

Sartorius Stedim Biotech S.A.
Z.I. Les Paluds
Avenue de Jouques, CS 91051
13781 Aubagne Cedex, France

Phone +33 (0)4.42.84.5600
Fax +33 (0)4.42.84.5619

Web : www.sartorius-stedim.com

Publication on Regulated Agreements
Pursuant to Article L225-40-2 of the French Commercial Code

CONCLUSION OF AN AGREEMENT WITH SARTORIUS AG
(Approval of the Board of Directors on February 5, 2021)

General Assistance and Administrative Services Agreement between Sartorius Stedim Biotech S.A. and Sartorius AG:

- Purpose of the Agreement: Recharge from Sartorius AG to Sartorius Stedim Biotech of Managers Directors services performed to the benefit of Sartorius Stedim Biotech.
- Duration: 1 year (2021)
- Identity of the related party: Sartorius AG
- Interested persons: Joachim Kreuzburg and René Fáber
- Nature of the relationship between the related party and the contracting party: main shareholder
- Date of the Regulated Agreement: 04/01/2021
- Amount of the transaction:
 - o Amount recharged for Mr. Kreuzburg's services (budget): 883 200 Euros
 - o Amount recharged for Mr. Fáber's services (budget): 511 600 Euros
- Interest of the regulated agreement for Sartorius Stedim Biotech: appropriate allocation of Managers Directors' respective costs between all concerned entities within the Sartorius Group.
- Termination : 31/12/2021