

Sartorius Stedim Biotech S.A. - 13781 Aubagne France

Sartorius Stedim Biotech S.A.
Z.I. Les Paluds
Avenue de Jouques, B.P. 1051
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 10 322 717,68 €
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 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 |
|--------------------------------------|--|--|
| June 4 th , 2009 | 16 951 888 | Gross total of voting rights: 18 966 518 |
| | | Net total of voting rights: 18 966 518 |

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