

Voting Rights Announcement: SARTORIUS AG

Release according to Article 26, Section 1 of the WpHG [the German Securities Trading Act] with the objective of Europe-wide distribution

1. Details of issuer

SARTORIUS AG Weender Landstraße 94 - 108 37075 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:

07 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 21, 22 WpHG)

ISIN	absolute		in %	
	direct (Sec. 21 WpHG)	indirect (Sec. 22 WpHG)	direct (Sec. 21 WpHG)	indirect (Sec. 22 WpHG)
DE0007165607	0	18754160	0 %	50.09 %
Total	18,754,160		50.09 %	

b.1. Instruments according to Sec. 25 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. 25 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. 22 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 7, 2017.

Language: English
Company: SARTORIUS AG
Weender Landstraße 94 - 108
37075 Göttingen
Germany
Internet: www.sartorius.com

End of News

25.09.2017