

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
July 31st, 2007	16 887 988	Gross total of voting rights: 18 902 368
		Net total of voting rights: 18 902 368

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