

Sartorius Stedim Biotech S.A. 13781 Aubagne France

Sartorius Stedim Biotech S.A.
Z.I. des Paluds
Avenue Jouques, B.P. 1051
13781 Aubagne, 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 10 301 672,68 €
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
BP 1051
Z I des 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
October 31st, 2007	16 897 988	Gross total of voting rights: 18 912 368
		Net total of voting rights: 18 912 368

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