



sartorius

Sartorius AG
2012 Financial Statements

2012

Contents

01 Financial Statements and Notes

- 4 Balance Sheet
- 5 Income Statement
- 6 Notes to the Individual Balance Sheet Items
- 10 Notes to the Income Statement
- 13 Remuneration Report
- 19 Declaration of the Executive Board
- 20 Independent Auditors' Report

02 Supplementary Information

- 22 Major Affiliates and Ownership in 2012
- 24 Executive Board and Supervisory Board | Positions Held
- 28 About This Publication

Forward-looking Statements Contain Risks

This annual report contains statements concerning the future performance of Sartorius AG. 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 origi-

nal 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.

Financial Statements and Notes

01

Balance Sheet

Assets	Notes		Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
A. Assets				
I. Property, plant and equipment	(4)		5,659	5,551
II. Financial assets	(5)		468,288	466,824
			473,947	472,375
B. Current assets				
I. Trade and other receivables	(6)		51,202	4,129
II. Cash on hand, deposits in banks			471	65
			51,673	4,194
C. Prepaid expenses	(7)		1,532	1,452
			527,152	478,021

Equity and Liabilities	Notes	€ in K	Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
A. Equity				
I. Subscribed capital	(8)	18,720		18,720
Nominal value of treasury shares		-1,673		-1,673
Issued capital			17,047	17,047
II. Capital reserves			101,397	101,397
III. Earnings reserves	(9)		10,867	10,867
IV. Retained profit incl. net profit	(9)		158,992	170,058
			288,303	299,369
B. Provisions	(10)		22,490	24,377
C. Liabilities	(11)		216,359	154,025
D. Deferred taxes	(26)		0	250
			527,152	478,021

Income Statement

	Notes	2012 € in K	€ in K	2011 € in K
1. Sales revenue	(15)	6,667		4,135
2. Other operating income	(16)	476		362
			7,143	4,497
3. Employee benefits expense	(17)	3,948		2,709
4. Depreciation and amortization expense	(18)	235		224
5. Other operating expenses	(19)	4,592		3,232
6. Income from investments	(20)	9,770		8,793
7. Profit reported due to a profit and loss transfer agreement	(21)	305		772
			1,301	3,400
8. Earnings before interest and taxes			8,444	7,897
9. Interest and similar income	(22)	1		0
10. Interest and similar expenses	(22)	5,952		5,828
			-5,951	-5,828
11. Profit before tax			2,493	2,069
12. Extraordinary expenses Extraordinary income	(23)		0	155,787
13. Income tax expense Income from income tax	(25) (26)	-250		4,380
14. Other taxes		1		2
			-249	4,382
15. Net profit			2,742	153,474
16. Profit brought forward			156,250	16,584
17. Retained profit			158,992	170,058

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 accounting and valuation methods of the previous reporting year were retained.

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. 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, 2012	7,481	0	239	251	7,971
Investments	268	0	4	70	342
Disposals	0	0	43	0	43
Transfers	251	0	0	-251	0
Gross book values as of Dec. 31, 2012	8,000	0	200	70	8,270
Depreciation as of Jan. 1, 2012	2,198	0	222	0	2,420
Depreciation in 2012	228	0	7	0	235
Disposals	0	0	43	0	43
Transfers	0	0	0		
Depreciation as of Dec. 31, 2012	2,426	0	186	0	2,612
Net book values as of Dec. 31, 2012	5,574	0	14	70	5,658
Net book values as of Dec. 31, 2011	5,283	0	17	251	5,551

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, provided that these are compatible with the German commercial principles of proper accounting. Where diminution in value is anticipated over the long term, unscheduled impairment losses are recorded.

3. Currency Translation

Liabilities in foreign currencies whose remaining term does not exceed one year are measured at the average spot rate valid on the reporting date. All other liabilities in foreign currencies are measured at their conversion rate upon the invoice date or the higher average spot rate on the reporting date.

Receivables in foreign currencies whose remaining term does not exceed one year are measured at the average spot rate valid on the reporting date. All other receivables in foreign currencies are measured at their conversion rate upon the invoice date or the higher average spot rate on the reporting date.

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

Buildings	15 to 50 years
Office equipment	3 to 13 years

5. 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, 2012	467,581	397	467,978
Investments	1,507	0	1,507
Disposals	43	0	43
Transfers	0	0	0
Gross book values as of Dec. 31, 2012	469,045	397	469,442
Impairment losses as of Jan. 1, 2012	757	397	1,154
Impairment losses in 2012	0	0	0
Disposals	0	0	0
Additions	0	0	0
Impairment losses as of Dec. 31, 2012	757	397	1,154
Net book values as of Dec. 31, 2012	468,288	0	468,288
Net book values as of Dec. 31, 2011	466,824	0	466,824

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.

Please refer to "Major Affiliates and Ownership."

6. Trade and Other Receivables

	Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
Receivables from affiliated companies	49,532	2,142
Other assets	1,670	1,987
	51,202	4,129
Of which due in more than one year:		
Other assets	1,144	867

Trade and other receivables were reported so that all discernible risks are covered. General credit risks are sufficiently covered by lump-sum valuation allowances. Receivables arising from reinsurance policies that are not exempt from attachment by all creditors are recognized in the actuarial reserves according to the company's business plan.

7. Prepaid Expenses

	Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
Loan discounts	1,532	1,451
Other prepaid expenses	0	1
	1,532	1,452

Loan discounts are released over the term of the respective finance loans.

8. 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.

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 (8.9%) of the capital stock. In fiscal 2012, no treasury shares were purchased.

The calculated par value of the treasury shares of €1,672,927.00 was deducted from the capital stock according to Section 272, Subsection 1a, of the German Commercial Code (HGB).

9. Earnings Reserves and Retained Profits

The earnings reserves remained unchanged at €10,868 K.

In fiscal 2012, the earnings reserves developed as follows:		€ in K
At Jan. 1, 2012		170,058
– Dividends paid in 2012		–13,808
= Profit carried forward		156,250
+ Retained profit incl. net profit for 2012		2,742
At Dec. 31, 2012		158,992

10. Provisions

	Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
Provisions for retirement benefits and similar obligations	16,085	15,725
Tax provisions	71	4,248
Other provisions	6,334	4,404
	22,490	24,377

Provisions for retirement benefits are set up according to actuarial principles in line with the valuation assumptions and methods pursuant to Section 249, Subsection 1, of the German Commercial Code (HGB) in conjunction with Section 253, Subsection 1, of HGB.

	Beginning of the reporting year	End of the reporting year
Accounting standards applied	HGB amended by BilMoG*	
Assessment standards applied	Projected Unit Credit Method (PUCM)	
Assumptions on mortality and invalidity	Ref. interest rate table "2005G" from Klaus Heubeck	
Annual discount rate acc. to RückAbzinsV** (Simplification rule: interest rate for remaining period of 15 years)	5.16% (as of Dec. 31, 2011)	5.13% (as of Dec. 31, 2012)
Expected annual increase in old-age pensions	2.00%	2.00%
Expected annual increase in income		
– Indiv. agreed pension plans	3.00%	3.00%
– Gen. in-house retirem. plan	–	–
Annual increase in fixed amounts considered	–	–
Fluctuation	–	–
Calculated retirement age	Preretirement age	

Insolvency-protected investments in reinsurance policies to hedge obligations for early retirement benefits that were reported at a present value of €897 K (2011: €843 K) and correspond to the acquisition costs were offset against the present value of the respective retirement benefits of €1,269 K (2011: €1,132 K).

Provisions for employee retirement benefits and similar obligations developed as follows during fiscal 2012:

	€ in K
Gross provisions as of January 1, 2012	16,567,851
Interest expense	823,204
Effect of the change in the actuarial interest rate	123,867
Other change	–533,880
Gross provisions as of December 31, 2012	16,981,042
Reinsurance reserves for hedging obligations acc. to Sec. 246, Sub. 2, of HGB, as of Dec. 31, 2012 (present value of pledged reinsurance policies)	896,685
Net provisions as of December 31, 2012	16,084,357

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 "Versorgungswerk" 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 €8.4 million of the provisions set up on December 31, 2012, is allocated to Sartorius AG's general in-house retirement plan, and €8.6 million to cover individually agreed pension plans.

The other provisions are recognized at the settlement amount for long-term provisions as are dictated by prudent business judgment. In assessing this settlement amount, all discernible risks from incomplete contracts and uncertain liabilities are considered. The other provisions with a term of less than one year are not discounted.

The other provisions include the following amounts:

	Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
Employee benefits expense	3,308	2,739
Accounts outstanding	650	100
Anticipated losses related to incomplete contracts	1,861	861
Other	515	704
	6,334	4,404

* BilMoG = German Accounting Law Modernization Act
 ** RückAbzinsV = German Regulation on the Discounting of Provisions

11. Liabilities

This item consists of the following:	Disclosed on balance sheet Dec. 31, 2012 € in K	Remaining term of more than five years € in K	Remaining term of up to one year € in K	Disclosed on balance sheet Dec. 31, 2011 € in K	Remaining term of more than five years € in K	Remaining term of up to one year € in K
Liabilities to banks	149,500	50,500	13,500	63,000	0	15,000
Trade payables	156	0	156	885	0	885
Payables to companies in which investments are held	65,238	0	65,238	88,872	0	88,872
Other liabilities	1,465	0	1,465	1,268	0	1,268
	216,359	50,500	80,359	154,025	0	106,025

The other liabilities consist of the following:	Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
Taxes	50	115
Liabilities relating to social security	0	0

Liabilities are reported as settlement amounts (amounts repayable).

12. Contingent Liabilities

	Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
Guarantees (suretyships) and similar contingent liabilities	15,085	15,981

Guarantees (suretyships) exist in the amount of €15,085 K granted to affiliated companies that are included in the consolidated financial statements.

The guarantees and/or suretyships in the amount of €15,085 K (2011: €15,981 K) given by Sartorius AG to affiliated companies for their local credit lines at banks are not to be recognized because there are no signs that these subsidiaries will fail to meet their financial obligations to banks on time.

13. Disclosures on Transactions Not Included in the Balance Sheet

In addition to provisions, liabilities and contingent liabilities, other financial obligations consist of the following:

	Dec. 31, 2012 € in K	Dec. 31, 2011 € in K
Rental and leasing contracts		
– due in fiscal 2012	50	
(previous year: due in fiscal 2011)		39
– due in any one fiscal year from 2013 to 2016	38	
(previous year: due in any one fiscal year from 2012 to 2015)		25
	88	64

14. Derivative Financial Instruments

To hedge the company's term loan of €55 million of its syndicated credit facility agreement – which was refinanced in April 2011 – against interest rate risk, rate swaps (interest rate hedges) have been concluded with a nominal volume of €45 million and for a term until September 2013. Moreover, we entered into forward interest swaps with a nominal volume of €30 million and for a term running from 2013 to 2016 in some instances. For these interest swaps, a reserve for anticipated losses related to incomplete contracts of €1,861 K (2011: €861 K) had to be set up by the reporting date.

Notes to the Income Statement

15. Sales Revenue

	2012 € in K	2011 € in K
Sales revenue, which is broken down by geographical market, was as follows:		
Germany	6,667	4,135
All other countries	0	0
Total	6,667	4,135

16. Other Operating Income

	2012 € in K	2011 € in K
Income from the release of provisions	150	102
Other income	326	260
Total	476	362

17. Employee Benefits Expense

	2012 € in K	2011 € in K
Wages and salaries	3,360	2,695
Social security	0	13
Retirement benefits expense	588	0
Total	3,948	2,708

In the reporting year, Sartorius AG as a holding company did not employ any staff. The employee benefits expense disclosed for fiscal 2012 refers to benefits for the Executive Board of Sartorius AG.

18. Depreciation Expense

	2012 € in K	2011 € in K
Depreciation of property, plant and equipment		
Scheduled	235	224

19. Other Operating Expenses

	2012 € in K	2011 € in K
Legal and consulting expenses	1,745	1,184
Costs for outside services	972	240
Supervisory Board remuneration	753	688
Insurance expenses	192	3
Ancillary staff costs	187	194
Travel, entertainment and representation costs	180	239
Fees and contributions	120	118
Car and leasing expenses	109	89
Maintenance costs	99	81
Advertising expenses	68	40
Office, mail and other administrative expenses	32	27
Other finance charges	16	8
Other expenses	120	321
Total	4,593	3,232

20. Income from Investments

	2012 € in K	2011 € in K
Income from investments	9,770	8,793

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

21. Profit Reported Due to a Profit and Loss Transfer Agreement

In fiscal 2011, Sartorius AG concluded a profit and loss transfer agreement with Sartorius Corporate Administration GmbH. The amount of €305 K reported for this contract in the year under review (2011: €772 K) is reflected in the company's earnings.

22. Interest

	2012 € in K	2011 € in K
Interest and similar income	1	0
– of which from affiliated companies	(0)	(0)
Interest and similar expenses	5,952	5,828
– of which from affiliated companies	(1,442)	(798)
	-5,951	-5,828

Interest expenses for fiscal 2012 include expenses of €868 K (2011: €891 K) that resulted from compounding provisions for employee retirement benefits and similar obligations.

23. Extraordinary Income | Extraordinary Expenses

In the extraordinary income posted for fiscal 2011, the company reports the result of disclosing hidden reserves from the spin-off of the operational business unit "static (lab) weighing technology" into Sartorius Weighing Technology GmbH.

24. 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 the previous year, under employee benefits expense.

For Sartorius AG, the income to be allocated to the other business years is €150 K (2011: €111 K). This concerns income from the release of provisions. Sartorius AG did not incur any material non-periodic expenses in fiscal 2012 (2011: €0 K).

25. Income Tax Expense

	2012 € in K	2011 € in K
Current income taxes	0	4,131
Deferred taxes	-250	250
Income taxes for the previous years	0	0
	-250	4,381

Extraordinary income accounts for the total amount of current income taxes reported.

26. Deferred Taxes

The temporary concept is used to calculate deferred taxes. Such deferred taxes are assessed for temporary differences between commercial accounting and tax accounting of assets, liabilities and prepaid expenses in Germany. This valuation not only includes the differences arising from Sartorius AG's own balance sheet items, but also those that exist for subsidiaries or business partnerships or other non-corporate entities in which Sartorius AG as a share owner holds an investment. Deferred taxes are assessed on the basis of the combined income tax rate for Sartorius AG, which is currently 30.0%. The combined income tax rate covers corporate income tax, industrial and commercial profits tax and the solidarity surcharge on income tax for reconstruction of eastern Germany. However, unlike in the former case, deferred taxes from temporary accounting differences for investments that have the legal form of a business partnership or a non-corporate entity are measured on the basis of a combined income tax rate, which covers only corporate income tax and the solidarity surcharge on income tax; this particular income tax rate is currently 15.83%. According to the option provided by Section 274, Subsection 1, sentence 2, of the German Commercial Code (HGB), deferred taxes are reported as amounts netted out. If an overall tax burden were to be yielded by these calculations, this would be disclosed as deferred tax liabilities. Conversely, if these calculations were to yield a tax benefit, the company would not exercise its right to elect to report this as a deferred tax asset. In the reporting year, assessment of the temporary differences yielded a deferred tax asset, which was not recognized.

Deferred tax assets result especially from the differences in the items disclosed for "Provisions for retirement benefits" and the "Other provisions." Furthermore, deferred tax assets are yielded for tax loss carry-forwards according to Section 15a of the German Income Tax Law (EStG), which may be used within the next five years. Deferred tax liabilities offsetting these deferred tax assets result primarily for the financial assets and concern accounting changes for business partnerships as well as different recognition amounts as a consequence of the spin-off in 2011.

Moreover, Sartorius AG has tax loss carry-forwards for corporate and commercial income tax as well as unused interest carry-forwards. As in the previous year, deferred tax assets are not considered for these carry-forwards in the reporting year because, from today's stance, use is not sufficiently probable for the remaining amounts carried forward.

27. Other Disclosures

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 7, 2012, 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:

	2012 € in K	2011 € in K
Auditing of the annual financial statements and of the consolidated annual financial statements	65	70
Other independent reporting services	93	341
Tax consultation incl. preparation of tax returns	99	195
Other services	90	14
	347	620

Disclosure acc. to Sec. 285, No. 21, of HGB

The following table covers all business with related companies and persons for fiscal 2012:

All amounts in € in K (thousand) for fiscal 2012	Sales transactions	Purchase transactions	Financing received	Financing provided	Other transactions
Affiliated companies	0	0	18,651	48,829	4,561
Companies in which investments are held	0	0	0	0	0
Persons in key positions	0	0	0	0	0

All amounts in € in K (thousand) for fiscal 2012	Sales transactions	Purchase transactions	Financing received	Financing provided	Other transactions
Affiliated companies	0	0	42,830	0	11,739
Companies in which investments are held	0	0	0	0	0
Persons in key positions	0	0	0	0	0

For information on the remuneration of the Supervisory Board members and of the Executive Board members, please refer to the following section.

Remuneration Report

Remuneration of the Supervisory Board Members

€ in K	2012	2011
Total remuneration	860	773
Fixed remuneration	300	290
Performance-related remuneration	283	217
Compensation for committee work	60	58
Meeting attendance fee	114	123
Remuneration from Sartorius Weighing Technology GmbH, Goettingen	20	2
Total remuneration for the Sartorius Stedim Biotech subgroup	83	83
Remuneration from Sartorius Stedim Biotech GmbH, Goettingen	38	38
Remuneration from Sartorius Stedim Biotech S.A., Aubagne	45	45

Prof. Dr. Dres. h.c. Arnold Picot (Chairman)

Total remuneration	255	230
Fixed remuneration	60	60
Performance-related remuneration	56	45
Compensation for committee work	18	20
Meeting attendance fee	18	20
Remuneration from Sartorius Weighing Technology GmbH, Goettingen	20	2
Total remuneration for the Sartorius Stedim Biotech subgroup	83	83
Remuneration from Sartorius Stedim Biotech GmbH, Goettingen	38	38
Remuneration from Sartorius Stedim Biotech S.A., Aubagne	45	45

Gerd-Uwe Boguslawski¹⁾ (Vice Chairman)

Total remuneration	103	102
Fixed remuneration	40	40
Performance-related remuneration	37	30
Compensation for committee work	12	12
Meeting attendance fee	14	20

Dr. Dirk Basting

Total remuneration	45	41
Fixed remuneration	20	20
Performance-related remuneration	19	15
Meeting attendance fee	6	6

Annette Becker¹⁾

Total remuneration	45	41
Fixed remuneration	20	20
Performance-related remuneration	19	15
Meeting attendance fee	6	6

€ in K	2012	2011
Christiane Benner¹⁾ (until April 19, 2012)		
Total remuneration	13	41
Fixed remuneration	6	20
Performance-related remuneration	6	15
Meeting attendance fee	1	6

Uwe Bretthauer¹⁾

Total remuneration	69	67
Fixed remuneration	20	20
Performance-related remuneration	19	15
Compensation for committee work	12	12
Meeting attendance fee	18	20

Michael Dohrmann¹⁾

Total remuneration	45	40
Fixed remuneration	20	20
Performance-related remuneration	19	15
Meeting attendance fee	6	5

Dr. Lothar Kappich

Total remuneration	45	48
Fixed remuneration	20	20
Performance-related remuneration	19	15
Compensation for committee work	0	3
Meeting attendance fee	6	10

Petra Kirchoff

Total remuneration	45	41
Fixed remuneration	20	20
Performance-related remuneration	19	15
Meeting attendance fee	6	6

Karoline Kleinschmidt¹⁾ (as of April 19, 2012)

Total remuneration	32	0
Fixed remuneration	14	0
Performance-related remuneration	13	0
Meeting attendance fee	5	0

Prof. Dr. Gerd Krieger

Total remuneration	56	47
Fixed remuneration	20	20
Performance-related remuneration	19	15
Compensation for committee work	6	3
Meeting attendance fee	11	9

¹⁾ 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 "Deutscher Gewerkschaftsbund."

€ in K	2012	2011
Prof. Dr. rer. nat. Dr.-Ing. Heribert Offermanns (until April 19, 2012)		
Total remuneration	13	41
Fixed remuneration	6	20
Performance-related remuneration	6	15
Meeting attendance fee	1	6

Prof. Dr. Thomas Scheper (as of April 19, 2012)		
Total remuneration	32	0
Fixed remuneration	14	0
Performance-related remuneration	13	0
Meeting attendance fee	5	0

Dr. Michael Schulenburg (until May 20, 2012)		
Total remuneration	0	29
Fixed remuneration	0	8
Performance-related remuneration	0	6
Compensation for committee work	0	7
Meeting attendance fee	0	8

Remuneration of the Executive Board Members

€ in K	2012	2011
Total remuneration		
Cash benefits for the fiscal year	2,679	2,458
Fixed remuneration for the fiscal year	1,246	1,186
Annually paid variable remuneration	558	659
Variable remuneration with multi-year components paid (fair value)	875	613
Change in the fair values of the variable multi-year components of remuneration from previous years	611	223
Pension commitments		
Additions to provisions for retirement benefits	283	120

Dr. Joachim Kreuzburg		
Cash benefits for the fiscal year	1,302	1,182
Fixed remuneration for the fiscal year	603	568
Annually paid variable remuneration	272	318
Variable remuneration with multi-year components paid (fair value)	427	296
Change in the fair values of the variable multi-year components of remuneration from previous years	366	156
Pension commitments		
Additions to provisions for retirement benefits	172	120

€ in K	2012	2011
Prof. Dr. Klaus Rüdiger Trützschler (as of November 29, 2011)		
Total remuneration	62	5
Fixed remuneration	20	2
Performance-related remuneration	19	1
Compensation for committee work	12	1
Meeting attendance fee	11	1

Beyond their Supervisory Board remuneration, the employee representatives who are employees within the Sartorius Group receive compensation that is not related to their service on the Supervisory Board.

€ in K	2012	2011
Jörg Pfirrmann		
Cash benefits for the fiscal year	575	541
Fixed remuneration for the fiscal year	269	262
Annually paid variable remuneration	119	145
Variable remuneration with multi-year components paid (fair value)	187	135
Change in the fair values of the variable multi-year components of remuneration from previous years	103	27
Pension commitments		
Additions to provisions for retirement benefits	111	0

Reinhard Vogt		
Cash benefits for the fiscal year	801	735
Fixed remuneration for the fiscal year	374	356
Annually paid variable remuneration	166	197
Variable remuneration with multi-year components paid (fair value)	261	182
Change in the fair values of the variable multi-year components of remuneration from previous years	142	39

Variable Remuneration Determined Using Multi-year Assessment

€ in K	Fair value when granted as of Jan. 1 of the particular year	Fair value at year end on Dec. 31, 2011	Fair value at year end on Dec. 31, 2012	Change in value in fiscal 2012
Dr. Joachim Kreuzburg				
Granted for fiscal 2010	138	149	153	4
Granted for fiscal 2011	138	140	142	2
Granted for fiscal 2012	146	0	148	2
Total	421	289	444	9
Jörg Pfirrmann				
Granted for fiscal 2010	53	57	59	2
Granted for fiscal 2011	63	64	64	1
Granted for fiscal 2012	64	0	65	1
Total	179	121	189	4
Reinhard Vogt				
Granted for fiscal 2010	75	81	84	2
Granted for fiscal 2011	85	86	88	1
Granted for fiscal 2012	90	0	91	1
Total	250	168	262	5

Components with a Long-term Incentive Effect

	Number of phantom stock units	Price on assignment* in €	Fair value when granted on Jan. 1 of the particular year in € in K	Fair value at year-end on Dec. 31, 2011 in € in K	Fair value at year-end on Dec. 31, 2012 in € in K	Paid in fiscal 2012 in € in K	Change in value in fiscal 2012 in € in K	Status
Dr. Joachim Kreuzburg								
Tranche for 2006	3,871	21.31	83	128	0	128	0	Paid in 2012
Tranche for 2007	3,593	34.79	125	119	249	0	130	Exercisable
Tranche for 2009	16,851	8.16	138	344	0	344	0	Paid in 2012
Tranche for 2010	8,715	15.78	138	271	334	0	63	Not exercisable
Tranche for 2011	5,165	26.62	138	156	323	0	167	Not exercisable
Sum of the tranches from the previous years	38,195		622	1,018	906	472	360	
Tranche for 2012	4,416	33.12	146	0	279	0	133	Not exercisable
Total sum of tranches	42,611		768	1,018	1,185	472	493	
Jörg Pfirrmann								
Tranche for 2009	2,703	8.16	22	55	0	55	0	Paid in 2012
Tranche for 2010	3,334	15.78	53	105	128	0	23	Not exercisable
Tranche for 2011	2,348	26.62	63	71	147	0	76	Not exercisable
Sum of the tranches from the previous years	8,385		138	231	275	55	99	
Tranche for 2012	1,937	33.12	64	0	122	0	58	Not exercisable
Total sum of tranches	10,322		202	231	397	55	157	
Reinhard Vogt								
Tranche for 2009	4,054	8.16	33	83	0	83	0	Paid in 2012
Tranche for 2010	4,754	15.78	75	148	182	0	34	Not exercisable
Tranche for 2011	3,193	26.62	85	96	200	0	104	Not exercisable
Sum of the tranches from the previous years	12,001		193	327	382	83	138	
Tranche for 2012	2,699	33.12	89	0	170	0	81	Not exercisable
Total sum of tranches	14,700		282	327	552	83	219	

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 plan replaces the former remuneration components that had been paid out annually until then. 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

minimum 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, which is included in the Group Management Report. Please refer to this information in the 2012 Annual Report for the Sartorius Group.

* Average share price over the last 20 days of trading in the preceding fiscal year

Remuneration of Former Managing Directors

	2012 € in K	2011 € in K
Remuneration of former managing directors and members of the Executive Board and their surviving dependents	386	390
Retirement benefits and pension obligations to former managing directors and members of the Executive Board as well as their surviving dependents	7,045	6,692

Number of Employees

No staff members were employed in the reporting year.

Proposal for Appropriation of Profits

The Executive Board will submit a proposal to the Annual Shareholders' Meeting to appropriate the retained profit of €158,991,646.29 reported by Sartorius AG for the year ended December 31, 2012, as follows:

	€
Payment of a dividend of €0.94 per ordinary share	8,016,372.64
Payment of a dividend of €0.96 per preference share	8,178,256.32
Unappropriated profit carried forward	142,797,017.33
	158,991,646.29

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 as of December 31, 2012, according to Section 21, Subsection 1 or 1a, of the German Securities Trading Act (WpHG):

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	30%	Apr. 6, 2011		30.01%	All acc. to Subsec. 1, No. 1

Goettingen, February 18, 2013

Sartorius Aktiengesellschaft
The Executive Board

Declaration of the Executive Board

We declare to the best of our knowledge that the year-end financial statements for fiscal 2012 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 its 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 18, 2013

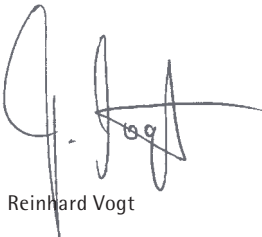
Sartorius Aktiengesellschaft
The Executive Board



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, Goettingen, Germany, and the management report integrated into the Group Management Report for the fiscal year from January 1 to December 31, 2012. Accounting and preparation of the annual financial statements and of the management report according to the German commercial accounting standards are the responsibility of the Executive Board of Sartorius AG. Our responsibility is to express an opinion on the company's annual financial statements, including its accounting records, and on its management report, based on our audit.

We conducted our annual 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 year-end financial statements and provides an overall true and fair view of the company's situation, and accurately presents the opportunities and risks of its future development.

Hanover, February 18, 2013

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



Henning Scharpenberg
Auditor



Dr. Meyer
Auditor

Supplementary Information

02

Major Affiliates and Ownership in 2012

	Sartorius AG's ownership in %	Equity at Dec. 31, 2012 € in K	Net profit at Dec. 31, 2012 € in K
Europe			
Sartorius Stedim Biotech S.A., Aubagne, France along with its subsidiaries:	74.4	117,742	26,197
Sartorius Stedim Belgium N.V., Vilvoorde, Belgium*)	100.0	1,065	407
Distribo GmbH, Goettingen, Germany***)	26.0	535	84
Sartorius Stedim Nordic A/S, Taastrup, Denmark*)	100.0	1,192	985
Sartorius Stedim Biotech GmbH, Goettingen, Germany*)	100.0	93,906	34,278
Sartorius Stedim Systems GmbH, Guxhagen, Germany*)	100.0	9,870	0
Sartorius Stedim Plastics GmbH, Goettingen, Germany*)	100.0	3,204	1,270
Sartorius Stedim UK Ltd., Epsom, U.K.*)	100.0	2,184	732
Sartorius Stedim Lab Ltd., Louth, U.K.*)	100.0	1,950	1,340
Sartorius Stedim France S.A.S., Aubagne, France*)	100.0	7,241	2,007
Sartorius Stedim Aseptics S.A., Lourdes, France*)	100.0	5,725	2,352
Sartorius Stedim Ireland Ltd., Dublin, Ireland*)	100.0	401	295
Sartorius Stedim Italy S.p.A., Florence, Italy*)	100.0	4,836	1,087
Sartorius Stedim Netherlands B.V., Nieuwegein, Netherlands*)	100.0	1,541	495
Sartorius Stedim Austria GmbH, Vienna, Austria*)	100.0	1,123	770
Sartorius Stedim Poland Sp. z.o.o., Kostrzyn, Poland*) ***)	100.0	108	-393
OOO Sartorius ICR, St. Petersburg, Russia*) **)	100.0	121	125
Sartorius Stedim Switzerland AG, Tagelswangen, Switzerland*)	100.0	5,850	1,888
Sartorius Stedim Spain S.A., Madrid, Spain*)	100.0	258	140
Sartorius Stedim Hungaria Kft., Budapest, Hungary*) ***)	100.0	294	9
North America			
Sartorius Stedim Filters Inc., Yauco, Puerto Rico*)	100.0	21,580	1,703
Sartorius Stedim North America Inc., Bohemia, New York, USA*)	100.0	59,340	17,101
Asia Pacific			
Sartorius Stedim Australia Pty. Ltd., Dandenong South, Victoria, Australia*)	100.0	1,351	722
Sartorius Stedim Biotech (Beijing) Co. Ltd., Beijing, China*)	100.0	2,075	231
Sartorius Stedim India Pvt. Ltd., Bangalore, India*)	100.0	3,013	241
Sartorius Stedim Japan K.K., Tokyo, Japan*)	100.0	3,397	1,071
Sartorius Korea Biotech Co. Ltd., Seoul, South Korea*)	49.0	4,398	2,399
Sartorius Stedim Malaysia Sdn. Bhd., Kuala Lumpur, Malaysia*)	100.0	833	759
Sartorius Stedim Singapore Pte. Ltd., Singapore*)	100.0	1,639	943
Other Remaining Markets			
Sartorius Stedim SUS S.A.R.L., M'Hamdia, Tunisia*)	100.0	4,539	1,041

The companies identified by an *) are indirectly held by Sartorius Stedim Biotech S.A., France.

The equity investments in the companies that are identified by **) are indirectly held.

As the financial statements of the subsidiaries identified by ***) were not available at the time our consolidated financial statements were prepared, the information from the annual financial statements from 2011 was considered.

	Sartorius AG's ownership in %	Equity at Dec. 31, 2012 € in K	Net profit at Dec. 31, 2012 € in K
Europe			
Sartorius Belgium N.V., Vilvoorde, Belgium**)	100.0	-579	-21
Sartorius Nordic A/S, Taastrup, Denmark**)	100.0	128	-164
Sartorius Weighing Technology GmbH, Goettingen, Germany**)	100.0	43,826	-3,788
Denver Instrument GmbH, Goettingen, Germany**)	100.0	1,794	1,257
Sartorius Mechatronics C&D GmbH & Co. KG, Aachen, Germany, including Sartorius-Verwaltungs-GmbH	100.0	-491	1,239
Sartorius Mechatronics T&H Hamburg GmbH, Hamburg, Germany	100.0	24,080	2,954
Sartorius Corporate Administration GmbH, Goettingen, Germany	100.0	639	0
SI Weende-Verwaltungs-GmbH, Goettingen, Germany	100.0	23	-2
SIV Weende GmbH & Co. KG, Goettingen, Germany	100.0	2,962	232
SI Grone 1-Verwaltungs-GmbH, Goettingen, Germany**)	100.0	24	-1
SIV Grone 1 GmbH & Co. KG, Goettingen, Germany**)	100.0	5,327	-124
SI Grone 2-Verwaltungs-GmbH, Goettingen, Germany	100.0	25	0
SIV Grone 2 GmbH & Co. KG, Goettingen, Germany	100.0	10	0
Sartorius Lab Holding GmbH, Goettingen, Germany	100.0	149,702	120
Sartorius UK Ltd., Epsom, U.K.**)	100.0	729	-1,410
Biohit Ltd., Torquay, U.K.**)	100.0	249	-169
Sartorius Biohit Liquid Handling Oy, Helsinki, Finland**)	100.0	5,617	-383
Sartorius France S.A.S., Aubagne, France**)	100.0	752	-927
VL Finance S.A.S., Aubagne, France	100.0	54,863	2,268
Biohit S.A.S., Dourdan, France**)	100.0	1,817	187
Sartorius Ireland Ltd., Dublin, Ireland**)	100.0	97	-206
Sartorius Italy S.r.l., Florence, Italy**)	100.0	787	-242
Sartorius Netherlands B.V., Nieuwegein, Netherlands**)	100.0	324	180
Sartorius Austria GmbH, Vienna, Austria**)	100.0	3,191	516
Sartorius Poland Sp. z o.o., Kostrzyn, Poland**)	100.0	339	-158
ZAO Sartogossm, St. Petersburg, Russia**)	51.0	675	193
Biohit OOO, St. Petersburg, Russia**)	100.0	761	362
Sartorius Mechatronics Switzerland AG, Dietikon, Switzerland**)	100.0	-285	-76
Sartorius Spain S.A., Madrid, Spain**)	100.0	-1,143	69
Sartorius Hungaria Kft., Budapest, Hungary**)	100.0	852	204
North America			
Sartorius North America Inc., Bohemia, New York, USA**)	100.0	24,508	0
Sartorius Corporation, Bohemia, New York, USA**)	100.0	-3,399	-9,633
Sartorius TCC Company, Arvada, Colorado, USA**)	100.0	6,955	167
Sartorius Canada Inc., Mississauga, Canada**)	100.0	72	-70
Asia Pacific			
Sartorius Australia Pty. Ltd., Dandenong South, Victoria, Australia**)	100.0	612	247
Denver Instrument (Beijing) Co. Ltd., Beijing, China**)	100.0	2,939	102
Sartorius Scientific Instruments (Beijing) Co. Ltd., Beijing, China**)	100.0	20,202	3,846
Biohit Biotech (Suzhou) Co. Ltd., Shanghai, China**)	100.0	881	-45
Sartorius Hong Kong Ltd., Kowloon, Hong Kong**)	100.0	3,049	41
Sartorius Mechatronics India Pvt. Ltd., Bangalore, India**)	100.0	4,962	10
Sartorius Weighing India Pvt. Ltd., Bangalore, India**)	100.0	2,957	-126
Biohit Biotech Systems (India) Pvt. Ltd., Chennai, India**)	100.0	18	-14
Sartorius Japan K.K., Tokyo, Japan**)	100.0	8,098	1,804
Sartorius Malaysia Sdn. Bhd., Kuala Lumpur, Malaysia**)	100.0	287	165
Sartorius Mechatronics Philippines Inc., Makati City, Philippines**)	100.0	389	41
Sartorius Singapore Pte. Ltd., Singapore**)	100.0	1,118	207
Sartorius Korea Ltd., Seoul, South Korea**)	100.0	4,096	541
Sartorius (Thailand) Co. Ltd., Bangkok, Thailand**)	49.0	672	189
Other Remaining Markets			
Sartorius Argentina S.A., Buenos Aires, Argentina**)	99.0	836	263
Sartorius do Brasil Ltda., São Paulo, Brazil**)	100.0	-1,166	96
Sartorius de México S.A. de C.V., Naucalpan, Mexico**)	99.0	299	513

Executive Board and Supervisory Board

During Fiscal 2012¹⁾

Executive Board

Dr. rer. pol. Joachim Kreuzburg

Dipl.-Ingenieur (Graduate Engineer)
CEO and Chairman
Executive for Labor Relations until July 23, 2012
Strategy, Operations, Legal Issues, Compliance
and Corporate Communications
Born April 22, 1965
Resident of Hanover, Germany
Member since November 11, 2002
Sprecher (Spokesman) from May 1, 2003,
to November 10, 2005
Chairman since November 11, 2005
Executive for Labor Relations from July 24, 2009,
to July 23, 2012
Appointed until November 10, 2015

Jörg Pfirrmann

Dipl.-Ökonom (Graduate Economist)
Executive for Labor Relations since July 24, 2012
Finance, Human Resources, IT and
General Administration
Born November 30, 1972
Resident of Noerten-Hardenberg, Germany
Member since Friday, July 24, 2009
Executive for Labor Relations since July 24, 2012
Appointed until July 23, 2017

Reinhard Vogt

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

Supervisory Board

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

Dipl.-Kaufmann (Graduate in Business,
Administration), university professor
Chairman
Executive Director of the Institute of Information,
Organization and Management,
Faculty of Economics of the
Ludwig Maximilian University in Munich, Germany
Resident of Gauting, Germany

Gerd-Uwe Boguslawski

Dipl.-Sozialwirt (Graduate Social Manager)
Vice Chairman
German Metalworkers' Union (IG Metall) in the southern
Lower Saxony/Harz region in Northeim, Germany
Resident of Hoeckelheim, Germany

Dr. Dirk Basting

Dipl.-Chemiker (Graduate Chemical Engineer)
Resident of Fort Lauderdale,
Florida, USA

Annette Becker

Personalfachkauffrau (HR Specialist)
Chairwoman of the Employees' Council of
Sartorius Corporate Administration GmbH
in Goettingen, Germany
Resident of Uslar, Germany

Uwe Bretthauer

Dipl.-Ingenieur (Graduate Engineer)
Chairman of the Employees' Council of
Sartorius Weighing Technology GmbH,
Chairman of the Group Employees' Council of
Sartorius AG in Goettingen, Germany
Resident of Goettingen, Germany

Michael Dohrmann

Feinmechaniker (Precision Engineer)
Chairman of the Employees' Council
of Sartorius Stedim Biotech GmbH
in Goettingen, Germany
Resident of Reinhausen, Germany

Dr. Lothar Kappich

Dipl.-Ökonom (Graduate Economist)
Managing Director of
ECE Projektmanagement GmbH & Co. KG
in Hamburg, Germany
Resident of Hamburg, Germany

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

Petra Kirchhoff

Dipl.-Volkswirtin (Graduate Political Economist)
 Vice President of Corporate Communications and
 Investor Relations
 Sartorius Corporate Administration GmbH
 in Goettingen, Germany
 Resident of Goettingen, Germany

Karoline Kleinschmidt

Dipl.-Sozialwirtin (Graduate Social Economist)
 Member since April 19, 2012
 Secretary of the German Metalworkers' Union
 (IG Metall) of the district management of Lower Saxony
 and Sachsen-Anhalt in Hanover, Germany
 Resident of Hanover, Germany

Prof. Dr. Gerd Krieger

Lawyer
 Honorary Professor at the
 Heinrich-Heine University in Duesseldorf, Germany
 Resident of Duesseldorf, Germany

Prof. Dr. Thomas Scheper

Dipl.-Chemiker (Graduate Chemical Engineer)
 Member since April 19, 2012
 University professor and head of the Institute of
 Technical Chemistry, Leibniz University
 in Hanover, Germany
 Resident of Hanover, Germany

Prof. Dr. Klaus Rüdiger Trützschler

Dipl.-Wirtschaftsmathematiker (Graduate Business
 Mathematician) and
 Dipl.-Mathematiker (Graduate Mathematician)
 Honorary Professor at the
 Westphalian Wilhelm University in Muenster, Germany
 Resident of Essen, Germany

Christiane Benner

Dipl.-Soziologin (Graduate Sociologist)
 Member until April 19, 2012
 Executive member of the managing committee of
 the German Metalworkers' Union (IG Metall)
 in Frankfurt am Main, Germany
 Resident of Frankfurt am Main, Germany

Prof. Dr. rer. nat. Dr.-Ing.**E.h. Heribert Offermanns**

Dipl.-Chemiker (Graduate Chemical Engineer)
 Member until April 19, 2012
 Honorary Professor at the Johann Wolfgang Goethe
 University of Frankfurt am Main, Germany
 Resident of Hanau, Germany

Committees of the Supervisory Board**Executive Task Committee**

Prof. Dr. Dres. h.c. Arnold Picot
 (Chairman)
 Gerd-Uwe Boguslawski
 Uwe Bretthauer
 Prof. Dr. Gerd Krieger

Audit Committee

Prof. Dr. Klaus Rüdiger Trützschler
 (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
 Prof. Dr. Gerd Krieger

Nomination Committee

Prof. Dr. Gerd Krieger
 Prof. Dr. Dres. h.c. Arnold Picot
 Dr. Lothar Kappich

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

Positions Held by the Members of the Executive Board¹⁾

as of December 31, 2012

Dr. rer. pol. Joachim Kreuzburg

Président-Directeur Général (CEO) of:

- Sartorius Stedim Biotech S.A., France²⁾

On the Supervisory Board of:

- Sartorius Stedim Biotech GmbH, Germany, Vice Chairman²⁾
- Sartorius Weighing Technology GmbH, Germany, Vice Chairman²⁾

On the Board of Directors of:

- Sartorius Stedim North America, Inc., USA²⁾
- Sartorius Stedim Filters, Inc., Puerto Rico²⁾
- Sartorius Japan K.K., Japan²⁾
- Sartorius Stedim Japan K.K., Japan²⁾
- Denver Instrument (Beijing) Co. Ltd., China²⁾
- Sartorius Scientific Instruments (Beijing) Co. Ltd., China²⁾
- Sartorius Hong Kong Ltd., China²⁾
- Sartorius Stedim Lab Ltd., U.K.²⁾
- Sartorius Biohit Liquid Handling Oy, Finland²⁾

On the Regionalbeirat (Regional Advisory Board) of:

- Commerzbank AG, Hamburg, Germany³⁾

On the Beirat (Advisory Board) of:

- Hameln Group GmbH, Germany³⁾

On the Wirtschaftsbeirat (Economic Advisory Board) of:

- Norddeutsche Landesbank, Germany³⁾

Jörg Pfirrmann

On the Verwaltungsrat (Administrative Board) of:

- Sartorius Mechatronics Switzerland AG, Switzerland²⁾

On the Board of Directors of:

- Sartorius North America, Inc., USA²⁾
- Sartorius Corporation, USA²⁾
- Sartorius TCC Company, USA²⁾
- Sartorius Stedim UK Ltd., U.K.²⁾
- Sartorius Stedim Ireland Ltd., Ireland²⁾
- Sartorius Stedim Biotech (Beijing) Co. Ltd., China²⁾

On the Comité Exécutif (Executive Committee) of:

- Sartorius Stedim France S.A.S., France²⁾

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

- Sartorius Stedim Italy S.p.A., Italy²⁾

On the Consejo de Administración

(Board of Directors) of:

- Sartorius Stedim Spain S.A., Spain²⁾

On the Supervisory Board of:

- Sartorius Stedim Nordic A/S, Denmark²⁾
- Sartorius Nordic A/S, Denmark²⁾

On the Unternehmerbeirat

(Employers' Advisory Board) of:

- Gothaer Versicherungsbank WVaG³⁾

Reinhard Vogt

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

- Sartorius Stedim Biotech S.A., France²⁾

On the Board of Directors of:

- Sartorius Stedim North America, Inc., USA²⁾
- Denver Instrument (Beijing) Co. Ltd., China²⁾
- Sartorius Scientific Instruments (Beijing) Co. Ltd., China²⁾
- Sartorius Stedim Biotech (Beijing) Co. Ltd., China²⁾
- Sartorius Stedim India Pvt. Ltd, India²⁾
- Sartorius Stedim Malaysia Sdn. Bhd., Malaysia²⁾
- Sartorius Japan K.K.²⁾
- Sartorius Stedim Japan K.K.²⁾
- Sartorius Hong Kong Ltd., China²⁾
- Sartorius Korea Ltd., South Korea²⁾
- Sartorius Stedim Australia Pty. Ltd., Australia²⁾

On the Verwaltungsrat (Administrative Board) of:

- Sartorius Stedim Switzerland AG, Switzerland, Chairman²⁾

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

²⁾ Positions held within the Group

³⁾ External positions held by members of the Executive Board as of December 31, 2012

Positions Held by the Members of the Supervisory Board¹⁾

as of December 31, 2012

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

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

- Sartorius Stedim Biotech S.A., France²⁾

On the Supervisory Board of:

- Sartorius Stedim Biotech GmbH, Germany, Chairman²⁾
- Sartorius Weighing Technology GmbH, Germany, Chairman²⁾
- Takkt AG, Germany³⁾
- Wissenschaftliches Institut für Infrastruktur und Kommunikationsdienste GmbH (Scientific Institute for Communication Services) and WIK-Consult GmbH, Germany³⁾

Gerd-Uwe Boguslawski

On the Supervisory Board of:

- Demag Cranes & Components GmbH, Germany³⁾
- Demag Cranes AG, Germany³⁾

Dr. Dirk Basting

None

Annette Becker

None

Uwe Bretthauer

None

Michael Dohrmann

None

Dr. Lothar Kappich

None

Petra Kirchhoff

None

Karoline Kleinschmidt

None

Prof. Dr. Gerd Krieger

On the Supervisory Board of:

- ARAG Lebensversicherungs-AG, Germany³⁾
- ARAG Krankenversicherungs-AG, Germany³⁾

Prof. Dr. Thomas Scheper

None

Prof. Dr. Klaus Rüdiger Trützschler

On the Supervisory Board of:

- Takkt AG, Germany, Chairman³⁾
- Wuppermann AG, Germany, Chairman³⁾
- Zwiesel Kristallglas AG, Germany, Chairman³⁾
- Bilfinger Berger AG, Germany³⁾
- Deutsche Bank AG, Germany³⁾

On the Verwaltungsrat (Administrative Board) of:

- Wilh. Werhahn KG, Germany³⁾

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

²⁾ Positions held within the Group

³⁾ External positions held by members of the Supervisory Board as of December 31, 2012

About This Publication

Published by

Sartorius AG
Group Corporate Communications
37070 Goettingen, Germany

Editorial Deadline

March 5, 2013

Published on

March 7, 2013

Original German version
published on March 7, 2013

Sartorius AG
Weender Landstrasse 94-108
37075 Goettingen, Germany

Phone: +49.551.308.0
Fax: +49.551.308.3289

info@sartorius.com
www.sartorius.com

turning science **into solutions**