

AlphaLISA®

Bovine Interferon Gamma (bIFN γ) Detection Kit

Product number: AL535 HV **Lot Number:** 3335650

Material provided:

Kit Format: AL535HV: 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.

AL535C: 500 assay points AL535F: 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: AL535S: 0.3 μ g (one vial contains an amount sufficient for performing 10 AlphaLISA standard curves)

Manufacturing date: November 12, 2024 **Document version:** 1

Product Information

Application: This kit is designed for the quantitative determination of bovine IFN-gamma (bIFN γ) in buffer and cell culture media using a homogeneous AlphaLISA assay (no wash steps). The assay shows negligible cross-reactivity with other bovine cytokines.

Kit contents: The kit contains 5 components: AlphaLISA Anti- bIFN γ Acceptor beads, Streptavidin-coated Donor beads, Biotinylated Anti- bIFN γ Antibody, Lyophilized bTNF γ and 10X AlphaLISA Bovine Immunoassay Buffer.

Sensitivity: Lower Detection Limit (LDL): 10 pg/mL
 Lower Limit of Quantification (LLOQ): 37 pg/mL
 EC₅₀: 9.43 \pm 6.5 ng/mL

Dynamic Range: 1 - 300000 pg/mL

Storage: Store kit in the dark at +4C. 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 bovine IFN γ analyte is stable for at least 18 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.

EC ₅₀ :	9.498 ng/mL
LDL:	0.040 pg/mL
LLOQ:	0.127 pg/mL
Min counts:	1082 counts
Max counts:	77744 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.

AL535-R Rev01

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

