

AlphaLISA®

Human TNF α and TNFRII Binding Kit

Product number: AL3121C **Lot Number:** 3240902

Material provided:

Kit Format: AL3121C: 500 assay points AL3121F: 5000 assay points

The number of assay points is based on an assay volume of 20 μ L in 384-well assay plates using kit components at the recommended concentrations.

Manufacturing date: 12/7/2023 **Document version:** 1

Product Information

- Application:** This kit is designed to assess inhibitors of human TNF α and TNFRII binding, using a homogeneous AlphaLISA assay (no wash steps). This assay can facilitate the design and development of antibody therapeutics by using competitive binding to human TNF α and TNFRII binding
- Kit contents:** The kit contains 5 components: anti-6xHis AlphaLISA Acceptor beads, Streptavidin-coated Donor beads, Biotinylated human TNF α , His tagged human TNFRII and AlphaLISA Immunoassay buffer.
- Storage:** The kit components must be stored at 4 $^{\circ}$ C in the dark. Reconstituted proteins can be stored at –20 $^{\circ}$ C for 3 months.
- Stability:** This kit is stable for at least 6 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

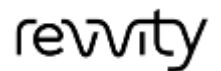
Quality Control

Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals and IC50 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.

Maximum signal: 244477 counts
Minimum signal: 135 counts
IC50: 0.95 nM
SB: 181

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.

AL3121-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.