

## Financial Schedule

Annual Shareholders' Meeting in Aubagne, France	March 24, 2020
Payment of dividends <sup>1)</sup>	March 31, 2020
Publication of first-quarter figures for 2020	April 21, 2020
Publication of first-half figures for 2020	July 21, 2020
Publication of nine-month figures for 2020	October 20, 2020
Publication of preliminary figures for fiscal 2020	January 2021
Annual Shareholders' Meeting in Aubagne, France	March, 2021
Publication of first-quarter figures for 2021	April 2021

<sup>1)</sup> Subject to approval by the Annual Shareholders' Meeting

### Contacts

**Petra Kirchhoff**  
Head of  
Corporate Communications & IR

Phone: 0551.308.1686  
petra.kirchhoff@sartorius.com

**Ben Orzelek**  
Head of Investor Relations

Phone: 0551.308.1668  
ben.orzelek@sartorius.com

### About This Publication

**Published by**  
Sartorius AG  
Corporate Communications  
37070 Goettingen, Germany

**Editorial Deadline**  
February 17, 2020

**Published on**  
February 18, 2020

**Financial Reporting System firesys**  
firesys GmbH  
Frankfurt | Main, Germany

**Photography**  
Peter Ginter  
Lohmar, Germany  
Frank Stefan Kimmel  
Goettingen

This is a translation of the original  
French-language  
"Document d'Enregistrement  
Universel 2019".