

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

**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 222-12-5 of the general regulations of the AMF

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

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 voting rights
February 28, 2017	92,180,190	162,041,944