

SARTORIUS STEDIM BIOTECH S.A. – Annual General Shareholders' Meeting –4 April 2017 – 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	82.984.391
Number of voices present, represented or with postal voting	152.593.391
Number of shares present	69.574.406
Number of voices present	139.148.376
Number of shares with postal voting	13.373.835
Number of voices with postal voting	13.373.865

Number of shares with proxy	36.150
Number of voices with proxy	71.150
Number of shareholders present, represented or with postal voting	227
Quorum	90,02%

	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 2016 and discharge to all directors	82.984.391	90,02%	152.267.736	99,79%	325.655	0,21%	0	0,00%	approved
2	Approval of the consolidated financial statements for the year ended 31 December 2016	82.984.391	90,02%	150.715.573	98,77%	1.877.818	1,23%	0	0,00%	approved
3	Assignment of the financial result for the year ended 31 December 2016 and setting of the dividend	82.984.391	90,02%	152.593.391	100,00%	0	0,00%	0	0,00%	approved
4	Approval of related party agreements as referred in articles L.225-38 and subsequent of the commercial code	14.533.985	61,25%	8.217.675	52,37%	7.474.916	47,63%	0	0,00%	approved
	<i>Being specified that shares who are parties didn't vote</i>									
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	14.533.985	61,25%	6.850.589	43,65%	8.842.002	56,35%	0	0,00%	rejected
	<i>Being specified that shares who are parties didn't vote</i>									
6	Setting of the annual Directors' fees for the members of the Board of Directors	82.984.391	90,02%	152.478.859	99,92%	114.532	0,08%	0	0,00%	approved
7	Approval of the elements of compensation due or granted for the 2016 financial year to Mr Joachim Kreuzburg, Chief Executive Officer	82.984.391	90,02%	144.086.087	94,42%	8.507.304	5,58%	0	0,00%	approved
8	Approval of the elements of compensation due or granted for the 2016 financial year to Mr Volker Niebel, Executive Vice President	82.984.391	90,02%	144.502.153	94,70%	8.091.238	5,30%	0	0,00%	approved
9	(Approval of the elements of compensation due or granted for the 2016 financial year to Mr Oscar-Werner Reif, Executive Vice President	82.984.391	90,02%	144.502.153	94,70%	8.091.238	5,30%	0	0,00%	approved
10	Approval of the elements of compensation due or granted for the 2016 financial year to Mr Reinhard Vogt, Executive Vice President)	82.984.391	90,02%	144.502.153	94,70%	8.091.238	5,30%	0	0,00%	approved
11	Proxy to carry out formalities)	82.984.391	90,02%	152.593.391	100,00%	0	0,00%	0	0,00%	approved