

Aubagne, France, March 27, 2023

The Annual Combined General Shareholders' meeting of Sartorius Stedim Biotech met on March 27, 2023 at the registered office in Aubagne, under the chairmanship of Mrs. Anne-Marie Graffin, member of the Board of Directors, associated with the videoconferencing of Mr. Joachim Kreuzburg, Chairman of the Board of Directors and Chief Executive Officer of the Company.

The 3251 shareholders who expressed their votes represented 92.89 % of shares carrying a voting right, adopted all resolutions submitted to them.

Mr. Joachim Kreuzburg presented the figures of the successful fiscal year 2022, as well as the Group and the Company's strategy and its future perspectives to the shareholders.

Complete results of the vote of the resolutions are available on the Company's website :

<https://www.sartorius.com/en/company/investor-relations/sartorius-stedim-biotech-sa-investor-relations/shareholders-meeting>