

1312519 25.03.2022 CET/CEST

Englisch

DGAP Voting Rights Announcement: SARTORIUS AG

Release according to Article 40, Section 1 of the WpHG [the German Securities Trading Act] with the objective of Europe-wide distribution

25.03.2022 / 16:10 CET/CEST

Dissemination of a Voting Rights Announcement transmitted by DGAP - a service of EQS Group AG.

The issuer is solely responsible for the content of this announcement.

Notification of Major Holdings

1. Details of issuer

Name:	SARTORIUS AG
Street:	Otto-Brenner-Straße 20
Postal code:	37079
City:	Göttingen Germany
Legal Entity Identifier (LEI):	529900EQV2DY4FOAMU38

2. Reason for notification

<input checked="" type="checkbox"/>	Acquisition/disposal of shares with voting rights
<input type="checkbox"/>	Acquisition/disposal of instruments
<input type="checkbox"/>	Change of breakdown of voting rights

Other reason:

3. Details of person subject to the notification obligation

Natural person (first name, surname): Alice N. Schwartz
Date of birth: 02 Jul 1926

4. Names of shareholder(s)

holding directly 3% or more voting rights, if different from 3.

Bio-Rad Laboratories GmbH

5. Date on which threshold was crossed or reached:

01 Oct 2013

6. Total positions

	% of voting rights attached to shares (total of 7.a.)	% of voting rights through instruments (total of 7.b.1 + 7.b.2)	Total of both in % (7.a. + 7.b.)	Total number of voting rights pursuant to Sec. 41 WpHG
New	33.28 %	0.00 %	33.28 %	9,360,000
Previous notification	n/a %	n/a %	n/a %	/

7. Details on total positions

a. Voting rights attached to shares (Sec. 33, 34 WpHG)

ISIN	Absolute		In %	
	Direct (Sec. 33 WpHG)	Indirect (Sec. 34 WpHG)	Direct (Sec. 33 WpHG)	Indirect (Sec. 34 WpHG)
DE0007165607	0	3,115,201	0.00 %	33.28 %
Total	3,115,201		33.28 %	

b.1. Instruments according to Sec. 38 (1) no. 1 WpHG

Type of instrument	Expiration or maturity date	Exercise or conversion period	Voting rights absolute	Voting rights in %
			0	0.00 %
		Total	0	0.00 %

b.2. Instruments according to Sec. 38 (1) no. 2 WpHG

Type of instrument	Expiration or maturity date	Exercise or conversion period	Cash or physical settlement	Voting rights absolute	Voting rights in %
				0	0.00 %
			Total	0.00	%

8. Information in relation to the person subject to the notification obligation

Person subject to the notification obligation is not controlled nor does it control any other undertaking(s) that directly or indirectly hold(s) an interest in the (underlying) issuer (1.).

X	Full chain of controlled undertakings starting with the ultimate controlling natural person or legal entity:
---	--

Name	% of voting rights (if at least 3% or more)	% of voting rights through instruments (if at least 5% or more)	Total of both (if at least 5% or more)
Alice N. Schwartz	%	%	%
David Schwartz Non-Exempt Marital Trust	%	%	%
Blue Raven Partners, L.P.	%	%	%
Bio-Rad Laboratories, Inc.	%	%	%
Bio-Rad Luxembourg S.à r.l.	%	%	%
Bio-Rad France Holding SAS	%	%	%
Bio-Rad Laboratories GmbH	33.28 %	%	33.28 %

9. In case of proxy voting according to Sec. 34 para. 3 WpHG
(only in case of attribution of voting rights in accordance with Sec. 34 para. 1 sent. 1 No. 6 WpHG)

Date of general meeting:

Holding total positions after general meeting (6.) after annual general meeting:

Proportion of voting rights	Proportion of instruments	Total of both
%	%	%

10. Other explanatory remarks:

Voting rights with a notification requirement have been held directly by Bio-Rad Laboratories GmbH since October 3, 2003 and were gradually increased in subsequent years, most recently to 30.01% on April 1, 2011. The voting rights were indirectly attributable to Alice N Schwartz at all times, but no notifications were made. On October 1, 2013, the Group was restructured, which is reflected in this one-time notification.

Date

24 March 2022

25.03.2022 CET/CEST The DGAP Distribution Services include Regulatory Announcements, Financial/Corporate News and Press Releases.
Archive at www.dgap.de

Language: English
Company: SARTORIUS AG
Otto-Brenner-Straße 20
37079 Göttingen
Germany
Internet: www.sartorius.com

End of News

DGAP News Service

