

AlphaPlex™-545

Human Interferon-gamma (IFN- γ) Detection Kit

Product number: AP217TB-HV **Lot Number:** 3274021

Material provided:

Kit Format:

AP217TB-HV: 100 assay points

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

AP217TB-C: 500 assay points AP217TB-F: 5000 assay points

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

Analyte Format:AL217S: 0.3 μ g (one vial contains an amount of human IFN- γ sufficient for performing 10 standard curves)**Manufacturing date:**

March 13, 2024

Document version:

1

Product Information

Application:

This kit is designed for the quantitative determination of human IFN- γ in serum, buffered solution or cell culture medium using a homogenous AlphaPlex 545 assay (no wash steps).

Kit contents:

The kit contains 5 components: AlphaPlex 545 Acceptor beads coated with an Anti-Analyte Antibody, Streptavidin-coated Donor beads, Biotinylated Anti-Analyte Antibody, lyophilized analyte and 10X AlphaLISA HiBlock Buffer.

Sensitivity:

Lower Detection Limit (LDL): 12 pg/mL

Lower Limit of Quantification (LLOQ): 48 pg/mL

EC₅₀: 50 \pm 30 ng/mL**Dynamic Range:**

10 - 300000 pg/mL

Storage:

Store kit in the dark at +4°C. Store reconstituted analyte at -20°C.

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. Note: Once reconstituted, the human Interferon-gamma (IFN- γ) analyte is stable for at least 3 months when stored at -20°C.

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.

Maximum signal: 83753 counts

Minimum signal: 124 counts

EC50: 115.600 ng/mL

LDL: 0.029 pg/mL

LLOQ: 0.098 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.

AP217TB-R Rev01

The Revvity logo is displayed in a lowercase, sans-serif font.

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.