

Financial Schedule

Annual Shareholders' Meeting in Goettingen, Germany	March 26, 2020
Payment of dividends ¹⁾	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 press conference in Goettingen, Germany	February 2021
Annual Shareholders' Meeting in Goettingen, Germany	March, 2021
Publication of first-quarter figures for 2020	April 2021

¹⁾ 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.