

**REQUEST FOR DOCUMENTS AND INFORMATION****COMBINED ORDINARY AND EXTRAORDINARY ANNUAL SHAREHOLDERS' MEETING 7 APRIL 2015**

To send back to:

**SARTORIUS STEDIM BIOTECH S.A.**  
**To the attention of Mr Olivier Guitard**  
**ZI Les Paluds - Avenue de Jouques**  
**CS 91051**  
**13781 Aubagne Cedex**

I, the undersigned

Name: .....

Forename: .....

Registered name: .....

Represented by: .....

Acting as: .....

Home address / Registered office: .....

.....

Owner of ..... registered shares,

And/or ..... bearer shares, registered in the accounts of (bank, financial institution) .....,

Asks to receive documents and information, concerning the Combined Ordinary and Extraordinary Annual Shareholders' Meeting of 7 April 2015, referred to in the provisions of Article R.225-88 of the French commercial code :

- To my address indicated above,
- To the following postal address: .....

Done at ..... on .....

(Signature)

The documents and information referred to in the Articles R.225-81 and R.225-83 of the French commercial code are available on the website of the company.  
According to the Article R.225-88 subparagraph 3 of the French commercial code, shareholders owners of registered shares can, by a single request, receive from the company documents and information at each of the next annual shareholders' meetings. To this facility, check this box. :