

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 223-6 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

<b>Date of information</b>	<b>Total number of shares making up the issued capital</b>	<b>Total number of theoretical voting rights including treasury shares</b>	<b>Total number of net voting rights</b>
August 31, 2019	92,180,190	160,981,410	160,979,030