

Sartorius Stedim Biotech S.A. - 13781 Aubagne France

Sartorius Stedim Biotech SA  
 Z.I. Les Paluds  
 Avenue de Jouques, CS 91051  
 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 10 395 806.66 €  
 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  
 CS 91051  
 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 11 <sup>th</sup> , 2013	17,042,306	Gross total of voting rights: 28,665,315
		Net total *of voting rights: 26,966,605

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