

SARTORIUS STEDIM BIOTECH S.A. - Combined Annual Shareholders' Meeting – 8 April 2014 - VOTES RESULTS

Number of shares of the issued capital the day of the meeting	17 042 306
Number of shares with voting rights	15 343 596
Number of shares present, represented or with postal voting	13 288 744
Number of voices present, represented or with postal voting	24 910 046
Number of shares present	11 676 508
Number of voices present	23 297 810
Number of shares with postal voting	1 612 236
Number of voices with postal voting	1 612 236
Number of shares with proxy	0
Number of voices with proxy	0
Number of shareholders present, represented or with postal voting	131
Quorum	86,61%

RESOLUTIONS	SHARES PRESENT	% OF THE ISSUED CAPITAL	FOR		AGAINST		ABSTENTION		RESULT
			VOICES	%	VOICES	%	VOICES	%	
1 Financial statements approval	13 288 744	77,98%	24 906 546	99,99%	3 500	0,01%	0	0,00%	adopted
2 Allocation of net income	13 288 744	77,98%	24 889 556	99,92%	20 490	0,08%	0	0,00%	adopted
3 Approval of the regulated agreements	13 288 744	77,98%	24 910 046	100,00%	0	0,00%	0	0,00%	adopted
4 Approval of the 2013 Directors' fees	13 288 744	77,98%	24 897 270	99,95%	12 776	0,05%	0	0,00%	adopted
5 Advisory vote on the elements of compensation due or granted for the 2013 financial year to Mr Joachim Kreuzburg, Chairman of the Board of Directors and Chief Executive Officer	13 288 744	77,98%	23 869 384	95,82%	1 040 662	4,18%	0	0,00%	adopted
6 Advisory vote on the elements of compensation due or granted for the 2013 financial year to Mr Reinhard Vogt, Executive Vice President	13 288 744	77,98%	23 869 384	95,82%	1 040 662	4,18%	0	0,00%	adopted
7 Advisory vote on the elements of compensation due or granted for the 2013 financial year to Mr Volker Niebel, Executive Vice President	13 288 744	77,98%	23 914 878	96,00%	995 168	4,00%	0	0,00%	adopted
8 Advisory vote on the elements of compensation due or granted for the 2013 financial year to Mr Oscar-Werner Reif, Executive Vice President	13 288 744	77,98%	23 869 384	95,82%	1 040 662	4,18%	0	0,00%	adopted
9 Amendment of article 4 of Title I of the Company's bylaws further to the realization of the transfer of activity on management of fluids to the benefit of our subsidiary, Sartorius Stedim FMT	13 288 744	77,98%	24 910 046	100,00%	0	0,00%	0	0,00%	adopted
10 Amendment of article 8.5 of Title III of the Company's bylaws	13 288 744	77,98%	24 910 046	100,00%	0	0,00%	0	0,00%	adopted
11 Amendment of article 11 of Title III of the Company's bylaws	13 288 744	77,98%	24 075 744	96,65%	834 302	3,35%	0	0,00%	adopted
12 Share capital increase through the capitalization of reserves, earnings or premium; determination of the conditions of such operation; authorization granted to the board of directors to carry out this capital increase	13 288 744	77,98%	24 910 046	100,00%	0	0,00%	0	0,00%	adopted
13 Authorization granted to the board of directors to reduce the share capital through the cancellation of shares purchased under a share buyback program	13 288 744	77,98%	24 877 046	99,87%	33 000	0,13%	0	0,00%	adopted

RESOLUTIONS		SHARES PRESENT	% OF THE ISSUED CAPITAL	FOR		AGAINST		ABSTENTION		RESULT
				VOICES	%	VOICES	%	VOICES	%	
14	Delegation of authority granted to the board of directors to issue shares and/or securities giving access to the share capital of the Company and/or securities giving the right to the allotment of debt instruments, with preferential subscription rights of the shareholders	13 288 744	77,98%	24 910 046	100,00%	0	0,00%	0	0,00%	adopted
15	Delegation of authority granted to the board of directors to issue shares and/or securities giving access to the share capital of the Company and/or securities giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders - through public offerings	13 288 744	77,98%	23 959 886	96,19%	950 160	3,81%	0	0,00%	adopted
16	Delegation of authority granted to the board of directors to issue shares and/or securities giving access to the share capital of the Company and/or securities giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders - through private placements as set forth in article L.411-2 II of the French Monetary and Financial Code	13 288 744	77,98%	23 886 044	95,89%	1 024 002	4,11%	0	0,00%	adopted
17	Authorization granted to the board of directors in the event of an issuance of shares and/or securities giving access to the share capital of the Company through public offerings or private placements as set forth in article L.411-2 II of the French Monetary and Financial Code, without preferential subscription rights of the shareholders, to set the issuance price under the conditions defined by the shareholders' meeting, up to a maximum of ten per cent (10%) of the share capital per year	13 288 744	77,98%	23 951 088	96,15%	958 958	3,85%	0	0,00%	adopted
18	Delegation of authority granted to the board of directors to increase the number of shares and/or securities giving access to the share capital of the Company to be issued in the event of a share capital increase with or without preferential subscription rights of the shareholders	13 288 744	77,98%	23 951 088	96,15%	958 958	3,85%	0	0,00%	adopted
19	Delegation of authority granted to the board of directors to issue shares and/or securities giving access to the share capital of the Company, as consideration for contributions in kind consisting of a company's shares and/or securities giving access to the share capital up to a maximum of ten per cent (10%) of the share capital, outside public exchange offers initiated by the Company	13 288 744	77,98%	24 838 053	99,71%	71 993	0,29%	0	0,00%	adopted
20	Delegation of authority granted to the board of directors to issue shares and/or securities giving access to the share capital of the Company, as consideration for securities tendered through public exchange offers initiated by the Company, without preferential subscription right of the shareholders	13 288 744	77,98%	23 989 463	96,30%	920 583	3,70%	0	0,00%	adopted
21	Delegation of authority granted to the board of directors to increase the share capital through the capitalization of reserves, earnings or premiums or any other sum upon which capitalization would be permitted	13 288 744	77,98%	24 910 046	100,00%	0	0,00%	0	0,00%	adopted

RESOLUTIONS		SHARES PRESENT	% OF THE ISSUED CAPITAL	FOR		AGAINST		ABSTENTION		RESULT
				VOICES	%	VOICES	%	VOICES	%	
22	Delegation of authority granted to the board of directors to issue shares and/or securities giving access to the share capital of the Company, reserved for members enrolled in an employer-sponsored company savings plan, without preferential subscription rights of the shareholders	13 288 744	77,98%	679 860	2,73%	24 230 186	97,27%	0	0,00%	rejected
23	Proxy for formalities	13 288 744	77,98%	24 910 046	100,00%	0	0,00%	0	0,00%	adopted