

AlphaLISA®

**Immunodeficiency Virus type-1 p24 protein (HIV p24) (high sensitivity)****Detection Kit****Product number:** AL291 C **Lot Number:** 3196824**Material provided:****Kit Format:** AL291C: 500 assay points AL291F: 5000 assay points

The number of assay points is based on an assay volume of 50 µL in 96- or 384-well assay plates using components at the recommended concentrations.

**Analyte Format:** AL291S: 0.1 µg (one vial contains an amount of human HIV p24 sufficient for performing 10 AlphaLISA standard curves)**Manufacturing date:** September 19, 2023 **Document version:** 1**Product Information****Application:** This kit is designed for the quantitative determination of HIV p24 in serum, buffered solution or cell culture medium using a homogeneous AlphaLISA assay (no wash steps).**Kit contents:** The kit contains 5 components: AlphaLISA Acceptor beads coated with an Anti-Analyte Antibody, Streptavidin-coated Donor beads, Biotinylated Anti-Analyte Antibody, lyophilized analyte and 10X AlphaLISA Immunoassay Buffer. . ICD solutions are **not** provided with this kit.**Sensitivity:** Lower Detection Limit (LDL): 1.8 pg/mL**Dynamic Range:** 1.8 – 30 000 pg/mL**Storage:** Store kit in the dark at +4°C. Store reconstituted analyte at -20°C.**Stability:** This product is stable for at least 12 months from the manufacturing date when stored in its original packaging and the recommended storage conditions. Note: Once reconstituted, the HIV p24 analyte is stable for at least 75 days at -20°C**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: 299,932 counts

Minimum signal: 294 counts

EC<sub>50</sub>: 3.826 ng/mL

LDL: 0.657 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.

AL291-R Rev01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000 [www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.