



**ANNUAL GENERAL SHAREHOLDER'S MEETING  
OF 26 MARCH 2019**

The Annual Combined General Shareholders' meeting of Sartorius Stedim Biotech met the 26<sup>th</sup> of March 2019 at the registered office in Aubagne, under the chairmanship of Mr Joachim Kreuzburg, Chairman and Chief Executive Officer of the Company.

The 405 shareholders who attended the meeting represented 89.13 % of shares carrying a voting right, adopted all resolutions submitted to them, except the fourth and fifth resolutions.

The complete results of the vote of the resolutions are available on the Company's website <http://www.sartorius.com/sartorius/en/EUR/company/investor-relations/sartorius-stedim-biotech-sa/shareholders-meeting/2019/>.

The shareholders' meeting was the occasion for Mr Joachim Kreuzburg to remind the good results of the Company and the Group.

He also presented to the shareholders the Group and the Company strategy and perspectives of development.

\*\*\*\*\*