



Disclosure of trading in own shares

Sartorius Stedim Biotech
 S.A. with capital of 10,369,626.68 €
 RCS of Marseille B 314 093 352
 SIRET 314 093 352 00073 – APE 252 H
 TVA FR 49 314 093 352

Sartorius Stedim Biotech S.A.
 Z.I. Les Paluds
 Avenue de Jouques
 B.P. 1051
 13781 Aubagne - France
www.sartorius-stedim.com

Goettingen, May 28, 2010

Pursuant to applicable law on share repurchases, Sartorius Stedim Biotech declares the following purchases of its own shares:

Investor Relations:

Andreas Wiederhold
 Phone: +49 (0) 551 308-1668
andreas.wiederhold@sartorius.com
sartorius.com

Trading session	Number of shares purchased	Weighted average price (in euros)	Amount (in euros)
May 24, 2010	543	31.97	17,361.18
May 25, 2010	2,842	30.47	86,603.98
Total	3,385	30.71	103,965.16

No derivative financial instruments were used to purchase these shares.

Contact: Andreas Wiederhold; Team Leader Treasury & Investor Relations ;
 Phone: +49 (0)551.308-1668; Fax: +49 (0)551.308-3153 ;
 e-mail: andreas.wiederhold@sartorius.com ; www.sartorius-stedim.com