



# AlphaLISA<sup>®</sup> Human Interferon-gamma (IFN- $\gamma$ ) Biotin-Free Detection Kit

**Product number:** AL327 HV/C/F

Research Use Only. Not for use in diagnostic procedures.

## Product Information

**Application:** This kit is designed for the quantitative determination of Human Interferon-gamma (IFN- $\gamma$ ) in cell culture supernatants using a homogeneous AlphaLISA assay (no wash steps).

**Sensitivity:** Lower Detection Limit (LDL): 2.4 pg/mL  
Lower Limit of Quantification (LLOQ): 4.2 pg/mL  
EC<sub>50</sub>: 4.8 ng/mL

**Dynamic range:** 2 -10 000 p/mL (Figure 1).

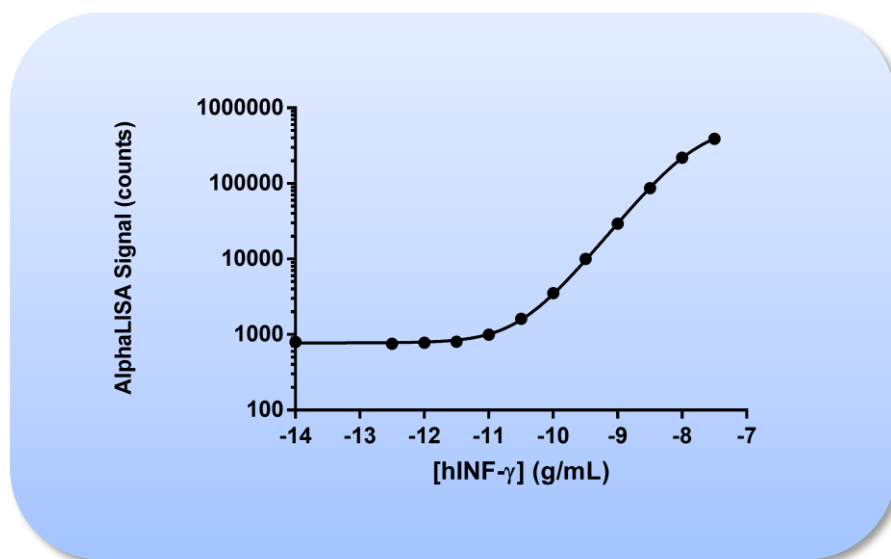


Figure 1. Typical sensitivity curves in AlphaLISA Immunoassay Buffer. The data was generated using a white Optiplate<sup>™</sup>-384 microplate and the EnVision<sup>®</sup> Multilabel Plate Reader with Alpha option 2103.

**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 hINF- $\gamma$  analyte is stable for at least 18 months when stored at -20°C.

## Analyte of Interest

Interferons (IFNs) activity has been discovered due to their antiviral effects. In humans, there are three families of IFNs: IFN type I (IFN- $\alpha$ ,  $\beta$ ,  $\omega$ ,  $\epsilon$ , and  $\kappa$ ), IFN type II (one single representative, IFN- $\gamma$ ), and IFN type III (IFN- $\lambda$ 1-3). Antigens and mitogens stimulate in Natural Killer (NK) and activated helper T lymphocytes (Th1) the production of IFN- $\gamma$ . Human IFN- $\gamma$  is a 140 amino acids polypeptide that shows multiple effects; it induces the production of cytokines, upregulates the expression of class I and II MHC antigens, and leukocyte adhesion molecules. It also activates macrophages, enhances the secretion of immunoglobulins by B cells, and potentiates Th1 cell expansion. Response to IFN- $\gamma$  is mediated by the heterodimeric IFN- $\gamma$  Receptor, triggering a signaling cascade involving JAK1, JAK2, and STAT1. Importantly, IFNs have been proved to be effective in the treatment of several viral infections and cancers.

## Description of the AlphaLISA Assay

AlphaLISA technology allows the detection of molecules of interest in buffer, cell culture media, serum and plasma in a highly sensitive, quantitative, reproducible and user-friendly mode. In this AlphaLISA assay, a DIG-labeled Anti-Analyte Antibody binds to the anti-DIG Alpha Donor beads, while another Anti-Analyte Antibody is conjugated to AlphaLISA Acceptor beads. In the presence of the analyte, the beads come into close proximity. The excitation of the Donor beads provokes the release of singlet oxygen molecules that triggers a cascade of energy transfer in the Acceptor beads, resulting in a sharp peak of light emission at 615 nm (Figure 2).

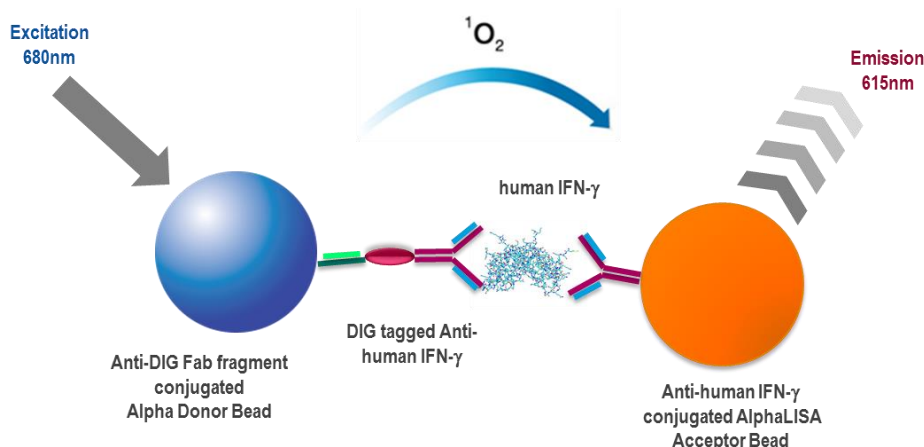


Figure 2. AlphaLISA Assay principle

## Precautions

- Anti-Digoxigenin Fab Fragment Donor beads are light-sensitive. All the other assay reagents can be used under normal light conditions. All Alpha assays using the Donor beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (LEE 090 filters (preferred) or Roscolux filters #389 from Rosco) can be applied to light fixtures.
- All blood components and biological materials should be handled as potentially hazardous. The analyte included in this kit is from a Mouse source.
- Some analytes are present in saliva. Take precautionary measures to avoid contamination of the reagent solutions.
- The DIG labeled anti-analyte antibody can be toxic. Contact with skin or inhalation should be avoided.

## Kit Content: Reagents and Materials

Kit components	AL327HV (100 assay points <sup>***</sup> )	AL327C (500 assay points <sup>***</sup> )	AL327F (5000 assay points <sup>***</sup> )
AlphaLISA Anti-hIFN- $\gamma$ Acceptor beads stored in PBS, 0.05% Proclin-300, pH 7.2	25 $\mu$ L @ 5 mg/mL (1 brown tube, <u>white</u> cap)	50 $\mu$ L @ 5 mg/mL (1 brown tube, <u>white</u> cap)	500 $\mu$ L @ 5 mg/mL (1 brown tube, <u>white</u> cap)
Anti-Digoxigenin Fab Fragment Donor beads stored in 25 mM HEPES, 100 mM NaCl, 0.05% Proclin-300, pH 7.4	100 $\mu$ L @ 5 mg/mL (1 brown tube, <u>black</u> cap)	200 $\mu$ L @ 5 mg/mL (1 brown tube, <u>black</u> cap)	2 mL @ 5 mg/mL (1 brown tube, <u>black</u> cap)
DIG labeled Anti-hIFN- $\gamma$ stored in PBS, 0.1% Tween-20, 0.05% NaN <sub>3</sub> , pH 7.4	25 $\mu$ L @ 500 nM (1 tube, <u>black</u> cap)	50 $\mu$ L @ 500 nM (1 tube, <u>black</u> cap)	500 $\mu$ L @ 500 nM (1 tube, <u>black</u> cap)
AlphaLISA hIFN- $\gamma$ (0.3 $\mu$ g), lyophilized analyte	1 tube, <u>clear</u> cap	1 tube, <u>clear</u> cap	1 tube, <u>clear</u> cap
AlphaLISA Immunoassay Buffer (10X)	2 mL, 1 small bottle	10 mL, 1 small bottle	100 mL, 1 large bottle

\* Reconstitute Human IFN- $\gamma$  in 100  $\mu$ L Milli-Q<sup>®</sup> grade H<sub>2</sub>O. The reconstituted analyte should be used within 60 minutes or aliquoted into screw-capped polypropylene vials and stored at -20°C for further experiments. Avoid multiple freeze-thaw cycles. It has been demonstrated that reconstituted h is stable for at least 18 months at -20°C. One vial contains an amount of human IFN- $\gamma$  sufficient for performing 10 standard curves. Additional vials can be ordered separately (cat # AL217S).

\*\* Extra buffer can be ordered separately (cat # AL000C: 10 mL, cat # AL000F: 100 mL).

\*\*\* The number of assay points is based on an assay volume of 100  $\mu$ L in 96-well plates (AL327HV) or 50  $\mu$ L in 96- or 384-well assay plates using the kit components at the recommended concentrations.

Sodium azide should **not** be added to the stock reagents. High concentrations of sodium azide (> 0.001 % final in the assay) might decrease the AlphaLISA signal.

### Specific additional required reagents and materials:

The following materials are recommended:

Item	Suggested source	Catalog #
TopSeal™-A Adhesive Sealing Film	Revvity Inc.	6050195
EnVision®-Alpha Reader	Revvity Inc.	-

## Recommendations

### General recommendations:

- The volume indicated on each tube is guaranteed for single pipetting. Multiple pipetting of the reagents may reduce the theoretical amount left in the tube. To minimize loss when pipetting beads, it is preferable not to pre-wet the tip.
- Centrifuge all tubes (including lyophilized analyte) before use to improve recovery of content (2000g, 10-15 sec). Re-suspend all reagents by vortexing before use.
- Use Milli-Q® grade H<sub>2</sub>O (18 MΩ•cm) to dilute 10X AlphaLISA Immunoassay Buffer to reconstitute the lyophilized analyte.
- When diluting the standard or samples, change tips between each standard or sample dilution. When loading reagents in the assay microplate, change tips between each standard or sample addition and after each set of reagents.
- When reagents are added to the microplate, make sure the liquids are at the bottom of the well.
- Small volumes may be prone to evaporation. It is recommended to cover microplates with TopSeal-A Adhesive Sealing Films to reduce evaporation during incubation. Microplates can be read with the TopSeal-A Film.
- The AlphaLISA signal is detected with an EnVision Multilabel Reader equipped with the Alpha option using the AlphaScreen standard settings (e.g. Total Measurement Time: 550 ms, Laser 680 nm Excitation Time: 180 ms, Mirror: D640as, Emission Filter: M570w, Center Wavelength 570 nm, Bandwidth 100 nm, Transmittance 75%).
- AlphaLISA signal will vary with temperature and incubation time. For consistent results, identical incubation times and temperature should be used for each plate.
- The standard curves shown in this technical data sheet are provided for information only. A standard curve must be generated for each experiment. The standard curve should be performed in the Immunoassay buffer for serum and/or plasma samples.

## Assay Procedure

IMPORTANT: PLEASE READ THE RECOMMENDATIONS BELOW BEFORE USE

- The manual described below is an example for generating one standard curve in a 50 µL final assay volume (48 wells, triplicate determinations). The manuals also include testing samples in 354 wells. If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly, as shown in the table below. These calculations do not include excess reagent to account for losses during transfer of solutions or dead volumes.
- The standard dilution manual is provided for information only. As needed, the number of replicates or the range of concentrations covered can be modified.
- Use of four background points in triplicate (12 wells) is recommended when LDL/LLOQ is calculated. One background point in triplicate (3 wells) can be used when LDL/LLOQ is not calculated.

Format	# of data points	Volume				Plate recommendation
		Final	Sample	AlphaLISA beads / Anti-DIG Antibody MIX	Donor beads	
AL327HV	100	100 µL	10 µL	40 µL	50 µL	White OptiPlate-96 (cat # 6005290) White ½ AreaPlate-96 (cat # 6005560)
AL327C	250	100 µL	10 µL	40 µL	50 µL	White OptiPlate-96 (cat # 6005290) White ½ AreaPlate-96 (cat # 6005560)
	500	50 µL	5 µL	20 µL	25 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290) Light gray AlphaPlate™-384 (cat # 6005350)
	1 250	20 µL	2 µL	8 µL	10 µL	Light gray AlphaPlate-384 (cat # 6005350) ProxiPlate™-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	2 500	10 µL	1 µL	4 µL	5 µL	Light gray AlphaPlate-1536 (cat # 6004350)
AL327F	5 000	50 µL	5 µL	20 µL	25 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290) Light gray AlphaPlate-384 (cat # 6005350)
	12 500	20 µL	2 µL	8 µL	10 µL	Light gray AlphaPlate-384 (cat # 6005350) ProxiPlate-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	25 000	10 µL	1 µL	4 µL	5 µL	Light gray AlphaPlate-1536 (cat # 6004350)

The manual described below is for 500 assay points including one standard curve (48 wells) and samples (452 wells).

If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly.

- 1) Preparation of 1X AlphaLISA Immunoassay Buffer:  
Add 3 mL of 10X AlphaLISA Immunoassay Buffer to 27 mL H<sub>2</sub>O.
- 2) Preparation of Anti- hIFN- $\gamma$  analyte standard dilutions:
  - a) Reconstitute lyophilized hIFN- $\gamma$  (0.3  $\mu$ g) in 100  $\mu$ L of H<sub>2</sub>O.
  - b) Prepare standard dilutions as follows in 1X AlphaLISA Immunoassay Buffer (change tip between each standard dilution):

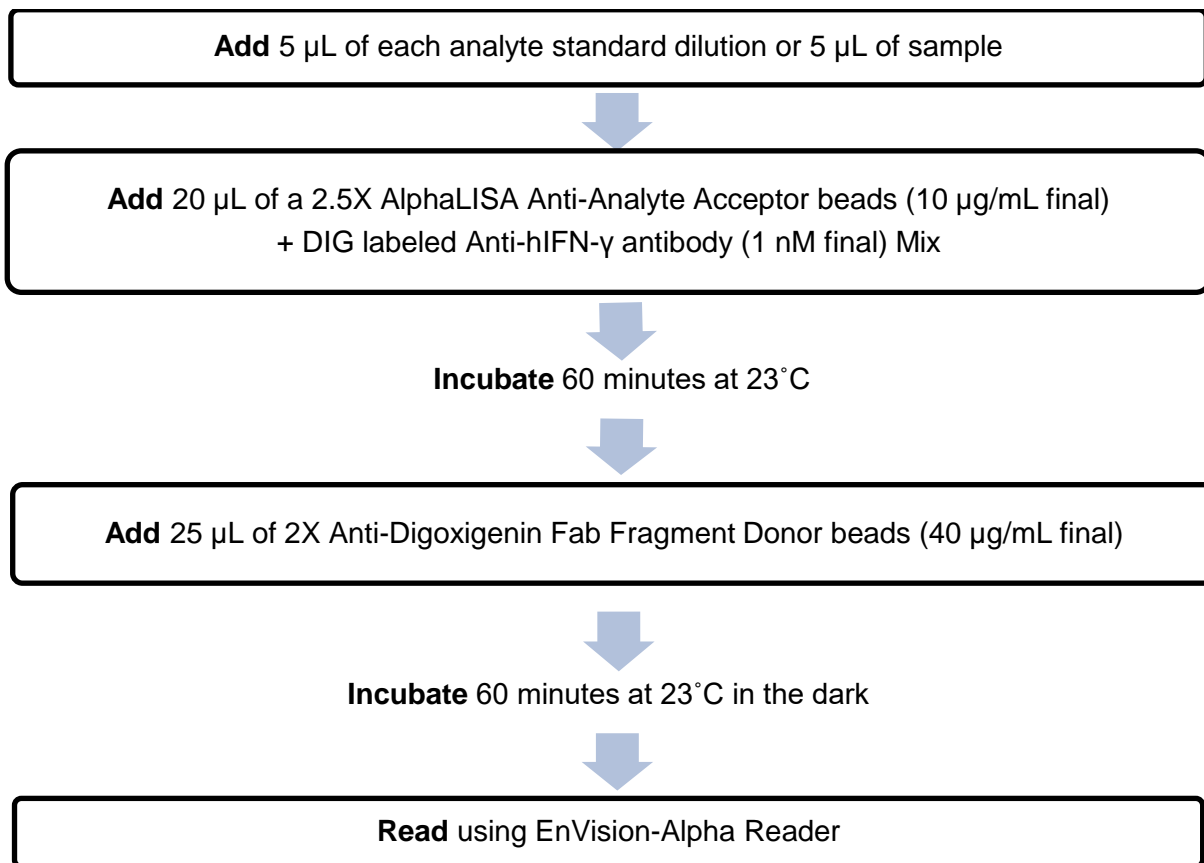
Tube	Vol. of hIFN- $\gamma$ ( $\mu$ L)	Vol. of diluent ( $\mu$ L) *	[hIFN- $\gamma$ ] in standard curve		Final [hIFN- $\gamma$ ] in well
			(g/mL in 5 $\mu$ L)	(pg/mL in 5 $\mu$ L)	(g/mL in 50 $\mu$ L)
A	10 $\mu$ L of provided hIFN- $\gamma$	90	3.00E-07	300000	3.00E-08
B	60 $\mu$ L of tube A	120	1.00E-07	100000	1.00E-08
C	60 $\mu$ L of tube B	140	3.00E-08	30000	3.00E-09
D	60 $\mu$ L of tube C	120	1.00E-08	10000	1.00E-09
E	60 $\mu$ L of tube D	140	3.00E-09	3000	3.00E-10
F	60 $\mu$ L of tube E	120	1.00E-09	1000	1.00E-10
G	60 $\mu$ L of tube F	140	3.00E-10	300	3.00E-11
H	60 $\mu$ L of tube G	120	1.00E-10	100	1.00E-11
I	60 $\mu$ L of tube H	140	3.00E-11	30	3.00E-12
J	60 $\mu$ L of tube I	120	1.00E-11	10	1.00E-12
K	60 $\mu$ L of tube J	140	3.00E-12	3	3.00E-13
L	60 $\mu$ L of tube K	120	1.00E-12	1	1.00E-13
M ** (background)	0	100	0	0	0
N ** (background)	0	100	0	0	0
O ** (background)	0	100	0	0	0
P ** (background)	0	100	0	0	0

\* Dilute standards in diluent (e.g. 1X AlphaLISA Immunoassay Buffer).  
At low concentrations of analyte, a significant amount of analyte can bind to the vial. Therefore, load the analyte standard dilutions in the assay microplate within 60 minutes of preparation.

\*\* Four background points in triplicate (12 wells) are used when LDL is calculated. If LDL does not need to be calculated, one background point in triplicate can be used (3 wells).

- 3) Preparation of 2.5X AlphaLISA Anti-hIFN- $\gamma$  Acceptor beads (25  $\mu$ g/mL) + DIG labeled Anti-hIFN- $\gamma$  Antibody ( 2.5nM) Mix:
  - a. Add 50  $\mu$ L of 5 mg/mL AlphaLISA Anti hIFN- $\gamma$  Acceptor beads and 50  $\mu$ L of 500nM Anti hIFN- $\gamma$  Antibody to 9900  $\mu$ L of 1X AlphaLISA Immunoassay Buffer.
  - b. Prepare just before use.
- 4) Preparation of 2X Anti-Digoxigenin Fab Fragment Donor beads (80  $\mu$ g/mL):
  - a. Keep the beads under subdued laboratory lighting.
  - b. Add 200  $\mu$ L of 5 mg/mL Anti-Digoxigenin Fab Fragment Donor beads to 12300  $\mu$ L of 1X AlphaLISA Immunoassay Buffer.
  - c. Prepare just before use.

5) In a white Optiplate (384 wells):



## Data Analysis

- Calculate the average count value for the background wells.
- Generate a standard curve by plotting the AlphaLISA counts versus the concentration of analyte. A log scale can be used for either or both axes. No additional data transformation is required.
- Analyze data according to a nonlinear regression using the 4-parameter logistic equation (sigmoidal dose-response curve with variable slope) and a  $1/Y^2$  data weighting (the values at maximal concentrations of analyte after the hook point should be removed for correct analysis).
- The LDL is calculated by interpolating the average background counts (12 wells without analyte) + 3 x standard deviation value (average background counts + (3xSD)) on the standard curve.
- The LLOQ as measured here is calculated by interpolating the average background counts (12 wells without analyte) + 10 x standard deviation value (average background counts + (10xSD)) on the standard curve. Alternatively, the true LLOQ can be determined by spiking known concentrations of analyte in the matrix and measuring the percent recovery, and then determining the minimal amount of spiked analyte that can be quantified within a given limit (usually +/- 20% or 30% of the real concentration).
- Read from the standard curve the concentration of analyte contained in the samples.
- If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.

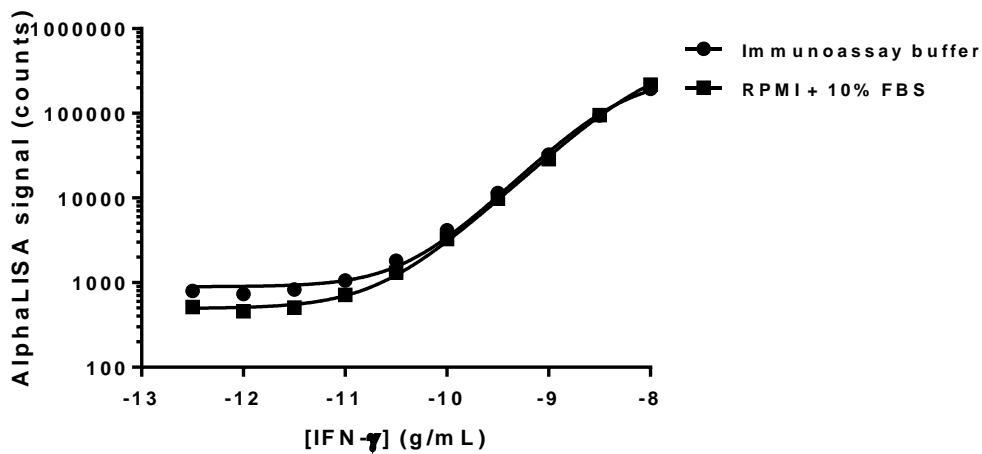
## Assay Performance Characteristics

AlphaLISA assay performance described below was determined using the 2 step manual.  
**When using cell culture media it is recommended to use 10% FBS.**

- Assay Sensitivity:

The LDL was calculated as described above. The values correspond to the lowest concentration of analyte that can be detected in a volume of 5  $\mu$ L using the recommended assay conditions.

LDL (pg/mL)	Buffer/Media	# of experiments
2.4	Immunoassay Buffer	12
3	RPMI with 10% FBS	6



\* Note that LDL can be decreased (i.e. sensitivity increased) by increasing the volume of analyte in the assay (e.g. use 10  $\mu$ L of analyte in a final assay volume of 50  $\mu$ L).

- Assay Precision:

The following assay precision data were calculated from the three independent assays using two different kit lots. In each lot, the analytes were prepared in AlphaLISA Immunoassay Buffer, or RPMI supplemented with 10% FBS. Each assay consisted of one standard curve comprising 12 data points (each in triplicate) and 12 background wells (no analytes). The assays were performed in 384-well format using AlphaLISA Immunoassay Buffer.

- Intra-assay precision:

The intra-assay precision was determined using a total of 16 independent determinations in triplicate. Shown is CV%.

hIFN- $\gamma$	Immunoassay Buffer	RPMI
CV%	10%	12%



- Inter-assay precision:

The inter-assay precision was determined using a total of 3 independent determinations with 9 measurements for 4ng/mL sample. Shown are CV%.

hIFN- $\gamma$ (1 ng/mL)	Immunoassay Buffer	RPMI
CV%	10%	13%

- Spike Recovery:

Four known concentrations of analyte were spiked in Immunoassay Buffer and cell culture media containing 10% FBS. All samples, including non-spiked Immunoassay Buffers and culture media were measured in the assay. The average recovery from three independent measurements is reported.

Spiked hIFN- $\gamma$ (ng/mL)	% Recovery	
	Immunoassay Buffer	RPMI
10	75%	102%
1	117%	121%
0.1	128%	130%

- Specificity:

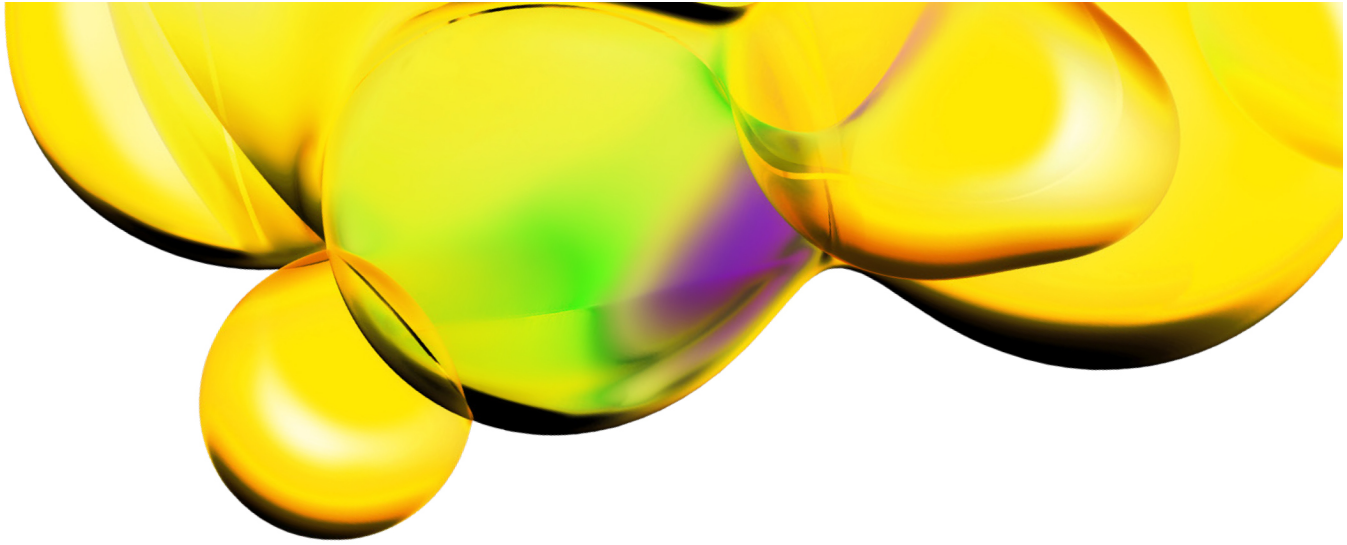
Cross-reactivity of the AlphaLISA Anti hIFN- $\gamma$  Kit was tested using the following proteins at 0.3  $\mu$ g/mL in AlphaLISA Immunoassay Buffer.

Protein	% Cross-reactivity
Mouse IFN- $\gamma$	0
Rat IFN- $\gamma$	0
Bovine IFN- $\gamma$	0
Rhesus macaque IFN- $\gamma$	0

## Troubleshooting Guide

You will find detailed recommendations for common situations you might encounter with your AlphaLISA Assay kit at: [www.revivity.com](http://www.revivity.com)

**FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. 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

[www.revvity.com](http://www.revvity.com)

revvity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.