




Authorised Economic Operators (AEO) Certificate

Certificate number

GB AEO 00350/19

| | |
|--|---|
| <p>1 Holder of the AEO Certificate</p> <p>Name: Sera Laboratories International Limited</p> <p>Eori No: GB 702816456000</p> <p>VAT ID: GB 702816456</p> <p>Legal Registration Number: 03572986</p> | <p>2 Issuing authority</p> <p>UK AEO Central Site HMRC</p> <p><i>[Signature]</i></p> <p> HM Revenue & Customs</p> |
|--|---|

The holder referred to at 1 above is an:

Authorised Economic Operator

Customs simplifications

Security and safety

Customs simplifications/security and safety

3 Date the certificate is effective from

2019-08-25