

Virtual General Meeting of Sartorius Aktiengesellschaft on June 26, 2020

Power of Attorney Granted to a Third Party

ID number confirming registration (AGM ticket): _____ Last name | Company: _____

Number of shares: _____ First name: _____

Phone number*: _____ Email address*: _____

Please return **no later than June 25, 2020, at 6:00 p.m. (CEST)** (deadline for receipt by the company) to:

Sartorius Aktiengesellschaft
c/o Better Orange IR & HV AG
Haidelweg 48
81241 Munich, Germany

or by **email**: sartorius@better-orange.de
or by **fax**: +49 (0)89 889 690 655

I | we hereby grant power of attorney to Mr. | Ms

Last name of proxy: _____

First name of proxy: _____

Proxy's place of residence: _____

if necessary by revoking a power of attorney already granted at an earlier time, to represent me | us in exercising my | our shareholder rights with respect to the virtual general meeting of Sartorius Aktiengesellschaft on June 26, 2020, and to exercise my | our voting rights by disclosing my | our name(s). This power of attorney includes the right of granting substitute power of attorney. I | we shall provide my | our individual access data for the password-protected Internet service platform to said proxies and draw their attention to the special condition that they are permitted to disclose this access data to third parties only within the scope of granting substitute power of attorney to substitute proxies for this virtual general meeting.

Place

Date

Legible signature(s) of person(s) making this declaration