

To be returned to:

SARTORIUS STEDIM BIOTECH S.A.
Legal department
Z.I. Les Paluds -
Avenue de Jouques
13400 Aubagne

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 Annual Combined General Shareholders' Meeting of 26 March 2024, 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: