

Annual General Shareholders' Meeting on March 28, 2019



Power of Attorney

Entrance card no. _____

No. of shares _____

I/we (Depositor)

First Name

Last Name

Postal Code (ZIP)

Complete Address (Street, City, Country)

hereby authorize

Mr./Ms./Mrs. (Proxy)

First Name

Last Name

Postal Code (ZIP)

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 **March 28, 2019**, and to exercise my/our rights (in particular voting rights).

Place/Date

Depositor's Signature