

## Sartorius AG 2009 Financial Statements





**Sartorius AG**  
2009 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. Amounts are reported in thousands, abbreviated as K, unless otherwise specified.

# 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, 2009 € in K	Dec. 31, 2008 € in K
<b>A. Fixed assets</b>			
I. Intangible assets	(4)	685	989
II. Plant, property and equipment	(5)	16,663	17,956
III. Financial assets	(6)	371,167	368,158
		<b>388,515</b>	<b>387,103</b>
<b>B. Current assets</b>			
I. Inventories	(7)	12,493	17,249
II. Trade and other receivables	(8)	27,152	40,677
III. Securities	(9)	16,082	15,069
IV. Cash on hand, deposits in banks		321	2,044
		<b>56,048</b>	<b>75,039</b>
<b>C. Prepaid expenses</b>	(10)	<b>1,002</b>	<b>1,429</b>
		<b>445,565</b>	<b>463,571</b>

Equity and Liabilities	Notes	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
<b>A. Equity</b>			
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)	23,082	27,810
		<b>168,475</b>	<b>173,203</b>
<b>B. Provisions</b>	(13)	<b>32,179</b>	<b>128,328</b>
<b>C. Liabilities</b>	(14)	<b>244,911</b>	<b>162,040</b>
		<b>445,565</b>	<b>463,571</b>

# Income Statement

	Notes	2009 € in K	€ in K	2008 € in K
1. Sales revenue	(18)	101,029		129,687
2. Decrease (-)   Increase in inventories of finished goods and work in progress	(19)	-2,607		447
			<b>98,422</b>	130,134
3. Work performed by the enterprise and capitalized	(20)	290		113
4. Other operating income	(21)	10,950		14,535
			<b>11,240</b>	14,648
			<b>109,662</b>	144,782
5. Raw materials and supplies	(22)	37,075		48,404
6. Employee benefits expense	(23)	42,136		50,176
7. Depreciation and amortization expense	(24)	3,259		3,459
8. Other operating expenses	(25)	44,089		41,689
9. Income from investments	(26)	24,064		18,699
10. Impairment losses on financial assets	(27)	14		1,527
			<b>102,509</b>	126,556
<b>11. Earnings before interest and taxes</b>			<b>7,153</b>	18,226
12. Interest and similar income	(28)	292		914
13. Interest and similar expenses	(28)	5,123		11,142
			<b>-4,831</b>	-10,228
<b>14. Profit before tax</b>			<b>2,322</b>	7,998
15. Income tax expense	(29)	-87		233
16. Other taxes		148		143
			<b>61</b>	376
<b>17. Net profit</b>			<b>2,261</b>	7,622
18. Profit brought forward			<b>20,821</b>	20,188
<b>19. Retained profit</b>			<b>23,082</b>	27,810

# Notes to the Financial Statements

Cash Flow Statement	2009 € in K	2008 € in K
Net profit for the period	2,261	7,621
Depreciation and amortization of fixed assets	3,259	3,459
Decrease (-)   Increase in long-term provisions	-755	554
Other material non-cash income (-)   expenses (+)	-10,869	-10,779
<b>Cash earnings acc. to the DVFA   SG*</b>	<b>-6,104</b>	<b>855</b>
Decrease (-) in short-term provisions	-1,195	-2,281
Gains resulting from the disposal of fixed assets	271	-1
Decrease   Increase (-) in inventories	4,756	-631
Decrease   Increase (-) in short- and long-term receivables and other assets	13,952	-4,692
Increase   Decrease (-) in liabilities (excluding liabilities to banks)	14,305	-3,871
<b>Cash flows from operating activities</b>	<b>25,985</b>	<b>-10,621</b>
Proceeds from fixed asset disposals	35	116
Payments (-) for intangible assets	-325	-388
Payments (-) for property, plant and equipment	-1,643	-3,597
Payments (-) for financial assets	-107,352	-7,239
<b>Cash flows from investing activities</b>	<b>-109,285</b>	<b>-11,108</b>
Dividend payments (-)	-6,989	-11,421
Additions to   Repayment of (-) financial liabilities	88,566	30,306
<b>Cash flows from financing activities</b>	<b>81,577</b>	<b>18,885</b>
<b>Change in cash and cash equivalents</b>	<b>-1,723</b>	<b>-2,843</b>
<b>Cash and cash equivalents at beginning of period</b>	<b>2,044</b>	<b>4,887</b>
<b>Cash and cash equivalents at end of period</b>	<b>321</b>	<b>2,044</b>
	<b>Dec. 31, 2009 € in K</b>	<b>Dec. 31, 2008 € in K</b>
<b>Composition of cash and cash equivalents</b>		
Cash on hand, deposits in banks	321	2,044

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).

The same accounting and valuation methods were applied as in the previous year.

## 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.

## 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, 2009	3,489	30	3,519
Investments	326	0	326
Disposals	0	0	0
Transfers	34	-30	4
Gross book values as of Dec. 31, 2009	3,849	0	3,849
Amortization as of Jan. 1, 2009	2,530	0	2,530
Amortization in 2009	633	0	633
Disposals	0	0	0
Amortization as of Dec. 31, 2009	3,163	0	3,163
Net book values as of Dec. 31, 2009	686	0	<b>686</b>
Net book values as of Dec. 31, 2008	959	30	989

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.

## 3. Currency Translation

Short-term receivables and liabilities in foreign currencies were translated at the year-end rates on December 31, 2009.

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

	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
Property, plant and equipment developed as follows:					
Gross book values as of Jan. 1, 2009	26,100	12,240	25,593	484	64,417
Investments	0	75	1,434	133	1,642
Disposals	814	115	874	0	1,803
Transfers	-8	21	465	-482	-4
Gross book values as of Dec. 31, 2009	25,278	12,221	26,618	135	64,252
Depreciation as of Jan. 1, 2009	14,984	10,554	20,923	0	46,461
Depreciation in 2009	628	456	1,542	0	2,626
Disposals	738	115	645	0	1,498
Transfers	-8	0	8		
Depreciation as of Dec. 31, 2009	14,866	10,895	21,828	0	47,589
Net book values as of Dec. 31, 2009	10,412	1,326	4,790	135	<b>16,663</b>
Net book values as of Dec. 31, 2008	11,116	1,686	4,670	484	17,956

Property, plant and equipment are reported at acquisition cost or production cost and, if subject to depreciation, are depreciated as scheduled.

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

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

Buildings	15 to 50 years
Machinery	5 to 15 years
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, 2009	369,006	383	369,389
Investments	3,009	14	3,023
Disposals	0	0	0
Transfers	0	0	0
Gross book values as of Dec. 31, 2009	372,015	397	372,412
Impairment losses as of Jan. 1, 2009	848	383	1,231
Impairment losses in 2009	0	14	14
Disposals	0	0	0
Reversals	0	0	0
Impairment losses as of Dec. 31, 2009	848	397	1,245
Net book values as of Dec. 31, 2009	371,167	0	<b>371,167</b>
Net book values as of Dec. 31, 2008	368,158	0	368,158

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 Sartorius Stedim Biotech S.A. listed on the French stock exchange and headquartered in Aubagne, France. 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 the Stedim S.A. shares and had been the company's majority shareholder until then. Sartorius AG subsequently 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 and 2009, Sartorius AG acquired additional 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, 2009, this deal as a whole entailed acquisition costs for the shares in the two companies amounting to €309,523 K. Of this amount, shares in VL Finance S.A. accounted for €151,520 K, and shares in Sartorius Stedim Biotech S.A. €158,003 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 approximately €89 million at December 31, 2009, due to a decrease in the prices 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.

Sartorius AG's investments in financial assets essentially resulted from purchases of shares in Sartorius Stedim Biotech S.A., Aubagne, France. In the reporting year, Sartorius AG acquired 141,590 shares, which represent 0.7% of Sartorius Stedim Biotech S.A.'s share capital.

Please refer to "Major Affiliates and Ownership."

## 7. Inventories

	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
Raw materials and supplies	5,314	7,177
Work in progress	3,676	5,638
Finished goods and merchandise	3,200	4,186
Payments on account	303	248
	<b>12,493</b>	<b>17,249</b>

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 the risk of limited usability, inventories are marked down accordingly.

## 8. Trade and Other Receivables

	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
Trade receivables	3,763	7,797
Receivables from affiliated companies	20,733	23,664
Receivables from companies in which investments are held	0	1
Other assets	2,656	9,216
	<b>27,152</b>	<b>40,678</b>
Of which due in more than one year:		
Trade receivables	0	0
Other assets	1,332	1,198
	<b>1,332</b>	<b>1,198</b>

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.

The reduction in trade receivables is essentially due to implementation of a factoring program as of June 30, 2009. Within the scope of this program, liabilities of €3,115 K were sold to a factor as of December 31, 2009.

## 9. Securities

	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
Own (treasury) shares	16,082	15,069

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 2009.

Because of the increase in the prices of Sartorius shares, additions of €1,013 K were reported for treasury shares in fiscal 2009.

## 10. Prepaid Expenses

	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
Loan discounts	960	1,168
Other prepaid expenses	42	261
	<b>1,002</b>	<b>1,429</b>

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 €1.00.

## 12. Earnings Reserves and Retained Profits

	€ in K
The reserves for own (treasury) shares of Sartorius AG developed as follows in fiscal 2009:	
At Jan. 1, 2009	15,069
Addition acc. to § 272, Subsec. 4, of HGB*	1,013
At Dec. 31, 2009	<b>16,082</b>

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 in fiscal 2009 as follows:	
At Jan. 1, 2009	10,207
Reclassified into the reserves for own shares	1,013
At Dec. 31, 2009	<b>9,194</b>

The retained profit of of €23,082 K consists of the profit brought forward of €20,821 K and of the net profit of €2,261 K.

\* HGB = Handelsgesetzbuch = German Commercial Code

### 13. Provisions

	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
Provisions for retirement benefits and similar obligations	14,614	14,721
Tax provisions	341	494
Deferred tax provisions	0	250
Other provisions	17,224	112,863
	<b>32,179</b>	<b>128,328</b>

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 €8.7 million of the reserves set up on December 31, 2009, is allocated to Sartorius AG's general in-house retirement plan, and €6.0 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, 2009 € in K	Dec. 31, 2008 € in K
The other provisions include the following amounts:		
Price warrants	0	103,691
Restructuring provision	10,130	0
Employee benefits expense	2,773	3,966
Payments to employees on the early retirement plan for offsetting reduced work hours	1,819	2,432
Commissions, credit notes, bonuses	359	402
Warranties	285	220
Other	1,858	2,152
	<b>17,224</b>	<b>112,863</b>

## 14. Liabilities

This item consists of the following:	Balance sheet Dec. 31, 2009 € in K	Remaining term of more than five years € in K	Remaining term of up to one year € in K	Balance sheet Dec. 31, 2008 € in K	Remaining term of more than five years € in K	Remaining term of up to one year € in K
Liabilities to banks	162,123	0	4,784	73,557	0	11,640
Payments received on account of orders	197	0	197	398	0	398
Trade payables	5,582	0	5,582	5,348	0	5,348
Payables to affiliated companies	72,984	0	72,984	81,148	0	81,148
Payables to companies in which investments are held	39	0	39	0	0	0
Other liabilities	3,986	0	3,986	1,589	0	1,589
	<b>244,911</b>	<b>0</b>	<b>87,572</b>	<b>162,040</b>	<b>0</b>	<b>100,123</b>

Other liabilities relate to the following:	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
Taxes	646	801
Liabilities relating to social security	0	0

Liabilities are reported as the amounts repayable.

## 15. Contingent Liabilities

Guarantees (suretyships) and other contingent liabilities	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
	18,735	35,841

Guarantees (suretyships) exist in the amount of €18,607 K granted to affiliated companies that are included in the consolidated financial statements.

## 16. Other Financial Obligations

	Dec. 31, 2009 € in K	Dec. 31, 2008 € in K
In addition to provisions, liabilities and contingent liabilities, other financial obligations consist of the following:		
Rental and leasing contracts		
– due in fiscal 2010 (previous year: due in fiscal 2009)	1,271	1,241
– due in any one fiscal year from 2011 to 2014 (previous year: due in any one fiscal year from 2010 to 2013)	1,051	1,302
	<b>2,322</b>	<b>2,543</b>

There are no other financial obligations to affiliated companies.

## 17. Derivative Financial Instruments

Within the scope of its operations, the Sartorius Group uses derivative financial instruments to hedge against risks arising from changes in interest and foreign exchange rates. The Group does not use 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 for the Group 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 concluded forward exchange transactions during 2009 up to the reporting date. For existing forward exchange transactions, a reserve of €55 K was set up in the reporting year for anticipated losses related to incomplete contracts.

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, 2009, Sartorius AG had two interest caps with an interest rate ceiling of 4.5%, one against the 3-month EURIBOR and the other against the 6-month EURIBOR, 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 €0 K as of December 31, 2009.

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

	Nominal value Dec. 31, 2009 € in K	Nominal value Dec. 31, 2008 € in K
Forward exchange transactions	9,308	0
Currency option transactions	0	55,190
	<b>9,308</b>	<b>55,190</b>

	Fair value Dec. 31, 2009 € in K	Fair value Dec. 31, 2008 € in K
Forward exchange transactions	9,253	0
Currency option transactions	0	-3,640
	<b>9,253</b>	<b>-3,640</b>

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

# Notes to the Income Statement

## 18. Sales Revenue

	2009 € in K	2008 € in K
Sales revenue, which is broken down by geographical markets, is as follows:		
Germany	39,714	50,843
All other countries	61,315	78,844
Total	<u>101,029</u>	<u>129,687</u>

## 19. Decrease | Increase in Inventories of Finished Goods and Work in Progress

	2009 € in K	2008 € in K
Finished goods	-645	-330
Work in progress	-1,962	777
	<u>-2,607</u>	<u>447</u>

## 20. Work Performed by the Enterprise and Capitalized

	2009 € in K	2008 € in K
This item is apportioned as follows to the items under "Property, plant and equipment":		
Machinery	0	25
Construction in progress	59	55
Factory and office equipment	199	0
Tools and equipment	32	33
	<u>290</u>	<u>113</u>



## 21. Other Operating Income

	2009 € in K	2008 € in K
Income due to currency translation	2,505	5,336
Income from affiliated companies	4,070	4,585
Income from the release of provisions	1,012	1,644
Government grants, promotions and other subsidies	713	993
Income from reversal of write-downs	508	423
Reversal of write-downs on financial assets	1,013	0
Other income	1,129	1,554
	<b>10,950</b>	<b>14,535</b>

## 22. Raw Materials and Supplies

	2009 € in K	2008 € in K
Cost of raw materials and supplies of purchased merchandise	36,048	46,902
Cost of purchased services	1,027	1,502
	<b>37,075</b>	<b>48,404</b>

## 23. Employee Benefits Expense

	2009 € in K	2008 € in K
Wages and salaries	34,014	41,041
Social security	7,034	7,274
Expenses for retirement benefits and pensions	1,088	1,861
	<b>42,136</b>	<b>50,176</b>

Sartorius AG employed an average of 765 people in the reporting year (2008: 796).

## 24. Depreciation and Amortization Expense

	2009 € in K	2008 € in K
Amortization of intangible assets and depreciation of property, plant and equipment		
Scheduled	<b>3,259</b>	<b>3,459</b>

## 25. Other Operating Expenses

	2009 € in K	2008 € in K
Costs for outside services	12,281	15,890
Costs for restructuring	11,448	0
Commissions and license fees	4,672	5,215
Expenses due to currency translation	3,103	6,614
Car and leasing expenses	1,896	2,120
Legal and consulting expenses including patent costs	1,776	1,270
Freight and packaging costs	1,255	2,058
Advertising expenses	1,135	1,608
Itemized allowance for bad debts and lump-sum valuation adjustment on receivables	1,073	1,553
Travel and entertainment expenses	1,024	1,315
Maintenance costs	1,007	1,402
Rent and leasing expenses	612	560
Office, mail and other administrative expenses	318	458
Insurance expenses	160	166
Other finance charges	106	152
Other expenses	2,223	1,308
	<b>44,089</b>	<b>41,689</b>

## 26. Income from Investments

	2009 € in K	2008 € in K
Income from investments	24,064	18,699

As in the year-earlier reporting period, the entire amount of Sartorius AG's income from investments consists of dividend payments from subsidiaries.

## 27. Impairment Losses on Financial Assets

	2009 € in K	2008 € in K
Impairment losses on financial assets	14	514
Impairment losses on own (treasury) shares	0	1,013
	<b>14</b>	<b>1,527</b>

## 28. Interest

	2009 € in K	2008 € in K
Interest and similar income	292	914
– of which from affiliated companies	(241)	(593)
Interest and similar expenses	5,123	11,142
– of which to affiliated companies	(537)	(3,758)
	<u>-4,831</u>	<u>-10,228</u>

The improvement in the interest figure resulted particularly from the lowered interest rates for bank loans and from the decrease in the interest-bearing liabilities to affiliated companies.

## 29. Income Tax Expense

	2009 € in K	2008 € in K
Current income taxes	106	228
Deferred taxes	-250	0
Income taxes of the previous years	57	5
	<u>-87</u>	<u>233</u>

The "Other taxes" item essentially discloses property tax and motor vehicle tax.

## 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 10, 2009, and made available to the shareholders on the company's Internet site at [www.sartorius.com](http://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:

	2009 € in K	2008 € in K
Auditing of the annual financial statements and of the consolidated financial statements	122	122
Other independent reporting services	93	68
Tax consultation including preparation of tax returns	12	52
Other services	118	93
	<u>345</u>	<u>335</u>

## Total Remuneration of the Supervisory Board and of the Executive Board

	2009 € in K	2008 € in K
<b>Remuneration of the Supervisory Board Members</b>		
Total remuneration	573	523
- Fixed remuneration	293	300
- Performance-related remuneration	73	75
- Compensation for committee work	60	60
- Meeting attendance fee	147	88
<b>Prof. Dr. Dres. h.c. Arnold Picot (Chairman)</b>		
Total remuneration	119	108
- Fixed remuneration	60	60
- Performance-related remuneration	15	15
- Compensation for committee work	18	18
- Meeting attendance fee	26	15
<b>Gerd-Uwe Boguslawski* (Vice Chairman)</b>		
Total remuneration	88	77
- Fixed remuneration	40	40
- Performance-related remuneration	10	10
- Compensation for committee work	12	12
- Meeting attendance fee	26	15
<b>Dr. Dirk Basting</b>		
Total remuneration	30	29
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Meeting attendance fee	5	4
<b>Annette Becker*</b>		
Total remuneration	30	29
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Meeting attendance fee	5	4
<b>Christiane Benner*</b>		
Total remuneration	30	29
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Meeting attendance fee	5	4
<b>Uwe Bretthauer*</b>		
Total remuneration	63	52
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Compensation for committee work	12	12
- Meeting attendance fee	26	15

	2009 € in K	2008 € in K
<b>Michael Dohrmann*</b>		
Total remuneration	31	29
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Meeting attendance fee	6	4
<b>Dr. Lothar Kappich</b>		
Total remuneration	31	28
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Meeting attendance fee	6	3
<b>Prof. Dr. Gerd Krieger</b>		
Total remuneration	31	28
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Meeting attendance fee	6	3
<b>Prof. Dr. rer. nat. Dr.-Ing. Heribert Offermanns</b>		
Total remuneration	31	28
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Meeting attendance fee	6	3
<b>Dr. Michael Schulenburg</b>		
Total remuneration	68	57
- Fixed remuneration	20	20
- Performance-related remuneration	5	5
- Compensation for committee work	18	18
- Meeting attendance fee	25	14
<b>Manfred Werner**</b>		
Total remuneration	21	29
- Fixed remuneration	13	20
- Performance-related remuneration	3	5
- Meeting attendance fee	5	4

\* The employee representatives declared that they donate their Supervisory Board remuneration to the foundation Hans Böckler Stiftung according to the guidelines of the German Trade Union Association.

\*\* His position remained dormant in the period from August 27, 2009, to March 7, 2010.

	2009 € in K	2008 € in K
<b>Remuneration of the Executive Board Members</b>		
Total remuneration	1,390	1,362
- Performance-independent remuneration*	820	806
- Performance-related remuneration	570	556
- Phantom stock units paid out	0	0
- Fair value of the phantom stock units apportioned in the reporting year	413	55
- Change in the fair value of the phantom stock units apportioned in the previous years	192	-399
- Fair value of mid-term remuneration components	0	63
- Additions to provisions for employee benefits	66	82
<b>Dr. Joachim Kreuzburg</b>		
Total remuneration	726	789
- Performance-independent remuneration*	418	480
- Performance-related remuneration	308	309
- Phantom stock units paid out	0	0
- Fair value of the phantom stock units apportioned in the reporting year	242	36
- Change in the fair value of the phantom stock units apportioned in the previous years	121	-262
- Fair value of mid-term remuneration components	0	63
- Additions to provisions for employee benefits	31	42

\* In fiscal 2009, the Executive Board waived parts of its performance-independent remuneration.

\*\* up to his retirement

	2009 € in K	2008 € in K
<b>Jörg Pfirrmann (since July 24, 2009)</b>		
Total remuneration	144	
- Performance-independent remuneration*	84	
- Performance-related remuneration	60	
- Phantom stock units paid out	0	
- Fair value of the phantom stock units apportioned in the reporting year	39	
- Change in the fair value of the phantom stock units apportioned in the previous years	-	
<b>Reinhard Vogt (since July 24, 2009)</b>		
Total remuneration	225	
- Performance-independent remuneration*	136	
- Performance-related remuneration	89	
- Phantom stock units paid out	0	
- Fair value of the phantom stock units apportioned in the reporting year	58	
- Change in the fair value of the phantom stock units apportioned in the previous years	-	
<b>Dr. Günther Maaz (until July 24, 2009)</b>		
Total remuneration**	295	573
- Performance-independent remuneration*	182	326
- Performance-related remuneration	113	247
- Phantom stock units paid out	0	0
- Fair value of the phantom stock units apportioned in the reporting year	74	19
- Change in the fair value of the phantom stock units apportioned in the previous years	71	-137
- Additions to provisions for employee benefits	35	40

Components with a Long-term Incentive Effect	Number of phantom stock units	Allotment price* €	Fair value when granted on Jan. 1 of the particular year € in K	Fair value at year-end on Dec. 31, 2009 € in K	Paid out € in K	Exercisable
<b>Dr. Joachim Kreuzburg</b>						
- Tranche of phantom stock units for fiscal 2005	4,053	15.42	63	64	0	No
- Tranche of phantom stock units for fiscal 2006	3,871	21.31	83	61	0	No
- Tranche of phantom stock units for fiscal 2007	3,593	34.79	125	55	0	No
- Tranche of phantom stock units for fiscal 2008	4,754	28.92	138	71	0	No
- Tranche of phantom stock units for fiscal 2009	16,851	8.16	138	242	0	No
	<b>33,122</b>		<b>547</b>	<b>493</b>		
<b>Jörg Pfirrmann</b>						
- Tranche of phantom stock units for fiscal 2009	2,703	8.16	22	39	0	No
<b>Reinhard Vogt</b>						
- Tranche of phantom stock units for fiscal 2009	4,054	8.16	33	58	0	No
<b>Dr. Günther Maaz until July 24, 2009</b>						
- Tranche of phantom stock units for fiscal 2005	2,594	15.42	40	41	0	No
- Tranche of phantom stock units for fiscal 2006	2,580	21.31	55	41	0	No
- Tranche of phantom stock units for fiscal 2007	1,796	34.79	63	27	0	No
- Tranche of phantom stock units for fiscal 2008	2,507	28.92	73	37	0	No
- Tranche of phantom stock units for fiscal 2009	5,162	8.16	42	74	0	No
	<b>14,639</b>		<b>273</b>	<b>220</b>		

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 replaces the remuneration element that was dependent upon the share price and had been paid out annually until then. As of 2005, the 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 2009 Annual Report for the Sartorius Group.

	2009 € in K	2008 € in K
<b>Former managing directors' remuneration and pension obligations</b>		
Remuneration of former managing directors and members of the Executive Board as well as of their surviving dependents	1,084	256
Pension obligations to former managing directors and members of the Executive Board as well as to their surviving dependents	4,991	4,683

### 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 €1,673 K (2008: €2,229 K). This amount includes income from the reversal of valuation allowances and the release of provisions. Sartorius AG's non-periodic expenses are €12,674 K (2008: €458 K) and concern restructuring measures in particular.

### Number of Employees

	2009	2008
Average workforce employed during the fiscal year:		
Employees	765	796
	765	796

### Proposal for Appropriation of Profits

The Executive Board will submit a proposal to the Annual General Shareholders' Meeting to appropriate the retained profit of €23,081,595.25 reported

by Sartorius AG for the year ended December 31, 2009, 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	16,092,385.71
	23,081,595.25

### 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 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.

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

The following information was available according to Section 21, Subsection 1, or 1a, of the German Securities Trading Act (WpHG) as of December 31, 2009:

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 19, 2010

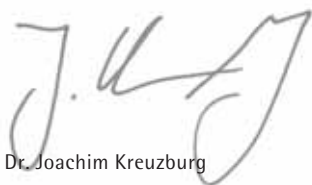
Sartorius Aktiengesellschaft  
The Executive Board

### Declaration of the Executive Board

To the best of our knowledge, we declare that the year-end financial statements for fiscal 2009 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 19, 2010

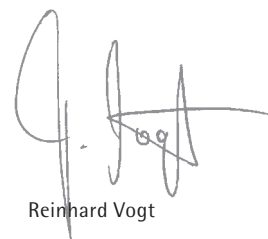
Sartorius Aktiengesellschaft  
The Executive Board



Dr. Joachim Kreuzburg



Jörg Pfirrmann



Reinhard Vogt

# 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, 2009. 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 annual 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 annual 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, annual 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 annual 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 19, 2010

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



Jürgen Reker  
Auditor



Dieter Tenambergen  
Auditor



# Major Affiliates and Ownership

	Ownership in percent %	Equity at Dec. 31, 2009 € in K	Net profit at Dec. 31, 2009 € in K
<b>Europe</b>			
<b>VL Finance S.A.S., Aubagne, France</b>	100.0	37,567	823
<b>Sartorius Stedim Biotech S.A., Aubagne, France along with its subsidiaries:</b>	71.1	84,320	14,159
Sartorius Stedim Aseptics S.A., Lourdes, France	100.0	3,611	1,787
Integrated Biosystems Sàrl, Fribourg, Switzerland	100.0	1,604	-152
Sartorius Stedim SUS S.A.R.L., M'Hamdia, Tunisia	100.0	787	596
<b>Sartorius Stedim Biotech GmbH, Goettingen, Germany along with its subsidiaries:</b>	100.0	71,353	16,266
Sartorius Stedim UK Ltd., Epsom, U.K.	100.0	930	1,007
Wave Biotech AG, Tagelswangen, Switzerland	100.0	1,567	98
Sartorius Stedim France S.A.S, Aubagne, France	100.0	3,017	1,996
Sartorius Stedim Italy S.p.A., Florence, Italy	100.0	3,625	646
Sartorius Stedim Spain S.A., Madrid, Spain	100.0	-645	117
Sartorius Stedim Hungaria Kft, Budapest, Hungary	100.0	94	93
Sartorius Stedim Poland Sp. z.o.o., Poznan, Poland	100.0	205	10
Sartorius Stedim Netherlands B.V., Nieuwegein, Netherlands	100.0	454	441
Sartorius Stedim Belgium N.V., Vilvoorde, Belgium	100.0	1,002	965
Sartorius Stedim Austria GmbH, Vienna, Austria	100.0	846	492
Sartorius Stedim Switzerland GmbH, Dietikon, Switzerland	100.0	40	-24
Sartorius Stedim Japan K.K., Tokyo, Japan	100.0	1,749	812
Sartorius Stedim Singapore Pte. Ltd., Singapore	100.0	725	275
Munktell & Filtrak GmbH, Baerenstein, Germany *)	49.0	3,997	743
Distribo GmbH, Goettingen, Germany *)	26.0	828	146
Sartorius Stedim Filters Inc., Yauco, Puerto Rico	100.0	5,282	3,503
Sartorius Stedim Plastics GmbH, Goettingen, Germany	100.0	2,348	1,060
Sartorius Technologies & Services GmbH, Goettingen, Germany	100.0	112	11
Sartorius Stedim Lab Ltd., Louth, U.K.	100.0	1,391	736
Sartorius Stedim Malaysia Sdn. Bhd., Kuala Lumpur, Malaysia	100.0	844	350
Sartorius Stedim Biotech (Beijing) Co. Ltd., Beijing, China	100.0	790	619
Sartorius Stedim Australia Pty. Ltd., East Oakleigh, Australia	100.0	1,343	992
Sartorius Stedim Nordic A/S, Taastrup, Denmark	100.0	208	-15
<b>Sartorius Stedim Systems GmbH, Melsungen, Germany along with its subsidiary:</b>	100.0	14,140	1,538
Sartorius Stedim India Pvt. Ltd., Bangalore, India	100.0	3,770	12
<b>Sartorius Stedim North America Inc., Edgewood, New York, USA along with its subsidiary:</b>	100.0	30,439	-5,587
Sartorius Stedim SUS Inc., Concord, California, USA	100.0	6,316	705
<b>Sartorius Mechatronics C&amp;D GmbH &amp; Co. KG, Aachen, Germany, including Sartorius-Verwaltungs-GmbH</b>	100.0	-3,312	-2,743
<b>Denver Instrument GmbH, Goettingen, Germany</b>	100.0	1,394	-59

	Ownership in percent %	Equity at Dec. 31, 2009 € in K	Net profit at Dec. 31, 2009 € in K
<b>Sartorius Corporate Administration GmbH, Goettingen, Germany</b>	100.0	129	-1,264
<b>Sartorius Mechatronics T&amp;H Hamburg GmbH, Hamburg, Germany along with its subsidiaries:</b>	100.0	17,902	-2,309
Sartorius Mechatronics India Pvt. Ltd., Bangalore, India	100.0	4,617	477
Sartorius Mechatronics Austria GmbH, Vienna, Austria	100.0	2,913	238
Sartorius Mechatronics France S.A.S, Palaiseau, France	100.0	668	-861
Sartorius Mechatronics Italy S.r.L., Florence, Italy	100.0	936	-208
Sartorius Mechatronics UK Ltd., Epsom, U.K.	100.0	1,856	-90
Sartorius Mechatronics Spain, S.A.U., Madrid, Spain	100.0	-912	-979
<b>Sartorius Mechatronics Netherlands B.V., Nieuwegein, Netherlands along with its subsidiary:</b>	100.0	-520	-1,094
GWT Global Weighing Technologies B.V. Netherlands, Nieuwegein, Netherlands	100.0	5	0
Sartorius Mechatronics Belgium B.V., Vilvoorde, Belgium	100.0	-71	-501
Sartorius Mechatronics Switzerland AG, Dietikon, Switzerland	100.0	-195	-198
Sartorius Mechatronics Sp. z.o.o., Poznan, Poland	100.0	334	-26
Sartorius Mechatronics Hungaria Kft., Budapest, Hungary	100.0	599	177
ZAO Sartogasm, St. Petersburg, Russia *)	51.0	-7	-114
OOO Sartorius ICR, St. Petersburg, Russia	100.0	0	0
<b>America</b>			
<b>Sartorius North America Inc., Edgewood, New York, USA along with its subsidiaries:</b>	100.0	22,448	0
Sartorius Mechatronics Corporation, Edgewood, New York, USA	100.0	7,550	-618
Sartorius Mechatronics Canada Inc., Mississauga, Canada	100.0	233	23
Sartorius TCC Company, Arvada, Colorado, USA	100.0	5,961	-1,085
Sartorius do Brasil Ltda., São Paulo, Brazil *)	100.0	-266	-92
Sartorius de México S.A. de C.V., Naucalpan, Mexico *)	99.0	109	-203
Sartorius Argentina S.A., Buenos Aires, Argentina *)	99.0	319	-20
<b>Asia   Pacific</b>			
Sartorius Mechatronics Australia Pty. Ltd., East Oakleigh, Australia	100.0	275	189
Sartorius Mechatronics Japan K.K., Tokyo, Japan	100.0	7,454	1,044
Sartorius Mechatronics Singapore Pte. Ltd., Singapore	100.0	1,109	139
<b>Sartorius Mechatronics Hong Kong Ltd., Kowloon, Hong Kong, China along with its subsidiary:</b>	100.0	2,984	589
Sartorius Mechatronics Korea Ltd., Seoul, South Korea	100.0	1,967	329
Beijing Sartorius Instrument & System Engineering Co. Ltd., Beijing, China	100.0	4,378	19
Sartorius Scientific Instruments (Beijing) Co. Ltd., Beijing, China	100.0	10,647	1,160
Sartorius Mechatronics Philippines Inc., Makati City, Philippines *)	100.0	268	69
Sartorius Mechatronics (Thailand) Co. Ltd., Bangkok, Thailand	49.0	118	37
PT. Sartorius Mechatronics Indonesia, Jakarta, Indonesia *)	95.0	19	5

As the financial statements of the subsidiaries identified by an \*) were not available at the time our annual financial statements were prepared, the information from the annual financial statements of 2008 were considered.

# Executive Board and Supervisory Board

During Fiscal 2009<sup>1)</sup>

## Executive Board

**Dr. rer. pol. Joachim Kreuzburg**  
Dipl.-Ingenieur (Graduate Engineer)  
CEO and Chairman  
Executive for Labor Relations  
Operations, Human Resources,  
Legal Issues, Internal Auditing and  
Corporate Communications  
Born April 22, 1965  
Hanover, Germany  
Member since November 11, 2002  
Sprecher (Spokesman) from May 1,  
2003, to November 10, 2005  
Chairman and CEO since  
November 11, 2005  
Executive for Labor Relations since  
July 24, 2009  
Appointed until November 10, 2015

**Jörg Pfirrmann**  
Dipl.-Ökonom (Graduate Economist)  
Finances, IT and  
General Administration  
Born November 30, 1972  
Nörten-Hardenberg, Germany  
Member since July 24, 2009  
Appointed until July 23, 2012

**Reinhard Vogt**  
Industriekaufmann (Industrial  
Business Manager)  
Marketing, Sales and Service  
Born August 4, 1955  
Dransfeld, Germany  
Member since July 24, 2009  
Appointed until July 23, 2014

**Dr. rer. nat. Günther Maaz**  
Dipl.-Physiker (Graduate Physicist)  
Member and Executive for Labor  
Relations until July 24, 2009  
Mechatronics Division,  
Human Resources and General  
Administration  
Born September 13, 1949  
Uslar, Germany  
Member since November 11, 2002  
Executive for Labor Relations since  
September 8, 2007

## Supervisory Board

**Prof. Dr. Dres. h.c. Arnold Picot**  
Dipl.-Kaufmann (Graduate in  
Business Administration)  
University professor  
Chairman of the Supervisory Board  
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  
1<sup>st</sup> Senior Officer of the German  
Metalworkers' Union  
in the 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)  
Gewerkschaftssekretärin  
(Union Secretary)  
German Metalworkers' Union  
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)  
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 in  
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  
Internal 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 the Members of the Executive Board<sup>1)</sup> as of December 31, 2009

**Dr. rer. pol. Joachim Kreuzburg**  
Président-Directeur Général (CEO)  
of:  
– Sartorius Stedim Biotech S.A.,  
France<sup>2)</sup>  
On the Supervisory Board of:  
– Sartorius Stedim Biotech GmbH,  
Germany; Vice Chairman<sup>2)</sup>  
On the Board of Directors of:  
– Sartorius Stedim North America,  
Inc., USA<sup>2)</sup>  
– Sartorius Stedim SUS, Inc., USA<sup>2)</sup>  
– Sartorius Stedim Filters, Inc.,  
Puerto Rico<sup>2)</sup>

- Sartorius Mechatronics Japan K.K., Japan<sup>2)</sup>
- Sartorius Stedim Japan K.K., Japan<sup>2)</sup>
- Beijing Sartorius Instrument & System Engineering Co. Ltd., China<sup>2)</sup>
- Sartorius Scientific Instruments (Beijing) Co. Ltd., China<sup>2)</sup>
- Sartorius Mechatronics Hong Kong Ltd., China<sup>2)</sup>
- Sartorius Stedim Lab Ltd., United Kingdom<sup>2)</sup>

On the Regionalbeirat (Regional Advisory Board) of:

- Commerzbank AG, Hamburg, Germany<sup>3)</sup>

On the Beirat (Advisory Board) of:

- Hameln Group GmbH, Germany<sup>3)</sup>

On the Wirtschaftsbeirat (Economic Advisory Board) of:

- Norddeutsche Landesbank, Germany<sup>3)</sup>

#### Jörg Pfirrmann

On the Verwaltungsrat (Administrative Board) of:

- Sartorius Mechatronics Switzerland AG, Switzerland<sup>2)</sup>

On the Board of Directors of:

- Sartorius North America, Inc., USA<sup>2)</sup>
- Sartorius Mechatronics Corporation, USA<sup>2)</sup>
- Sartorius TCC Company, USA<sup>2)</sup>
- Sartorius Stedim UK Ltd., United Kingdom<sup>2)</sup>

On the Comité Exécutif (Executive Committee) of:

- Sartorius Stedim France S.A.S., France<sup>2)</sup>

On the Consiglio di Amministrazione (Board of Directors) of:

- Sartorius Stedim Italy S.p.A., Italy<sup>2)</sup>

On the Consejo de Administracion (Board of Directors) of:

- Sartorius Stedim Spain S.A.<sup>2)</sup>

On the Supervisory Board of:

- Sartorius Stedim Nordic A/S<sup>2)</sup>

#### Reinhard Vogt

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

- Sartorius Stedim Biotech S.A., France<sup>2)</sup>

On the Board of Directors of:

- Sartorius Stedim North America, Inc., USA<sup>2)</sup>
- Sartorius Stedim SUS, Inc., USA<sup>2)</sup>
- Beijing Sartorius Instrument & System Engineering Co. Ltd., China<sup>2)</sup>
- Sartorius Scientific Instruments (Beijing) Co. Ltd., China<sup>2)</sup>
- Sartorius Stedim Biotech (Beijing) Co. Ltd., China<sup>2)</sup>
- Sartorius Stedim India Pvt. Ltd., India<sup>2)</sup>
- Sartorius Stedim Malaysia Sdn. Bhd., Malaysia<sup>2)</sup>
- Sartorius Mechatronics Hong Kong Ltd., China<sup>2)</sup>
- Sartorius Stedim Australia Pty. Ltd., Australia<sup>2)</sup>

On the Verwaltungsrat (Administrative Board) of:

- Wave Biotech AG, Switzerland; Chairman<sup>2)</sup>

#### Dr. rer. nat. Günther Maaz

On the Board of Directors of:

- Sartorius North America, Inc., USA, until July 24, 2009<sup>2)</sup>
- Sartorius Mechatronics Corporation, USA, until July 24, 2009
- Sartorius TCC Company, USA, until July 24, 2009
- Denver Instrument, Inc., USA, until July 24, 2009<sup>2)</sup>
- Sartorius Stedim Filters, Inc., Puerto Rico, until July 24, 2009<sup>2)</sup>
- Sartorius Mechatronics Hong Kong Ltd., China, until July 24, 2009<sup>2)</sup>
- Beijing Sartorius Instrument & System Engineering Co. Ltd., China; Vice Chairman until July 24, 2009<sup>2)</sup>
- Sartorius Scientific Instruments (Beijing) Co. Ltd., China, Vice Chairman until July 24, 2009<sup>2)</sup>
- Sartorius Mechatronics India Pvt. Ltd., India, July 24, 2009<sup>2)</sup>

#### Positions Held by the Supervisory Board<sup>1)</sup> as of December 31, 2009

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

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

- Sartorius Stedim Biotech S.A., France<sup>2)</sup>

On the Supervisory Board of:

- Sartorius Stedim Biotech GmbH, Germany; Chairman<sup>2)</sup>
- Takt AG, Germany<sup>3)</sup>
- Wissenschaftliches Institut für Infrastruktur und Kommunikationsdienste GmbH (Scientific Institute for Communication Services) and WIK-Consult GmbH, Germany<sup>3)</sup>

##### Gerd-Uwe Boguslawski

On the Supervisory Board of:

- Demag Cranes & Components GmbH, Germany<sup>3)</sup>
- Demag Cranes AG, Germany<sup>3)</sup>

##### Dr. Dirk Basting

None

##### Annette Becker

None

##### Christiane Benner

On the Supervisory Board of:

- T-Systems International GmbH, Germany<sup>3)</sup>

Uwe Bretthauer  
None

Michael Dohrmann  
None

Dr. Lothar Kappich  
None

##### Prof. Dr. Gerd Krieger

On the Supervisory Board of:

- ARAG Lebensversicherungs-AG, Germany<sup>3)</sup>
- ARAG Krankenversicherungs-AG, Germany<sup>3)</sup>

##### 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<sup>3)</sup>

##### Dr. Michael Schulenburg

On the Administrative Board of:

- TÜV Rheinland, Berlin, Brandenburg, Pfalz e.V., Cologne<sup>3)</sup>

On the Beirat (Advisory Board) of:

- Lohmann GmbH & Co. KG, Germany<sup>3)</sup>

On the Board of Directors of:

- Cognis Holding Luxembourg S.à.r.l., Luxembourg<sup>3)</sup>

On the Supervisory Board of:

- Cognis GmbH, Germany; Chairman<sup>3)</sup>

##### Manfred Werner

On the Consejo de Administración (Board of Directors) of:

- Sartorius Mechatronics Spain S.A., Spain; Presidente (President)<sup>2)</sup>
- Sartorius Stedim Spain S.A., Spain; Presidente (President)<sup>2)</sup>

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

- Sartorius Mechatronics Belgium N.V., Belgium<sup>2)</sup>
- Sartorius Stedim Belgium N.V., Belgium<sup>2)</sup>

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

<sup>2)</sup> Positions held within the Group

<sup>3)</sup> External positions held by the Supervisory Board or Executive Board members as at December 31, 2009

# About This Publication

**Published by**

Sartorius AG  
Group Corporate Communications  
37070 Goettingen, Germany

**Editorial Deadline**

March 8, 2010

**Published in**

March 2010

**Photography**

Peter Ginter | Lohmar

Original German version  
published on March 9, 2010.

Sartorius AG  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289

[www.sartorius.com](http://www.sartorius.com)