

Aubagne, France, March 26, 2024

The Annual Combined General Shareholders' meeting of Sartorius Stedim Biotech met on March 26, 2024 at its registered office in Aubagne, under the chairmanship of Mr. Kreuzburg, Chairman of the Board of Directors and in the presence of Mr René Faber, Chief Executive Officer of the Company.

The 2 882 shareholders who expressed their votes represented 93,45 % of shares carrying a voting right, adopted all resolutions submitted.

Mr. René Faber presented to the shareholders the numbers of the fiscal year 2023, as well as the Group and the Company's strategy, along with future developments.

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

[Shareholders' Meeting | Sartorius Stedim Biotech S.A.](#)