

Sartorius Stedim Biotech S.A. - 13781 Aubagne France

Sartorius Stedim Biotech SA
 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 10 386 682,28 €
 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
 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
March 7 th , 2012	17,028,448	Gross total of voting rights: 28,860,461
		Net total *of voting rights: 27,161,751

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