

**Notification of transactions made by persons discharging managerial responsibilities and persons closely related to them**

**1. Contact information for the person discharging managerial responsibilities or the closely related person**

Company name: Sartorius AG  
Legal form: AG (*Aktiengesellschaft*)

**2. Reason for notification**

The declarant is a closely related person referred to in Article 3.26) of EU Regulation 596/2014 of April 16, 2014 on market abuse.

Please specify the positions held within the issuer.

Positions: Joachim Kreuzburg is chairman and CEO of Sartorius Stedim Biotech SA and chairman of the management board of Sartorius AG.

This is an initial notification.

**3. Contact information of the issuer**

Company name: Sartorius Stedim Biotech S.A.  
LEI: 52990006IVXY7GCSSR39

**4. Information about the transaction**

Date of the transaction: 02/11/2020  
Place of the transaction: outside a trading venue  
Nature of the transaction: sale  
Description of the financial instrument: shares  
Identification code: FR0013154002

**Detailed information by transaction:**

Unit price: EUR 327.00  
Volume: 405,887 shares

**Aggregated information:**

Weighted average price: EUR 327.00  
Volume: 405,887 shares

The transaction is NOT related to the exercise of a stock option program or to a bonus or performance share grant.

## **5. Comments**

Dr. René Fáber is a member of the board of Sartorius Stedim Biotech SA and a member of the management board of Sartorius AG. Dr. Lothar Kappich is a member of the board of Sartorius Stedim Biotech SA and chairman of the supervisory board of Sartorius AG.