

AlphaLISA®

High Performance (HP) Mouse IL1 β Detection Kit**Product number:** AL592C **Lot Number:** 3337325**Material provided:****Kit Format:**

AL592HV: 100 assay points

The number of assay points based on an assay volume of 100 μ L in a 96-well assay plate using kit components at the recommended concentrations.

AL592C: 500 assay points AL592F: 5000 assay points

The number of assay points is based on an assay volume of 50 μ L in 384-well assay plates using kit components at the recommended concentrations.

Analyte Format:AL592S: 0.1 μ g (one vial contains an amount sufficient for performing 10 AlphaLISA standard curves)**Manufacturing date:**

September 16, 2024

Document version:

1

Product Information**Application:**

This kit is designed for the quantitative determination of mouse IL1 β using a homogeneous no wash AlphaLISA assay. This assay is also compatible with rat IL1 β but requires the use of a recombinant rat standard (not provided in the kit) to generate a standard curve.

Kit contents:

The kit contains 5 components: AlphaLISA Anti-mIL1 β Acceptor beads, Streptavidin-coated Donor beads, Biotinylated Anti- mIL1 β Antibody, Lyophilized mIL1 β and 10X AlphaLISA Immunoassay Buffer.

Sensitivity:

Lower Detection Limit (LDL): 0.79 pg/mL;

Lower Limit of Quantification (LLOQ): 2.60 pg/mL

EC₅₀: 9.70 ng/mL**Dynamic Range:**

0.79– 30 000 pg/mL

Storage:

Store kit in the dark at +4°C. For reconstituted analyte, aliquot and store at -20 °C. The aliquoted analyte at -20 °C is stable up to 28 days. Avoid freeze-thaw cycles.

Stability:

This kit is stable for at least 24 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, EC₅₀ 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.

| | |
|--------------------|---------------|
| EC ₅₀ : | 6.487 ng/mL |
| LDL: | 4.094 pg/mL |
| LLOQ: | 11.220 pg/mL |
| Min counts: | 231 counts |
| Max counts: | 318031 counts |

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.

AL592-R Rev01

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