

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

**TOTAL NUMBER OF SHARES
 AND
 TOTAL NUMBER OF VOTING RIGHTS
 MAKING UP THE ISSUED CAPITAL**

S.A. au capital de 10 384 303,28 €
 RCS de Marseille B 314 093 352
 SIRET 314 093 352 00073 – APE 252 H
 TVA FR 49 314 093 352

Deadline date for information	Total number of shares making up the issued capital	Total number of voting rights
March 14 th , 2011	17,023,448	Gross total of voting rights: 18,674,142
		Net total *of voting rights: 16,975,432

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