

SARTORIUS STEDIM BIOTECH S.A. – Annual General Shareholders' Meeting – 26 March 2019 – VOTING RESULTS

Number of shares of the issued capital the day of the meeting	92.180.190
Number of shares with voting rights	92.177.210
Number of shares present, represented or with postal voting	82.157.386
Number of total voting rights	160.973.737
Number of voices present, represented or with postal voting	150.770.982
Number of shares present	68.613.512
Number of voices present	137.180.874
Number of shares with postal voting	13.498.046
Number of voices with postal voting	13.499.376

Number of shares with proxy	22.914
Number of voices with proxy	45.366
Number of shareholders present, represented or with postal voting	405
Quorum	89,13%

	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 2018 and discharge to all directors	82.157.386	89,13%	149.721.716	99,30%	1.003.900	0,67%	0	0,00%	approved
2	Approval of the consolidated financial statements for the year ended 31 December 2018	82.157.386	89,13%	150.648.967	99,92%	76.649	0,05%	0	0,00%	approved
3	Assignment of the financial result for the year ended 31 December 2018	82.157.386	89,13%	150.694.871	99,95%	30.745	0,02%	0	0,00%	approved
4	Approval of regulated agreements covered by Article L.225-38 and subsequent of the French Commercial Code	13.706.986	57,77%	3.804.346	27,43%	9.740.396	70,23%	0	0,00%	rejected
5	Approval of regulated commitments covered by article L. 225-42-1 of the French Commercial Code related to Mr. Joachim Kreuzburg, Chief Executive Officer	13.706.986	57,77%	3.263.751	23,53%	10.280.991	74,12%	0	0,00%	rejected
6	Setting of the annual Directors' fees for the members of the Board of Directors	82.157.386	89,13%	150.695.730	99,95%	29.886	0,02%	0	0,00%	approved
7	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 2018 financial year	82.157.386	89,13%	140.864.901	93,43%	9.860.715	6,54%	0	0,00%	approved
8	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 2019 financial year	82.157.386	89,13%	140.867.799	93,43%	9.857.817	6,54%	0	0,00%	approved
9	Renewal of the term of Mr. Joachim Kreuzburg as Director	82.157.386	89,13%	145.213.558	96,31%	5.512.058	3,66%	0	0,00%	approved
10	Renewal of the term of Mr. Lothar Kappich as Director	82.157.386	89,13%	145.571.224	96,55%	5.154.392	3,42%	0	0,00%	approved
11	Renewal of the term of M. Henri Riey as Director	82.157.386	89,13%	147.502.494	97,83%	3.223.122	2,14%	0	0,00%	approved
12	Nomination of Mrs. Pascale Boissel as new member of the Board of Directors of the Company	82.157.386	89,13%	142.688.823	94,64%	8.036.793	5,33%	0	0,00%	approved
13	Nomination of Mr. René Fäber as new member of the Board of Directors of the Company	82.157.386	89,13%	148.397.527	98,43%	2.328.089	1,54%	0	0,00%	approved
14	Authorization granted to the Board of directors to enable the Company to trade in its own shares	82.157.386	89,13%	143.068.508	94,89%	7.657.108	5,08%	0	0,00%	approved
15	Proxy to carry out formalities	82.157.386	89,13%	150.696.815	99,95%	28.801	0,02%	0	0,00%	approved
16	Update of Bylaws; adoption of the new Bylaws	82.157.386	89,13%	140.215.111	93,00%	10.510.505	6,97%	0	0,00%	approved
17	Update of Bylaws for determine the procedures for appointing directors and directors representing employees; update of the article 15 of Bylaws	82.157.386	89,13%	150.645.376	99,92%	80.240	0,05%	0	0,00%	approved
18	Proxy to carry out formalities	82.157.386	89,13%	150.696.815	99,95%	28.801	0,02%	0	0,00%	approved