



## Disclosure of trading in own shares

Sartorius Stedim Biotech  
 S.A. with capital of 10,375,726.68 €  
 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, Dec. 15, 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)
Dec. 09, 2010	370,000	34.32	12,698,400.00
<b>Total</b>	<b>370,000</b>	<b>34.32</b>	<b>12,698,400.00</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)