

AlphaLISA®

High Performance Human Tumor Necrosis Factor alpha (TNF α) Biotin-Free Detection Kit

Product number: AL3158 C **Lot Number:** 3337330

Material provided:

Kit Format:

AL3158HV: 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.

AL3158C: 500 assay points AL3158F: 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: AL3158S: 0.1 μ g (one vial contains an amount sufficient for performing 10 AlphaLISA standard curves)

Manufacturing date: August 7, 2024 **Document version:** 1

Product Information

- Application:** This kit is designed for the quantitative determination of human TNF α using a homogeneous no wash AlphaLISA assay.
- Kit contents:** The kit contains 5 components: AlphaLISA Acceptor beads coated with Anti-hTNF α Antibody, Anti-Digoxigenin Fab Fragment Donor beads, Digoxigenin Anti-hTNF α Antibody, Lyophilized human TNF α standard analyte and 10X AlphaLISA Immunoassay Buffer.
- Sensitivity:** Lower Detection Limit (LDL): 7.88 pg/mL
Lower Limit of Quantification (LLOQ): 31.12 pg/mL
EC₅₀: 57.02 ng/mL
- Dynamic Range:** 7.88 – 100 000 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 : | 40.333 ng/mL |
| LDL: | 31.149 pg/mL |
| LLOQ: | 92.232 pg/mL |
| Min counts: | 293 counts |
| Max counts: | 73920 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.

AL3158-R Rev01



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

(800) 762-4000 www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.