and Income Statement

To enhance the clarity of the presentation, individual balance sheet and income statement items have been combined and reported separately in the Notes.

3. Currency Translation

Short-term receivables and liabilities in foreign currencies were translated at the year-end rates on December 31, 2008.

4. Intangible Assets

	Concessions, industrial property rights and similar rights as well as licenses for such rights € in K	Payments on account € in K	Total € in K
Intangible assets developed during the fiscal year as follows:			
Gross book values as of Jan. 1, 2008	3,097	21	3,118
Investments	384	4	388
Disposals	0	0	0
Transfers	8	5	13
Gross book values as of Dec. 31, 2008	3,489	30	3,519
Amortization as of Jan. 1, 2008	1,870	0	1,870
Amortization in 2007	660	0	660
Disposals	0	0	0
Amortization as of Dec. 31, 2008	2,530	0	2,530
Net book values as of Dec. 31, 2008	959	30	989
Net book values as of Dec. 31, 2007	1,227	21	1,248

Intangible assets acquired are stated at acquisition cost less the accumulated, regular amortization. Where diminution in value is anticipated over the long term, these assets are written off as impairment losses.

Intangible assets are subject to scheduled amortization according to the straight-line method over a useful life of two to five years.

5. Property, Plant and Equipment

Property, plant and equipment developed as follows:	Land and leasehold rights and improvements, including buildings on third-party land € in K	Technical equipment and machinery € in K	Other equipment, factory and office equipment € in K	Payments on account relating to plant and equipment and construction in progress € in K	Total € in K
Gross book values as of Jan. 1, 2008	26,053	11,514	23,194	391	61,152
Additions	47	752	2,355	443	3,597
Disposals	0	26	293	0	319
Transfers	0	0	337	-350	-13
Gross book values as of Dec. 31, 2008	26,100	12,240	25,593	484	64,417
Depreciation as of Jan. 1, 2008	14,357	10,002	19,614	0	43,973
Depreciation in 2008	627	577	1,595	0	2,799
Disposals	0	25	286	0	311
Depreciation as of Dec. 31, 2008	14,984	10,554	20,923	0	46,461
Net book values as of Dec. 31, 2008	11,116	1,686	4,670	484	17,956
Net book values as of Dec. 31, 2007	11,696	1,512	3,580	391	17,179

Property, plant and equipment are reported at acquisition cost or production cost and, if subject to depreciation, are depreciated as scheduled. The straight-line method was used for calculating the depreciation reported on Sartorius AG's annual financial statements up to and including 1994. For additions reported under property, plant and equipment, scheduled depreciation has been calculated since fiscal 1995 according to the declining balance method of depreciation, because in this business year, tax write-offs in accordance with Section 3 of the German law on subsidies for former eastern borderline areas ("Zonenrandförderungsgesetz") were eliminated along with this law.

German tax simplification methods are applied. Where diminution in value is anticipated over the long term, unscheduled impairment losses are recorded.

On account of the tax write-offs reported in previous fiscal years, the annual profit disclosed by Sartorius AG is €0.1 million higher than the amount that would have had to have been declared without these write-offs. The tax savings of the previous fiscal years yield a tax burden of approx. €0.1 million.

Amortization | depreciation of fixed assets is based on the following periods of useful life:

Buildings	15 to 50 years
Machinery	5 to 15 years
Factory and office equipment	3 to 13 years

6. Financial Assets

	Shares in affiliated companies € in K	Participating interests € in K	Total € in K
Financial assets changed as follows:			
Gross book values as of Jan. 1, 2008	311,560	477	312,037
Investments	57,446	14	57,460
Disposals	0	108	108
Transfers	0	0	0
Gross book values as of Dec. 31, 2008	369,006	383	369,389
Impairment losses as of Jan. 1, 2008	348	369	717
Impairment losses in 2008	500	14	514
Disposals	0	0	0
Reversals	0	0	0
Impairment losses as of Dec. 31, 2008	848	383	1,231
Net book values as of Dec. 31, 2008	368,158	0	368,158
Net book values as of Dec. 31, 2007	311,212	108	311,320

Financial assets are accounted for at acquisition cost or at par value, unless they have to be reported at a lower recoverable amount on the balance sheet date.

In fiscal 2007, Sartorius AG acquired a controlling stake in Stedim S.A. listed on the French stock exchange and headquartered in Aubagne, France (now known as Sartorius Stedim Biotech S.A.). In the first stage of this acquisition, Sartorius AG purchased all of the shares in VL Finance S.A.S., Aubagne, France, which owned 49.8% of Stedim S.A. shares and had been the company's majority shareholder until then. Sartorius AG then increased its holding to a 69.8% majority stake by transferring its shares in Sartorius Stedim Biotech GmbH, Goettingen, into Stedim S.A. in the form of a capital increase against the issue of new shares. Both of the stages described were completed on June 29, 2007. In fiscal 2008, Sartorius AG acquired an additional 1.5% of the shares in Sartorius Stedim Biotech S.A. Moreover, this gave rise to subsequent acquisition costs in connection with the share price warrants granted to non-affiliated shareholders. As of December 31, 2008, the deal as a whole entailed acquisition costs for the shares in the two companies amounting to €306,583 K. Of this amount, shares in VL Finance S.A. accounted for €151,520 K and shares in Sartorius Stedim Biotech S.A. €155,063 K.

The fair value of the shares in VL Finance S.A.S. derived from the market value of shares in Sartorius Stedim Biotech S.A. was €76 million at December 31, 2008, due to a decrease in the price of the Sartorius Stedim Biotech S.A. shares. No use was made of the mechanism for unscheduled write-downs on the shares, as the fair value of the total Stedim transaction, including the shares directly held in Sartorius Stedim Biotech S.A., did not go below the book value on the whole.

The acquisition of the equity in Sartorius Mechatronics Hungaria Kft., Budapest, Hungary, and the founding of Sartorius Mechatronics Poland Sp. z.o.o., Poznan, Poland, resulted in a further increase in the financial assets.

The impairment loss reported for fiscal 2008 concerns the book value of the participating interest of Sartorius Mechatronics C&D GmbH & Co. KG, Aachen, Germany, because of the continued impairment of shares in this company.

Please refer to "Major Affiliates and Ownership."

7. Inventories

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
Raw materials and supplies	7,177	6,955
Work in progress	5,638	4,860
Finished goods and merchandise	4,186	4,629
Payments on account	248	174
	17,249	16,618

Under the item "Inventories," raw materials and supplies, including merchandise, are reported at average acquisition cost; work in progress and finished goods, at manufacturing or production cost. Besides including direct costs that can be allocated to these materials, the acquisition cost covers the appropriate portion of production and materials handling overhead and normal depreciation | amortization, provided that the latter relates to production. Interest on loans is not capitalized.

The fixed valuation method is used to a certain extent for raw materials and supplies.

Lower stock market or market prices or lower attributable values are taken into account by write-offs. Where inventory risks exist, such as the risk of reduced shelf life as a result of storage periods or limited usability, inventories are marked down accordingly.

8. Trade and Other Receivables

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
Trade receivables	7,797	9,034
Receivables from affiliated companies	23,664	21,491
Receivables from companies in which investments are held	0	0
Other assets	9,216	6,846
	40,677	37,371
Of which due in more than one year:		
Trade receivables	0	0
Other assets	1,199	1,065
	1,199	1,065

Trade and other receivables were reported so that all discernible risks are covered. General credit risks are sufficiently covered by lump-sum valuation allowances. Liabilities arising from insurance for financial backing are recognized in the actuarial reserves according to the company's business plan.

9. Securities

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
Own (treasury) shares	15,069	16,082

As authorized at the Annual Shareholders' Meeting on June 21, 2000, Sartorius AG repurchased its own shares pursuant to Section 71, Subsection 1, No. 8, of the German Stock Corporation Law (AktG).

A reserve for the corresponding amount was set up for own shares according to Section 272, Subsection 4, Sentence 1, of the German Commercial Code (HGB). Other earnings reserves were used to establish this reserve pursuant to Section 272, Subsection 4, Sentence 3, of HGB.

Between October 27, 2000, and the reporting date, 831,944 ordinary shares were repurchased at an average price of €11.27 and 840,983 preference shares were repurchased at an average price of €7.98. This corresponds to €1,673 K or 8.9% of the company's issued capital. No treasury shares were purchased in fiscal 2008.

Due to the drop in the prices of Sartorius shares, treasury shares had to be written down for a total of €1,013 K.

10. Prepaid Expenses

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
Loan discounts	1,168	0
Other prepaid expenses	261	43
	1,429	43

The loan discount is released over the term of the finance loan.

11. Issued Capital

Sartorius AG's capital stock is divided into 9,360,000 bearer-type ordinary shares and 9,360,000 non-voting preference shares, each having a calculated par value of €1.

12. Earnings Reserves and Retained Profits

	€ in K
The reserves for own (treasury) shares of Sartorius AG developed as follows in the reporting year:	
At Jan. 1, 2008	16,082
Withdrawal acc. to Sec. 272, Subsec. 4, of HGB*	1,013
At Dec.31, 2008	<u>15,069</u>

The earnings allocated to the reserves equate to the amount reported as assets for own (treasury) shares. The corresponding amount is reported under the current marketable securities item.

	€ in K
The other earnings reserves developed as follows during the reporting year:	
At Jan. 1, 2008	9,194
Reclassified from the reserves for own shares	1,013
At Dec. 31, 2008	<u>10,207</u>

The retained profit of €27,810 K consists of the profit brought forward of €20,188 K and the net profit of €7,622 K.

* = Handelsgesetzbuch: German Commercial Code

13. Provisions

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
Provisions for retirement benefits and similar obligations	14,721	13,960
Tax provisions	494	475
Deferred tax provisions	250	250
Other provisions	112,863	62,456
	128,328	77,141

Provisions for retirement benefits were calculated according to actuarial principles for the going-concern value stipulated by Section 6a of the German Income Tax Law (EStG).

Besides allowing for Sartorius AG's general in-house retirement plan, these provisions are set aside to meet individually agreed pension plans for active and former Executive Board members and senior corporate officers. Since the termination of the General Pension Plan in 1983, the portion of the general provisions is only for employees who began their employment with the company before

January 1, 1983. A total of €8.8 million of the reserves set up on December 31, 2008, is allocated to Sartorius AG's general in-house retirement plan, and €5.9 million to cover individually agreed pension plans. The valuation of these provisions was based on an assumed 6% rate of interest and according to the reference charts "Richttafeln 2005 G" published by Dr. Klaus Heubeck.

In assessing the amounts for the remaining provisions, all discernible risks and uncertain liabilities have been considered.

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
The other provisions include the following amounts:		
Warrants	103,691	50,777
Employee benefits expense	3,966	6,218
Payments to employees on the early retirement plan for offsetting reduced work hours	2,432	2,638
Commissions, credit notes, bonuses	402	385
Warranties	220	287
Other	2,152	2,151
	112,863	62,456

14. Liabilities

This item consists of the following:	Balance sheet	Remaining term	Remaining term	Balance sheet	Remaining term	Remaining term
	Dec. 31, 2008 € in K	of more than five years € in K	of up to one year € in K	Dec. 31, 2007 € in K	of more than five years € in K	of up to one year € in K
Liabilities to banks	73,557	0	11,640	43,251	0	34,348
Payments received on account of orders	398	0	398	417	0	417
Trade payables	5,348	0	5,348	5,249	0	5,249
Payables to affiliated companies	81,148	0	81,148	100,085	0	100,085
Payables to companies in which investments are held	0	0	0	16	0	16
Other payables	1,589	0	1,589	1,587	0	1,587
	162,040	0	100,123	150,605	0	141,702

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
Other liabilities relate to the following:		
Taxes	801	831
Liabilities relating to social security	0	0

Liabilities are reported as the amounts repayable.

15. Contingent Liabilities

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
Guarantees (suretyships) and other contingent liabilities	35,841	158,125

Guarantees (suretyships) exist in the amount of €35,241 K granted to affiliated companies that are included in the consolidated financial statements.

16. Other Financial Obligations

	Dec. 31, 2008 € in K	Dec. 31, 2007 € in K
In addition to provisions, liabilities and contingent liabilities, other financial obligations consist of the following:		
Rental and leasing contracts		
– due in fiscal 2009 (previous year: due in fiscal 2008)	1,241	1,199
– due in any one fiscal year from 2010 to 2013 (previous year: due in any one fiscal year from 2009 to 2012)	1,302	1,448
Obligation from purchase contracts	0	0
Forward exchange transactions for hedging of commodity trade	0	7,473
	2,543	10,120

There are no other financial obligations to affiliated companies.

17. Derivative Financial Instruments

Within the scope of its operations, Sartorius AG uses derivative financial instruments to hedge against risks arising from changes in interest and foreign exchange rates. Sartorius AG does not hold or issue derivative financial instruments for speculative purposes. Banks of first-rate creditworthiness are the exclusive contracting parties of these financial transactions that Sartorius AG conducts centrally by strictly separating trading, settlement and control within defined limits.

To hedge against future risks arising from changes in the foreign exchange rates, Sartorius AG bought currency options during 2008 up to the reporting date. For existing forward exchange transactions, a reserve for anticipated losses related to incomplete contracts of €153 K had to be set up.

To reduce its exposure to risks of changes in interest rates, Sartorius AG uses interest derivatives in the form of interest caps. On the reporting date of December 31, 2008, Sartorius AG had two interest caps with an interest rate ceiling of 4.5% and three interest caps with a ceiling of 2.5%, each against the 6-month EURIBOR. The fair value, determined according to the marking-to-market method, of the interest cap reported under other receivables was €10 K as of December 31, 2008.

On the reporting date, Sartorius AG concluded forward exchange transactions as follows:

	Nominal value Dec. 31, 2008 € in K	Nominal value Dec. 31, 2007 € in K
Forward exchange transactions	0	7,473
Currency option transactions	55,190	57,752
	55,190	65,225
	Fair value Dec. 31, 2008 € in K	Fair value Dec. 31, 2007 € in K
Forward exchange transactions	0	799
Currency option transactions	–3,640	2,985
	–3,640	3,784

Fair values are determined according to the marking-to-market method.

Notes to the Income Statement

18. Sales Revenue

	2008 € in K	2007 € in K
Sales revenue, which is broken down by business area and geographical market, is as follows:		
Germany	50,843	61,151
All other countries	78,844	114,737
Total	129,687	175,888

19. Increase in Inventories of Finished Goods and Work in Progress

	2008 € in K	2007 € in K
Finished goods	-330	147
Work in progress	777	568
	447	715

20. Work Performed by the Enterprise and Capitalized

	2008 € in K	2007 € in K
This item is apportioned as follows to the items under "Property, plant and equipment":		
Machinery	25	30
Construction in progress	55	75
Factory and office equipment	0	13
Tools and equipment	33	51
	113	169

21. Other Operating Income

	2008 € in K	2007 € in K
Income due to currency translation	5,336	4,378
Income from affiliated companies	4,585	8,953
Income from the release of provisions	1,644	784
Government grants, promotions and other subsidies	993	1,640
Income from the reversal of write-downs	423	2,078
Income from investment disposals	0	17,526
Reversal of write-downs on financial assets	0	1,945
Other income	1,554	1,610
	14,535	38,914

22. Raw Materials and Supplies

	2008 € in K	2007 € in K
Cost of raw materials and supplies of purchased merchandise	46,902	63,669
Cost of purchased services	1,502	1,627
	48,404	65,296

23. Employee Benefits Expense

	2008 € in K	2007 € in K
Wages and salaries	41,040	55,038
Social security	7,275	9,127
Expenses for retirement benefits and pensions	1,861	1,541
	50,176	65,706

Sartorius AG employed an average of 796 people in the reporting year (2007: 975).

24. Depreciation and Amortization Expense

	2008 € in K	2007 € in K
Depreciation and amortization of intangible assets and property, plant and equipment:		
Scheduled	3,459	6,065

25. Other Operating Expenses

	2008 € in K	2007 € in K
Costs for outside services	15,890	19,001
Expenses due to currency translation	6,614	3,805
Commissions and license fees	5,215	7,520
Car and leasing expenses	2,120	2,627
Freight and packaging expenses	2,058	3,238
Advertising	1,608	2,122
Itemized allowance for bad debts and lump-sum valuation adjustment on receivables	1,553	950
Maintenance costs	1,402	1,875
Travel and entertainment expenses	1,315	2,161
Legal and consulting expenses including patent costs	1,270	3,497
Rent and leasing expenses	560	1,491
Office, mail and other administrative expenses	458	1,050
Insurance expenses	166	1,647
Other finance charges	152	193
Other expenses	1,308	3,068
	41,689	54,245

26. Income from Investments

	2008 € in K	2007 € in K
Income from investments	18,699	5,879

The entire amount of Sartorius AG's income from investments consists of dividend payments from subsidiaries.

27. Impairment Losses on Financial Assets and Marketable Securities

	2008 € in K	2007 € in K
Impairment losses on financial assets	514	0
Impairment losses on own (treasury) shares	1,013	0
	1,527	0

Impairment losses on financial assets were recognized particularly for shares of Sartorius Mechatronics C&D GmbH & Co. KG, Aachen, Germany.

28. Interest

	2008 € in K	2007 € in K
Interest and similar income	914	1,441
- of which from affiliated companies	(593)	(551)
Interest and similar expenses	11,142	8,248
- of which to affiliated companies	(3,758)	(1,978)
	-10,228	-6,807

Interest was impacted in particular by the loans taken out in conjunction with the acquisition of Stedim S.A. and the associated refinancing measures.

29. Income Tax Expense

	2008 € in K	2007 € in K
Current income taxes	228	249
Deferred taxes	0	30
Income taxes of the previous years	5	810
	233	1,089

Sartorius AG's income tax rate for the year ended was 3.0% (2007: 4.7%).

Real property tax and motor vehicle tax are essentially reported under the "Other taxes" item.

30. Other Information

Declaration acc. to Sec. 285, No. 16, of the German Commercial Code (HGB)

The declaration prescribed by Section 161 of the German Stock Corporation Law (AktG) was submitted on December 11, 2008, and made available to the shareholders on the company's website at www.sartorius.com.

Disclosure acc. to Sec. 285, No. 17, of HGB

The consultation expenses reported under other operating expenses include the fees for services rendered by the auditing company for the following:

	2008 € in K	2007 € in K
Auditing of the first-half financial report, the annual financial statements of the parent corporation and of the consolidated financial statements	190	122
Tax consultation incl. preparation of tax returns	52	83
Other services	93	623
	335	828

Total Remuneration of the Supervisory Board and the Executive Board

	2008 € in K	2007 € in K
Remuneration of the Supervisory Board Members		
– Total remuneration	384	459
– Fixed remuneration	150	150
– Performance-related remuneration	106	181
– Compensation for committee work	30	30
– Meeting attendance fee	98	98
– Compensation for individually performed services	0	216*

* lawyers' fees for Hengeler & Müller, whose partner Prof. Gerd Krieger is a member of the Supervisory Board

	2008 € in K	2007 € in K
Remuneration of the Executive Board Members		
Total		
– Total remuneration	1,362	1,665
– Performance-independent remuneration	806	875
– Performance-related remuneration	556	790
– Phantom stock units paid out	0	0
– Fair value of the phantom stock units held (see separate table)	205	513
– Fair value of mid-term remuneration components	63	63
– Addition to provisions for employee benefits	82	60

Dr. Joachim Kreuzburg

– Total remuneration	789	915
– Performance-independent remuneration	480	440
– Performance-related remuneration	309	475
– Phantom stock units paid out	0	0
– Fair value of the phantom stock units held (see separate table)	130	319
– Fair value of mid-term remuneration components	63	63
– Addition to provisions for employee benefits	42	15

Dr. Günther Maaz

– Total remuneration	573	498
– Performance-independent remuneration	326	285
– Performance-related remuneration	247	213
– Phantom stock units paid out	0	0
– Fair value of the phantom stock units held (see separate table)	75	194
– Addition to provisions for employee benefits	40	24

Olaf Grothey until Sept. 7, 2007

– Total remuneration*		252
– Performance-independent remuneration		150
– Performance-related remuneration		102
– Phantom stock units paid out		**
– Fair value of the phantom stock units held (see separate table)		0
– Addition to provisions for employee benefits		21

* up to the time the Board member left the company

** liquidated within the scope of an overall agreement at the time he left the company

	Number of phantom stock units	Fair value when granted on Jan. 1 of the particular year € in K	Fair value at year-end on Dec. 31, 2008 € in K	Paid out € in K
Components with a Long-term Incentive Effect				
Dr. Joachim Kreuzburg				
- Tranche of phantom stock units for 2005	4,053	63	33	0
- Tranche of phantom stock units for 2006	3,871	83	32	0
- Tranche of phantom stock units for 2007	3,593	125	28	0
- Tranche of phantom stock units for 2008	4,754	138	37	0
	16,271	409	130	
Dr. Günther Maaz				
- Tranche of phantom stock units for 2005	2,594	40	21	0
- Tranche of phantom stock units for 2006	2,580	55	21	0
- Tranche of phantom stock units for 2007	1,796	63	14	0
- Tranche of phantom stock units for 2008	2,507	73	19	0
	9,477	231	75	

A phantom stock plan established in fiscal 2005 is used as the variable component that serves as a long-term incentive and is subject to risk. This remuneration component depends on the development of the Sartorius share price over a period of at least three years and is payable only if this price exceeds an established appreciation or the development of a comparative index.

By having introduced this component, Sartorius AG followed the suggestion of the German Corporate Governance Code. The specific nature of this component is explained in the Remuneration Report printed in the Group Management Report. Please refer to this information in the 2008 Annual Report for the Sartorius Group.

	2008 € in K	2007 € in K
Remuneration of former managing directors		
Remuneration of former managing directors, Executive Board members and their surviving dependents	256	1,896
Pension obligations to former managing directors, members of the Executive Board and their surviving dependents	4,683	4,377

Non-periodic Income and Expenses

Non-periodic income and expenses are items that do influence current results, but concern changes in transactions of the past years. Essentially, these are reported in other operating income and in employee benefits expense.

For Sartorius AG, the income to be allocated to the other business years is €2,229 K (2007: €5,071 K). This amount includes income from the reversal of valuation allowances and the release of provisions. Sartorius AG's non-periodic expenses are €458 K (2007: €1,764 K) and concern severance pay, in particular.

Number of Employees

	2008	2007
Average workforce employed during the fiscal year:		
Wage earners	0	402
Salaried employees	796	573
	796	975

Since the introduction of the reformed collective wage agreement (ERA = Entgelt-Rahmen-Abkommen) at Sartorius AG in Germany, the

company no longer differentiates between wage earners and salaried employees and reports the number as one item.

Proposal for Appropriation of Profits

The Executive Board will submit a proposal to the Annual Shareholders' Meeting to appropriate the retained profit of €27,809,559.63 reported

by Sartorius AG for the year ended December 31, 2008, as follows:

	€
Payment of a dividend of €0.40 per ordinary share	3,411,222.40
Payment of a dividend of €0.42 per preference share	3,577,987.14
Unappropriated profit carried forward	20,820,350.09
	27,809,559.63

Disclosure acc. to Section 160, Subsection 1, No. 4, of the German Stock Corporation Law (AktG)

Subject to approval by the Supervisory Board, the Executive Board is authorized to sell treasury shares held by the corporation, including sales through channels other than the stock exchange or by tendering an offer to all shareholders in proportion to

their participation in the company, provided that these shares are offered within the scope of acquiring companies or shareholdings in companies in return. Under these circumstances, the pre-emptive rights of the shareholders are excluded.

Disclosure acc. to Section 160, Subsection 1, No. 8, of the German Stock Corporation Law (AktG)

The following information was available acc. to Section 21, Subsection 1 or 1a, of the German Securities Trading Act (WpHG) as of December 31, 2008:

Name Company	Residence Registered office	Threshold value (THV)		Percentage of own voting shares held (55.59% were additionally held by the community of heirs to Horst Sartorius's estate)	Percentage of voting shares held	Information on share allocation according to §22 of the German Securities Trading Act (WpHG)
		Which THV was attained, exceeded or which % went below THV	Day on which THV was attained, exceeded or on which % went below THV			
Mrs. C. Franken	Bovenden, Germany	50%	Apr. 1, 2002	3.66%	59.25%	None
Mrs. U. Baro	Munich, Germany	50%	Apr. 1, 2002	4.99%	60.58%	None
Mrs. K. Sartorius-Herbst	Northeim, Germany	50%	Apr. 1, 2002	7.54%	63.13%	None
Prof. Dr. A. Picot (Executor of Horst Sartorius's estate)	Gauting, Germany	50%	Apr. 1, 2002		55.59%	All acc. to Subsec. 1, Sent. 1, No. 6
Bio-Rad Laboratories Inc.	Hercules, California, USA	25%	Sept. 19, 2006		25.02%	All acc. to Subsec. 1, No. 1

Goettingen, February 23, 2009

Sartorius Aktiengesellschaft
The Executive Board

Declaration of the Executive Board

To the best of our knowledge, we declare that the financial statements for fiscal 2008 present a true and fair view of the actual net worth, financial situation and profitability of the company in accordance with the accounting standards used in preparing these statements. We also certify that the progress of the company's business, including its business performance and situation, are represented accurately in the Group Management Report in all material respects and present the most important opportunities and risks of the company's future development during the fiscal year.

Goettingen, February 2009

Sartorius Aktiengesellschaft
Executive Board



Dr. Joachim Kreuzburg



Dr. Günther Maaz

Independent Auditors' Report

We audited the year-end financial statements – consisting of the balance sheet, income statement and notes to the financial statements – including the accounting records of Sartorius Aktiengesellschaft (AG), Goettingen, Germany, and the management report summarized in the Group Management Report for the business year from January 1 through December 31, 2008. These statements and reports drawn up according to German commercial accounting standards are the responsibility of the Executive Board of Sartorius AG. Our responsibility is to express an opinion on Sartorius AG's financial statements, including its accounting records, and management report, based on our audit.

We conducted our audit in accordance with Section 317 of the German Commercial Code (HGB), taking into account the principles of proper auditing established by the German Institute of Independent Auditors, "Institut der Wirtschaftsprüfer." These principles require that we plan and perform the audit to obtain reasonable assurance that there are no misrepresentations and infractions that would have a material impact on the presentation of the net worth, financial position and earnings in the financial statements, in consideration of the accounting principles to be applied, or in the management report. In determining the audit focus, information on the business activities and the economic and legal background of the company as well as expectations concerning possible errors are considered. Within the scope of the audit, the effectiveness of the internal controlling system and evidence supporting the amounts and disclosures of the accounting records, financial statements and management report are predominantly examined on a test basis. This audit covers assessment of the accounting principles applied and the significant estimates made by the Executive

Board as well as the overall presentation of the year-end financial statements and the management report. We believe that our audit provides a reasonable basis for our opinion.

Our audit did not result in any objections.

According to our assessment based on the information we obtained during the audit, the financial statements drawn up by Sartorius Aktiengesellschaft, Goettingen, Germany, present fairly, in all material respects, the net worth, financial position and earnings of the joint stock corporation in conformity with generally accepted accounting principles. The management report is consistent with the financial statements and provides an overall true and fair view of the company's situation, and accurately presents the risks of its future development.

Hanover, February 23, 2009

Deloitte & Touche GmbH
Wirtschaftsprüfungsgesellschaft
(Name of the independent auditing company)



Jürgen Reker
Auditor



Dieter Tenambergen
Auditor

Major Affiliates and Ownership

	Ownership in %	Equity at Dec. 31, 2008 € in K	Net profit at Dec. 31, 2008 € in K
Europe			
VL Finance S.A.S., Aubagne, France	100.0	56,735	2,225
Sartorius Stedim Biotech S.A., Aubagne, France along with its subsidiaries:	71.1	74,750	5,655
Sartorius Stedim Aseptics S.A., Lourdes, France	100.0	2,824	1,416
Integrated Biosystems Sàrl, Fribourg, Switzerland	100.0	1,753	374
Sartorius Stedim SUS S.A.R.L., M'Hamdia, Tunisia	100.0	207	254
Sartorius Stedim Biotech GmbH, Goettingen, Germany, along with its subsidiaries:	100.0	65,087	11,333
Sartorius Stedim UK Ltd., Epsom, U.K.	100.0	1,491	1,365
Wave Biotech AG, Tagelswangen, Switzerland	80.0	1,464	91
Sartorius Stedim France S.A.S., Aubagne, France	100.0	2,320	1,034
Sartorius Stedim Italy S.p.A., Florence, Italy	100.0	2,994	588
Sartorius Stedim Spain S.A., Madrid, Spain	100.0	-762	225
Sartorius Stedim Poland sp. z.o.o., Poznan, Poland	100.0	No business activity in 2008	
Sartorius Stedim Hungaria Kft, Budapest, Hungary	100.0	92	90
Sartorius Stedim Netherlands B.V., Nieuwegein, Netherlands	100.0	13	-25
Sartorius Stedim Belgium N.V., Vilvoorde, Belgium	100.0	267	197
Sartorius Stedim Austria GmbH, Vienna, Austria	100.0	784	430
Sartorius Stedim Switzerland GmbH, Dietikon, Switzerland	100.0	65	8
Sartorius Stedim Japan K.K., Tokyo, Japan	100.0	1,005	475
Sartorius Stedim Singapore Pte. Ltd., Singapore	100.0	450	141
Munktell & Filtrak GmbH, Baerenstein, Germany*)	49.0	4,097	945
Distribo GmbH, Goettingen, Germany*)	26.0	782	172
Sartorius Stedim Filters Inc., Yauco, Puerto Rico	100.0	3,024	2,865
Sartorius Stedim Plastics GmbH, Goettingen, Germany	100.0	1,788	757
Sartorius Stedim F&B GmbH, Goettingen, Germany	100.0	4,588	684
Sartorius Technologies & Services GmbH, Goettingen, Germany	100.0	101	23
Sartorius Stedim Lab Ltd., Louth, U.K.	100.0	1,285	785
Sartorius Stedim Malaysia Sdn. Bhd., Kuala Lumpur, Malaysia	100.0	764	219
Sartorius Stedim Biotech (Beijing) Co. Ltd., Beijing, China	100.0	212	51
Sartorius Stedim Australia Pty. Ltd., East Oakleigh, Australia	100.0	1,179	1,331
Sartorius Stedim Nordic A/S, Taastrup, Denmark	100.0	222	300
Sartorius Stedim Systems GmbH, Melsungen, Germany, along with its subsidiary:	100.0	12,602	4,232
Sartorius Stedim India Pvt. Ltd., Bangalore, India	100.0	3,719	-131
Sartorius Stedim North America Inc., Edgewood, New York, USA, along with its subsidiaries:	100.0	36,945	1,779
Sartorius Stedim Systems Inc., Springfield, Missouri, USA	100.0	11,976	-2,354
Sartorius Stedim SUS Inc., Concord, California, USA	100.0	5,806	-337
Sartorius Mechatronics C&D GmbH & Co. KG, Aachen, Germany, including Sartorius-Verwaltungs-GmbH	100.0	-569	-653
Denver Instrument GmbH, Goettingen, Germany	100.0	1,453	178

	Ownership in %	Equity at Dec. 31, 2008 € in K	Net profit at Dec. 31, 2008 € in K
Sartorius Corporate Administration GmbH, Goettingen, Germany	100.0	1,393	156
Sartorius Mechatronics T&H Hamburg GmbH, Hamburg, Germany, along with its subsidiary:	100.0	20,211	3,155
Sartorius Mechatronics India Pvt. Ltd., Bangalore, India	100.0	4,094	662
Sartorius Mechatronics Austria GmbH, Vienna, Austria	100.0	3,122	447
Sartorius Mechatronics France S.A.S, Palaiseau, France	100.0	1,529	-18
Sartorius Mechatronics Italy S.r.L., Florence, Italy	100.0	1,151	177
Sartorius Mechatronics UK Ltd., Epsom, U.K.	100.0	1,806	396
Sartorius Mechatronics Spain, S.A.U., Madrid, Spain	100.0	67	-310
Sartorius Mechatronics Netherlands B.V., Nieuwegein, Netherlands, along with its subsidiary:	100.0	574	314
GWT Global Weighing Technologies B.V. Netherlands, Nieuwegein, Netherlands	100.0	5	-1
Sartorius Mechatronics Belgium B.V., Vilvoorde, Belgium	100.0	428	104
Sartorius Mechatronics Switzerland AG, Dietikon, Switzerland	100.0	7	201
Sartorius Mechatronics Poland sp. z.o.o., Poznan, Poland	100.0	343	-14
Sartorius Mechatronics Hungaria Kft., Budapest, Hungary	100.0	596	173
ZAO Sartogasm, St. Petersburg, Russia*)	51.0	748	54
OOO Sartorius ICR, St. Petersburg, Russia	100.0	0	0
America			
Sartorius North America Inc., Edgewood, New York, USA, along with its subsidiaries:	100.0	23,137	0
Sartorius Mechatronics Corporation, Edgewood, New York, USA	100.0	7,055	-285
Sartorius Mechatronics Canada Inc., Mississauga, Canada	100.0	183	17
Sartorius TCC Company, Arvada, Colorado, USA	100.0	7,224	121
Denver Instrument Inc., Denver, USA	100.0	560	16
Sartorius do Brasil Ltda., São Paulo, Brazil*)	100.0	-237	-76
Sartorius de México S.A. de C.V., Naucalpan, Mexico*)	99.0	336	41
Sartorius Argentina S.A., Buenos Aires, Argentina*)	99.0	353	100
Asia Pacific			
Sartorius Mechatronics Australia Pty. Ltd., East Oakleigh, Australia	100.0	52	61
Sartorius Mechatronics Japan K.K., Tokyo, Japan	100.0	7,403	1,082
Sartorius Mechatronics Singapore Pte. Ltd., Singapore	100.0	972	64
Sartorius Mechatronics Hong Kong Ltd., Kowloon, Hong Kong, along with its subsidiaries:	100.0	2,487	262
Sartorius Mechatronics Korea Ltd., Seoul, South Korea	100.0	1,549	253
Beijing Sartorius Instrument & System Engineering Co. Ltd., Beijing, China	100.0	4,494	-40
Sartorius Scientific Instruments (Beijing) Co. Ltd., Beijing, China	100.0	9,817	1,185
Sartorius Mechatronics Philippines Inc., Makati City, Philippines*)	100.0	195	26
Sartorius Mechatronics (Thailand) Co. Ltd., Bangkok, Thailand	49.0	80	-26
PT. Sartorius Mechatronics Indonesia, Jakarta, Indonesia*)	95.0	36	22

As the financial statements of the subsidiaries identified by an *) were not available at the time our financial statements were prepared, the information from the annual financial statements of 2007 were considered.

Executive Board and Supervisory Board

During Fiscal 2008 ¹⁾

Executive Board

Dr. rer. pol. Joachim Kreuzburg
Dipl.-Ingenieur (Graduate Engineer)
CEO and Chairman
Biotechnology Division and Finance
Born April 22, 1965
Hanover, Germany
Member since November 11, 2002
"Sprecher" (Spokesman) from
May 1, 2003, to Nov. 10, 2005
Chairman since Nov. 11, 2005
Appointed until November 10, 2010

Dr. rer. nat. Günther Maaz
Dipl.-Physiker (Graduate Physicist)
Executive for Labor Relations
Mechatronics Division
Human Resources and General
Administration
Born Sept. 13, 1949
Uslar, Germany
Member since November 11, 2002
Executive for Labor Relations since
Sept. 8, 2007
Appointed until November 10, 2010

Supervisory Board

Prof. Dr. Dres. h.c. Arnold Picot
Dipl.-Kaufmann (Graduate in
Business Administration) and
university professor
Chairman
CEO and Executive Director of the
Institute of Information,
Organization and Management,
Faculty of Economics at the
Ludwig Maximilian University
in Munich
Gauting, Germany

Gerd-Uwe Boguslawski
Dipl.-Sozialwirt (Graduate Social
Economist) Vice Chairman of the
Supervisory Board
1st senior local officer of the
German Metalworkers' Union in
southern Lower Saxony | Harz region
Northeim, Germany

Dr. Dirk Basting
Dipl.-Chemiker (Graduate Chemical
Engineer)
Fort Lauderdale, Florida, USA

Annette Becker
Kauffrau (Business Administrator)
Chairwoman of the Employees'
Council of Sartorius Corporate
Administration GmbH
Goettingen, Germany

Christiane Benner
Dipl.-Soziologin (Graduate
Sociologist) Gewerkschafts-
sekretärin (Union Secretary)
Executive Committee, German
Metalworkers' Union IG Metall
Frankfurt am Main, Germany

Uwe Bretthauer
Dipl.-Ingenieur (Graduate Engineer)
Chairman of the Employees'
Council of Sartorius AG
Goettingen, Germany

Michael Dohrmann
Feinmechaniker (Precision
Engineer)
Chairman of the Employees'
Council of Sartorius Stedim
Biotech GmbH
Goettingen, Germany

Dr. Lothar Kappich
Dipl.-Ökonom (Graduate
Economist)
Member since April 27, 2007
Managing Director of ECE
Projektmanagement GmbH & Co. KG
Hamburg, Germany

Prof. Dr. Gerd Krieger
Lawyer
Honorary Professor at the Heinrich-
Heine University in Duesseldorf
Duesseldorf, Germany

**Prof. Dr. rer. nat. Dr.-Ing. E.h.
Heribert Offermanns**
Dipl.-Chemiker (Graduate Chemical
Engineer)
Honorary Professor at the
Johann-Wolfgang-Goethe
University of Frankfurt am Main
Hanau, Germany

Dr. Michael Schulenburg
Dipl.-Ingenieur (Graduate Engineer)
Management Consultant
Mettmann, Germany

Manfred Werner
Dipl.-Ökonom (Graduate
Economist)
Senior Vice President of
Administration, Organization and
Auditing, Sartorius Corporate
Administration GmbH
Goettingen, Germany

Committees of the Supervisory Board

Executive Task Committee
Prof. Dr. Dres. h.c. Arnold Picot
(Chairman)
Gerd-Uwe Boguslawski
Uwe Bretthauer
Dr. Michael Schulenburg

Audit Committee
Dr. Michael Schulenburg (Chairman)
Gerd-Uwe Boguslawski
Uwe Bretthauer
Prof. Dr. Dres. h.c. Arnold Picot

Conciliation Committee
Prof. Dr. Dres. h.c. Arnold Picot
(Chairman)
Gerd-Uwe Boguslawski
Uwe Bretthauer
Dr. Michael Schulenburg

Nomination Committee
Prof. Dr. Gerd Krieger
Prof. Dr. Dres. h.c. Arnold Picot
Dr. Michael Schulenburg

Positions Held by Members of the Supervisory Board ¹⁾ as of December 31, 2008

Dr. rer. pol. Joachim Kreuzburg

Président-Directeur Général (CEO) of:

- Sartorius Stedim Biotech S.A., France ²⁾

On the Supervisory Board of:

- Sartorius Stedim Biotech GmbH, Germany, Vice Chairman ²⁾

On the Board of Directors of:

- Sartorius Stedim North America, Inc., USA ²⁾
- Sartorius Stedim SUS, Inc., USA ²⁾
- Sartorius Stedim Filters, Inc., Puerto Rico ²⁾
- Sartorius Mechatronics Japan K.K., Japan ²⁾
- Sartorius Stedim Japan K.K., Japan ²⁾
- Beijing Sartorius Instrument & System Engineering Co. Ltd., China ²⁾
- Sartorius Scientific Instruments (Beijing) Co. Ltd., China ²⁾
- Sartorius Stedim Lab Ltd., United Kingdom ²⁾

On the Landesbeirat (Regional Advisory Board) of:

- Commerzbank AG, Hamburg, Germany ³⁾

Dr. rer. nat. Günther Maaz

On the Board of Directors of:

- Sartorius North America, Inc., USA ²⁾
- Sartorius Mechatronics Corporation, USA ²⁾
- Sartorius TCC Company, USA ²⁾
- Denver Instrument, Inc., USA ²⁾
- Sartorius Stedim Filters, Inc., Puerto Rico ²⁾
- Sartorius Mechatronics Hong Kong Ltd., China ²⁾
- Beijing Sartorius Instrument & System Engineering Co. Ltd., China, Vice Chairman ²⁾
- Sartorius Scientific Instruments (Beijing) Co. Ltd., China, Vice Chairman ²⁾
- Sartorius Mechatronics India Pvt. Ltd., India ²⁾

Positions Held by Members of the Supervisory Board ¹⁾ as of December 31, 2008

Prof. Dr. Dres. h.c. Arnold Picot

On the Conseil d'Administration (Board of Directors) of:

- Sartorius Stedim Biotech S.A., France ²⁾

On the Supervisory Board of:

- Sartorius Stedim Biotech GmbH, Germany, Chairman ²⁾
- eteleon e-solutions AG, Germany, Vice Chairman ³⁾
- Takkt AG, Germany ³⁾
- Wissenschaftliches Institut für Infrastruktur und Kommunikationsdienste GmbH (Scientific Institute for Communication Services) and WIK-Consult GmbH, Germany ³⁾

Gerd-Uwe Boguslawski

On the Supervisory Board of:

- Novelis Deutschland GmbH, Germany ³⁾
- Demag Cranes & Components GmbH, Germany ³⁾
- Demag Cranes AG, Germany ³⁾

Dr. Dirk Basting

None

Annette Becker

None

Christiane Benner

On the Supervisory Board of:

- IAV GmbH, Ingenieurgesellschaft Auto und Verkehr, (auto and traffic engineering company), Germany ³⁾

Uwe Bretthauer

None

Michael Dohrmann

None

Dr. Lothar Kappich

None

Prof. Dr. Gerd Krieger

On the Supervisory Board of:

- ARAG Lebensversicherungs-AG, Germany ³⁾
- ARAG Krankenversicherungs-AG, Germany ³⁾

Prof. Dr. rer. nat. Dr.-Ing. E.h.

Heribert Offermanns

On the Supervisory Board of:

- Innovectis (Gesellschaft für Innovative Technologien und FuE-Dienstleistungen) GmbH, Germany, Vice Chairman ³⁾

Dr. Michael Schulenburg

On the Advisory Board of:

- Lohmann GmbH & Co. KG, Germany ³⁾
- Odenwald Faserplattenwerk GmbH, Germany ³⁾

On the Board of Directors of:

- Cognis Holding Luxembourg S.à.r.l., Luxembourg ³⁾

On the Supervisory Board of:

- Cognis GmbH, Germany, Chairman ³⁾

Manfred Werner

On the Consejo de Administración (Advisory Board) of:

- Sartorius Mechatronics Spain S.A., Spain, "Presidente" (President) ²⁾
- Sartorius Stedim Spain S.A., Spain, "Presidente" (President) ²⁾

On the Raad van Bestuur (Board of Management) of:

- Sartorius Mechatronics Belgium N.V., Belgium ²⁾
- Sartorius Stedim Belgium N.V., Belgium ²⁾

¹⁾ Information required pursuant to Section 285, No. 10, of the German Commercial Code (HGB)

²⁾ Positions held within the Group

³⁾ External positions held

About This Publication

Published by

Sartorius AG
Group Corporate Communications
37070 Goettingen, Germany

Editorial Deadline

March 10, 2009

Published on

March 11, 2009

Photography

Peter Ginter | Lohmar

Original German version published on
March 11, 2009

Sartorius AG
Weender Landstrasse 94-108
37075 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

www.sartorius.com