

Sartorius Stedim Biotech S.A. - 13400 Aubagne, France

Sartorius Stedim Biotech SA
ZI les Paluds Avenue de Jouques
13781 Aubagne Cedex

t:+33.(0)4.42.84.56.00
f:+33.(0)4.42.84.56 19
www.sartorius-stedim.com

S.A. au capital de 19,466,081.00€
RCS de Marseille B 314 093 352
SIRET 314 093 352 00073 - APE 6420Z
TVA FR 49 314 093 352

**DECLARATION RELATIVE TO
THE NUMBER OF SHARES AND VOTING RIGHTS
MAKING UP THE ISSUED CAPITAL**

Article L 233-8-II of the Code of commerce
and article 223-6 of the general regulations of the AMF

Designation of issuer

SARTORIUS STEDIM BIOTECH SA
CS 91051
Z. I. Les Paluds
Avenue de Jouques
13400 Aubagne

Date of information	Total number of shares making up the issued capital	Total number of theoretical voting rights including treasury shares	Total number of net voting rights
February 28, 2026	97,330,405	164,793,754	164,751,543