

Sartorius Stedim Biotech S.A. - 13781 Aubagne Cedex, France

Sartorius Stedim Biotech SA  
Z.I. Les Paluds  
Avenue de Jouques, CS 91051  
13781 Aubagne Cedex, France

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 18,436,038 €  
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  
13781 Aubagne Cedex

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
March 31, 2023	92,180,190	160,432,226	160,416,259