

BUREAU VERITAS
Certification



SERA LABORATORIES INTERNATIONAL LIMITED

UNIT 2 SITE J
ALBERT DRIVE
BURGESS HILL
WEST SUSSEX
RH15 9TN

Bureau Veritas Certification Holding SAS – UK Branch certify that the Management System of the above organisation has been audited and found to be in accordance with the requirements of the management system standards detailed below

BS EN ISO 9001:2015

Scope of certification

THE PURCHASE, REPACKAGING AND SUPPLY OF SERUM AND RELATED PRODUCTS.

Original cycle start date: **12 JANUARY 2011**

Expiry date of previous cycle: **04 FEBRUARY 2017**

Recertification Audit date: **26 JANUARY 2017**

Recertification cycle start date: **24 FEBRUARY 2017**

Subject to the continued satisfactory operation of the organisation's Management System, this certificate expires on: **04 FEBRUARY 2020**

Certificate No. UK008493 Version: 1 Revision Date 24 FEBRUARY 2017



*Certification body address: 5th Floor, 66 Prescot Street, London E1 8HG, United Kingdom.
Local office: 5th Floor, 66 Prescot Street, London E1 8HG, United Kingdom.*

0008

*Further clarifications regarding the scope of this certificate and the applicability of the management system requirements may be obtained by consulting the organisation.
To check this certificate validity please call: +44 (0) 207 550 8998*

