

Special Report of the Statutory Auditors on Related Party Agreements

This is a translation into English of the statutory auditors' Special report on related party agreements of the Company issued in French and it is provided solely for the convenience of English speaking users.

This report should be read in conjunction with, and construed in accordance with, French law and professional auditing standards applicable in France.

General meeting of shareholders to approve the financial statements for the year ended 31 December 2020.

To the Shareholders of SARTORIUS STEDIM BIOTECH S.A.,

As statutory auditors of your company, we present to you our report on related party agreements.

We are required to inform you, on the basis of the information provided to us, of the terms and conditions including the reasons justifying their benefit to the company, of the related party agreements that we have been informed of or that we may have identified in the performance of our engagement. We are not required to comment as to whether they are beneficial or appropriate or to ascertain the existence of any such agreements. It is your responsibility, in accordance with article R. 225-31 of the French commercial code ('Code de Commerce'), to evaluate the benefits resulting from these agreements prior to their approval.

In addition, we are required, where applicable, to inform you in accordance with article R. 225-31 of the French commercial code concerning the implementation, during the year, of the agreements already approved by the general meeting of shareholders.

We performed those procedures which we considered necessary to comply with professional guidance issued by the French national auditing body ('Compagnie nationale des commissaires aux comptes') relating to this type of engagement. These procedures consisted in verifying that the information provided to us is consistent with the documentation from which it has been extracted.

Related Party Agreements Submitted for Approval by the General Meeting of Shareholders

Related party agreements authorized and concluded during the previous accounting period

We hereby inform you that we have not been advised of any agreements authorized and concluded during the previous accounting period to be submitted to the general meeting of shareholders for their approval in accordance with article L. 225-38 of the French Commercial Code.

Related party agreements from prior years not approved by the General meeting of the shareholders

We hereby inform you that the following agreement, authorized and concluded during the year ended 31 December 2018, which was mentioned in our special report on related party agreements for the year ended 31 December 2019 and which was not approved by the general meeting of shareholders approving the financial statements for the year ended 31 December 2019.

- General Assistance and Administrative Services Agreement
- With the company, Sartorius AG (SAG) 73.8% shareholder of the company Sartorius Stedim Biotech S.A. (SSB S.A.)
- Persons concerned: Mr Joachim Kreuzburg (Chairman and Chief Executive Officer of SSB SA and Chief Executive Officer of the Executive Committee of SAG) and Mr René Fáber (Member of the Board of Directors of SSB S.A. and member of the Executive Committee of SAG)
- Nature and purpose: general assistance and administrative services agreement signed on 15 February 2018 with retrospective effect commencing 1 January 2015 for an indefinite duration. This agreement covers the recharging by SAG to SSB S.A. of a part of the remuneration of Mr Joachim Kreuzburg and Mr René Fáber in respect of the services they perform and provide within the company.
- Details: the recharge of the said services of the person concerned is calculated using an allocation based on work performed and time spent by the executives for the benefit of SSB S.A.

The amounts invoiced (excluding tax) by SAG to SSB S.A. for the years ended 31 December 2018, 2019 and 2020 is detailed below:

	Mr Joachim Kreuzburg	Mr René Fáber
Fiscal year 2020	€ 761,917	€ 608,400
Fiscal year 2019	€ 582,804	€ 410,004
Fiscal year 2018	€ 674,216	

Related Party Agreements Already Approved by the General Meeting of Shareholders

We hereby inform you that we have not been advised of any agreements already approved by the general meeting of shareholders and which continued during the previous financial year.

Marseille, 15 February 2021

The Statutory Auditors
French original signed by

KPMG Audit
A division of KPMG S.A.
John Evans
Partner

Deloitte & Associés
Philippe Battisti
Partner