



## 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](http://www.sartorius-stedim.com)

Goettingen, May 21, 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](mailto:andreas.wiederhold@sartorius.com)  
[sartorius.com](http://sartorius.com)

Trading session	Number of shares purchased	Weighted average price (in euros)	Amount (in euros)
May 17, 2010	1,273	35.55	45,248.91
May 18, 2010	2,215	35.30	78,191.27
May 19, 2010	800	34.50	27,600.00
May 20, 2010	2,446	33.91	82,954.87
May 21, 2010	1,369	32.25	44,147.79
<b>Total</b>	<b>8,103</b>	<b>34.33</b>	<b>278,142.84</b>

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](mailto:andreas.wiederhold@sartorius.com) ; [www.sartorius-stedim.com](http://www.sartorius-stedim.com)