

## Financial Schedule

Annual Shareholders' Meeting in Goettingen, Germany	March 28, 2019
Payment of dividends <sup>1)</sup>	April 2, 2019
Publication of first-quarter figures for 2019	April 18, 2019
Publication of first-half figures for 2019	July 19, 2019
Publication of nine-month figures for 2019	October 22, 2019
German Equity Forum in Frankfurt   Main, Germany	November 25-27, 2019
Publication of preliminary figures for fiscal 2019	January 2020
Annual press conference in Goettingen, Germany	February 2020
Annual Shareholders' Meeting in Goettingen, Germany	March 26, 2020
Publication of first-quarter figures for 2020	April 2020

<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

#### Katharina Tillmanns-Pelzer

Manager Corporate Responsibility

Phone: 0551.308.2893  
katharina.tillmanns@sartorius.com

### About This Publication

#### Published by

Sartorius AG  
Corporate Communications  
37070 Goettingen, Germany

#### Editorial Deadline

February 18, 2019

#### Published on

February 19, 2019

#### 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  
German-language annual report.