

Sartorius Stedim Biotech S.A. - 13781 Aubagne 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 10 386 682,28 €
RCS de Marseille B 314 093 352
SIRET 314 093 352 00073 – APE 252 H
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

Deadline date for information	Total number of shares making up the issued capital	Total number of voting rights
January 7 th , 2013	17,041,306	Gross total of voting rights: 28,664,315
		Net total *of voting rights: 26,965,605

* Net total = total number of voting rights attached to shares – private share voting right