
POWER OF ATTORNEY

Entrance card no.: _____

No. of shares: _____

I/we

First Name of Depositor

Last Name of Depositor

Depositor's Postal Code (ZIP)

Depositor's Complete Address (Street, City, Country)

hereby authorize Mr./Ms./Mrs.

Proxy's First Name

Proxy's Last Name

Proxy's Postal Code (ZIP)

Proxy's Complete Address (Street, City, Country)

to grant substitute power of attorney, and, by way of exemption from the restrictions set out in § 181 Alternative 2 of the German Civil Code (BGB), to represent me/us at the Annual General Shareholders' Meeting of the Sartorius Aktiengesellschaft on April 10, 2014, and to exercise my/our voting rights.

Place/Date

Depositor's Signature