

Aubagne, France, March 24, 2021

Annual Combined Shareholders' Meeting of 24 March 2021

The Annual Combined General Shareholders' meeting of Sartorius Stedim Biotech was convened on the 24th of March 2021, under the chairmanship of Mr. Joachim Kreuzburg, Chairman and Chief Executive Officer of the Company. In accordance with the French Ordinance N°. 2020-321 of March 25, 2020, modified by the French Ordinance N° 2020- 1497 of 2, December 2020, that have adapted the rules for shareholder meetings and conferences, the meeting was held without the physical attendance of shareholders and was fully live-streamed on the company's website at www.sartorius.com due to the coronavirus pandemic and the associated restrictions imposed on travel and in-person meetings.

The 2 220 shareholders who expressed their votes represented 88,82% of shares carrying a voting right, adopted all resolutions submitted to them, except the fourth resolution.

The complete results of the vote of the resolutions are available on the Company's website.

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

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