

AlphaLISA®

**SARS-CoV2 Spike S1 Protein Detection Kit****Product number:** AL3142 C **Lot Number:** 3315532**Material provided:****Kit Format:** AL3142C: 500 assay points AL3142F: 5000 assay points

The number of assay points is based on an assay volume of 20 µL in 96- or 384-well assay plates using components at the recommended concentrations.

**Analyte Format:** AL3142S: 10 µL at 0.25 mg/mL (one vial contains an amount sufficient for performing 10 AlphaLISA standard curves)**Manufacturing date:** 08/02/24 **Document version:** 1**Product Information****Application:** This kit is designed for the quantitative determination of SARS-Cov2 Spike S1 protein using a homogeneous no wash AlphaLISA assay.**Kit contents:** The kit contains six components: AlphaLISA Acceptor beads coated with anti-Spike S1 Antibody, Streptavidin-coated Donor beads, Biotinylated anti-Spike S1, frozen Spike S1 standard, 10X AlphaLISA Immunoassay Buffer, and 5X AlphaLISA Lysis Buffer.**Sensitivity:** Lower Detection Limit (LDL): 34 pg/mL  
Lower Limit of Quantification (LLOQ): 104 pg/mL  
EC<sub>50</sub>: 24.7 ng/mL**Dynamic Range:** 24.7 – 100 000 pg/mL**Storage:** Store kit in the dark at 4 °C, Store frozen analyte at -20 °C. The aliquoted analyte at -20 °C is stable up to 90 days. Avoid freeze-thaw cycles.**Stability:** This kit is stable for at least 6 months from the date of manufacture when stored in its original packaging and the recommended storage conditions.

## Quality Control

Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals, EC<sub>50</sub> and LDL were measured on the EnVision Multilabel Plate Reader with Alpha option. We certify that these results meet our quality release criteria. Maximum counts may vary between bead lots and the instrument used, with no impact on LDL measurement.

Maximum signal: 86204 counts

Minimum signal: 502 counts

EC50: 43.860 ng/mL

LDL: 0.004 pg/mL

LLOQ: 0.025 pg/mL

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

AL3142-R Rev01

The Revvity logo consists of the word "revvity" in a lowercase, sans-serif font. The letters are black and have a modern, clean appearance.