

## Financial Schedule

|                                                    |                  |
|----------------------------------------------------|------------------|
| Annual Shareholders' Meeting<br>in Aubagne, France | March 26, 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 |
| Publication of preliminary figures for fiscal 2019 | January 2020     |
| Annual Shareholders' Meeting<br>in Aubagne, France | March 24, 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

### About This Publication

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

**Editorial Deadline**  
February 15, 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  
French-language  
"Document de Référence 2018".