

Sartorius AG



WKN: 716560 ISIN: DE0007165607 Land: Deutschland

Nachricht vom 20.03.2018 | 14:34

**Correction of a release from 25.09.2017, 17:05 CET/CEST - Sartorius AG:
Release according to Article 26, Section 1 of the WpHG [the German
Securities Trading Act] with the objective of Europe-wide distribution**

SARTORIUS AG

20.03.2018 / 14:34

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.

Correction of a notification of Major Holding published on 25.09.2017

1. Details of issuer

SARTORIUS AG
Otto-Brenner-Straße 20
37079 Göttingen
Germany

2. Reason for notification

	Acquisition/disposal of shares with voting rights
	Acquisition/disposal of instruments
	Change of breakdown of voting rights
X	Other reason: Change in execution of last will

3. Details of person subject to the notification obligation

Name:	City and country of registered office:
Mr Dr. Lothar Kappich, Date of birth: 15 Feb 1957	

4. Names of shareholder(s)

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

Karin Sartorius-Herbst, Christine Franken, Ulrike Baro, Andreas Franken, Kai Christian Franken (in Erbengemeinschaft)

5. Date on which threshold was crossed or reached:

08 Sep 2017

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 of issuer
Resulting situation	50.09 %	0 %	50.09 %	37,440,000
Previous notification	n/a %	n/a %	n/a %	/

7. Notified details of the resulting situation

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

ISIN	absolute	in %

	direct (Sec. 33 WpHG)	indirect (Sec. 34 WpHG)	direct (Sec. 33 WpHG)	indirect (Sec. 34 WpHG)
DE0007165607	0	18754160	0 %	50.09 %
Total	18,754,160		50.09 %	

b.1. Instruments according to Sec. 38 para. 1 No. 1 WpHG

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

b.2. Instruments according to Sec. 38 para. 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 %
					%
			Total		%

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

<input checked="" type="checkbox"/>	Person subject to the notification obligation is not controlled and does itself not control any other undertaking(s) holding directly or indirectly an interest in the (underlying) issuer (1.).
	Full chain of controlled undertakings starting with the ultimate controlling natural person or legal entity:

Name	% of voting rights (if at least held 3% or more)	% of voting rights through instruments (if at least held 5% or more)	Total of both (if at least held 5% or more)

9. In case of proxy voting according to Sec. 34 para. 3 WpHG

Date of general meeting:	
Holding position after general meeting:	% (equals voting rights)

10. Other explanatory remarks:

Notification due to the death of the former executor Prof. Dr. Arnold Picot on July 9, 2017, and appointment of Dr. Kappich as his successor, effective as of September 8, 2017.
--

20.03.2018 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