

Aubagne, France, March 24, 2026

The Annual Combined General Shareholders' meeting of Sartorius Stedim Biotech met on March 24, 2026 at its registered office in Aubagne, under the chairmanship of M. Michael Grosse, Chairman of the Board of Directors and in the presence of M. René Faber, Chief Executive Officer of the Company.

The 2478 shareholders who expressed their votes represented 89,90% of shares carrying voting rights, adopted all resolutions submitted.

Mr. René Faber presented to the shareholders the results of the fiscal year 2025, as well as the Group and the Company's future developments.

Complete results of the vote of the resolutions are available on the Company's website: [Shareholders' Meeting | Sartorius Stedim Biotech S.A.](#)