

SARTORIUS STEDIM BIOTECH S.A. – Annual General Shareholders' Meeting – 3 avril 2018 - VOTING RESULTS

Number of shares of the issued capital the day of the meeting	92,180,190
Number of shares with voting rights	92,180,190
Number of shares present, represented or with postal voting	83,002,685
Number of voices present, represented or with postal voting	151,612,902
Number of shares present	69,073,858
Number of voices present	137,648,385
Number of shares with postal voting	13,893,292
Number of voices with postal voting	13,894,132

Number of shares with proxy	350,535
Number of voices with proxy	70,385
Number of shareholders present, represented or with postal voting	223
Quorum	90.04%

RESOLUTIONS	SHARES PRESENT	% OF THE ISSUED CAPITAL	FOR		AGAINST		ABSTENTION		RESULT
			VOICES	%	VOICES	%	VOICES	%	
1 Approval of Financial statements for the year ended 31 December 2017 and discharge to all directors	83,002,685	90.04%	151,454,489	99.90%	158,413	0.10%	0	0.00%	approved
2 Approval of the consolidated financial statements for the year ended 31 December 2017	83,002,685	90.04%	151,588,882	99.98%	24,020	0.02%	0	0.00%	approved
3 Assignment of the financial result for the year ended 31 December 2017	83,002,685	90.04%	151,588,382	99.98%	24,520	0.02%	0	0.00%	approved
4 Approval of regulated agreements covered by Article L.225-38 and subsequent of the French Commercial Code	14,552,285	90.04%	6,800,292	46.22%	7,911,810	53.78%	0	0.00%	rejected
5 Setting of the annual Directors' fees for the members of the Board of Directors	83,002,685	90.04%	151,588,382	99.98%	24,520	0.02%	0	0.00%	approved
6 Approval of the fixe, variable and extraordinary components of the remuneration and the benefits of all kinds due or awarded to the Chairman and Chief Executive Officer for the 2017 financial year	83,002,685	90.04%	144,074,915	95.03%	7,537,987	4.97%	0	0.00%	approved
7 Approval of the criteria and principles for determining, allocating and awarding components of the compensation and benefits of the Chairman and Chief Executive Officer for the 2018 financial year	83,002,685	90.04%	141,518,086	93.34%	10,094,816	6.67%	0	0.00%	approved
8 Renewal of the term of Mrs. Anne-Marie Graffin as Director	83,002,685	90.04%	151,555,721	99.96%	57,181	0.04%	0	0.00%	approved
9 Renewal of the term of Mrs. Susan Dexter as Director	83,002,685	90.04%	151,555,721	99.96%	57,181	0.04%	0	0.00%	approved
10 Ratification of the cooptation of Mr. Lothar Kappich as Directors by the Board of Directors	83,002,685	90.04%	147,948,607	97.58%	3,664,295	2.42%	0	0.00%	approved
11 Renewal of the term of Deloitte et Associés as one the Statutory Auditors	83,002,685	90.04%	151,511,381	99.93%	101,521	0.07%	0	0.00%	approved
12 Acknowledgment of the end of term and non renewal of one of the Alternate Auditors	83,002,685	90.04%	151,588,882	99.98%	24,020	0.02%	0	0.00%	approved
13 Authorization granted to the Board of directors to enable the Company to trade in its own shares	83,002,685	90.04%	143,661,996	94.76%	7,950,906	5.24%	0	0.00%	approved
14 Proxy to carry out formalities	83,002,685	90.04%	151,588,882	99.98%	24,020	0.02%	0	0.00%	approved
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, with preferential subscription rights of the shareholders	83,002,685	90.04%	144,581,092	95.36%	7,031,810	4.64%	0	0.00%	approved
16 Delegation of authority granted to the Board of directors to issue shares and/or securities giving or capable of 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	83,002,685	90.04%	143,170,999	94.43%	8,441,903	5.57%	0	0.00%	approved

17	Delegation of authority granted to the Board of directors to issue shares and/or securities giving or capable of giving access to the share capital of the Company and/or the issuance of securities giving the right to the allotment of debt instruments, without the preferential subscription rights of the shareholders – through private placements referred to in article L. 411-2 II of the French Monetary and Financial Code	83,002,685	90.04%	142,837,354	94.21%	8,775,548	5.79%	0	0.00%	approved
18	Delegation of authority to the Board of directors to increase the number of shares and/or securities giving or capable of giving access to the share capital of the Company to be issued in case of share capital increase, with or without preferential subscription rights of the shareholders	83,002,685	90.04%	143,196,290	94.45%	8,416,612	5.55%	0	0.00%	approved
19	Delegation of authority granted to the Board of directors to decide to issue shares and/or securities giving or capable of giving access to the share capital of the Company as consideration for contributions in kind in shares and/or securities giving or capable of giving access to capital, without preferential subscription rights of shareholders	83,002,685	90.04%	144,574,596	95.36%	7,038,306	4.64%	0	0.00%	approved
20	Delegation of authority granted to the Board of directors to increase the share capital of the Company through the capitalization of reserves, earnings or premiums or any other sum upon which capitalization would be permitted	83,002,685	90.04%	144,720,503	95.45%	6,892,399	4.55%	0	0.00%	approved
21	Delegation of authority granted to the Board of directors to issue shares and/or securities giving or capable of giving access to the share capital of the Company, reserved for members of company savings plan, without preferential subscription rights of the shareholders	83,002,685	90.04%	151,190,741	99.72%	422,161	0.28%	0	0.00%	approved