

Sartorius Stedim Biotech

Special Report of the Statutory Auditors on Related Party Agreements

Extract from the Universal Registration Document 2023

6.3 Special Report of the Statutory Auditors on Related Party Agreements

General meeting of Shareholders to approve the financial statements for the year ended 31 December 2023

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.

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 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, on the 9 February 2024

The Statutory Auditors

French original signed by

KPMG Audit
Department of KPMG S.A.

Deloitte & Associés

Nicolas BLASQUEZ

Philippe BATTISTI