



## **IFN- $\gamma$ (human), ELISA kit**

**Catalog No. ADI-900-136**

**96 Well Kit**

### Table of Contents

Description	Page	2
Introduction		2
Precautions		2
Materials Supplied		3
Storage		3
Materials Needed but Not Supplied		3
Sample Handling		4
Procedural Notes		4
Reagent Preparation		5
Assay Procedure		6
Calculation of Results		7
Typical Results		7
Typical Standard Curve		8
Calibration		8
Performance Characteristics		9
Sample Dilution Recommendations		9
References		10
Limited Warranty		12

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

## **Description**

The IFN- $\gamma$  (human), ELISA kit is a complete kit for the quantitative determination of human IFN- $\gamma$  in biological fluids. Please read the complete kit insert before performing this assay. The kit uses an antibody to human IFN- $\gamma$  immobilized on a microtiter plate to bind the human IFN- $\gamma$  in the standards or sample. A recombinant human IFN- $\gamma$  Standard is provided in the kit. A biotinylated antibody is added to the standards and samples. This labeled antibody binds to the human IFN- $\gamma$  captured on the plate. After a short incubation the excess standard, sample and antibody are washed out and streptavidin conjugated to Horseradish peroxidase is added, which binds to the biotinylated human IFN- $\gamma$  antibody. Excess conjugate is washed out and substrate is added. The substrate reacts with the conjugate bound to the IFN- $\gamma$  captured on the plate. After a short incubation, the enzyme reaction is stopped and the color generated is read at 450 nm. The measured optical density is directly proportional to the concentration of human IFN- $\gamma$  in either standards or samples. For further explanation of the principles and practices of immunoassays please see the excellent books by Chard<sup>1</sup> or Tijssen<sup>2</sup>.

## **Introduction**

Human Interferon-gamma (IFN- $\gamma$ ) is a polypeptide of 166 amino acids including the signal sequence of 23 amino acids. Two active forms of the protein have been found with different glycosylation sites. The first form is glycosylated only at position 25, while the second is glycosylated at positions 25 and 97.<sup>4,8</sup> IFN- $\gamma$  is alternatively known as antigen-induced interferon, immune interferon (IIF), Type-2 interferon, T interferon, mitogen-induced interferon, and pH2-labile interferon.<sup>8</sup> IFN- $\gamma$  plays an important role in immune response. It is secreted by Th1 cells which in turn activates macrophage through the CD40 surface ligand. This secretion also inhibits the development of Th2 cells.<sup>3</sup> Human and mouse IFN- $\gamma$  share only a 40% homology on the amino acid level.<sup>9</sup>

## **Precautions**

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

1. Stop Solution is a 0.18 M sulfuric acid solution. This solution is caustic; care should be taken in use.
2. The activity of the Horseradish peroxidase conjugate is affected by nucleophiles such as azide, cyanide and hydroxylamine.
3. We test this kit's performance with a variety of samples, however it is possible that high levels of interfering substances may cause variation in assay results.
4. The human IFN- $\gamma$  Standard provided, Catalog No. 80-1289, should be handled with care because of the known and unknown effects of IFN- $\gamma$ .

## **Materials Supplied**

- 1. human IFN- $\gamma$  Microtiter Plate, One Plate of 96 Wells, Catalog No. 80-1288**  
A strip microtiter plate coated with antibody specific to human IFN- $\gamma$ .
- 2. human IFN- $\gamma$  Antibody , 8 mL, Catalog No. 80-1291**  
A solution of biotinylated antibody to human IFN- $\gamma$ .
- 3. human IFN- $\gamma$  Standard Diluent, 12 mL, Catalog No. 80-1290**  
Phosphate buffered saline containing proteins and antibiotics.
- 4. human IFN- $\gamma$  Streptavidin-HRP Concentrate, 75  $\mu$ L, Catalog No. 80-1292**  
A concentrated solution of streptavidin conjugated to Horseradish peroxidase.
- 5. human IFN- $\gamma$  Streptavidin-HRP Dilution Buffer, 14 mL, Catalog No. 80-1293**
- 6. Wash Buffer Concentrate, 50 mL, Catalog No. 80-1253**  
Tris buffered saline containing detergents.
- 7. human IFN- $\gamma$  Standard, 2 vials, Catalog No. 80-1289**  
Two vials of lyophilized recombinant human IFN- $\gamma$ .
- 8. human IFN- $\gamma$  TMB Substrate, 13 mL, Catalog No. 80-1294**  
A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide. **Protect from prolonged exposure to light.**
- 9. human IFN- $\gamma$  Stop Solution, 13 mL, Catalog No. 80-1295**  
A 0.18 M solution of sulfuric acid in water. Keep tightly capped. Caution: **Caustic.**
- 10. human IFN- $\gamma$  Assay Layout Sheet, 1 each, Catalog No. 30-0212**
- 11. Plate Sealer, 3 each, Catalog No. 30-0012**

## **Storage**

All components of this kit are stable at 4°C until the kit's expiration date.

## **Materials Needed but Not Supplied**

1. Deionized or distilled water. No difference in assay results is seen with distilled water.
2. Precision pipets for volumes between 50  $\mu$ L and 1,000  $\mu$ L.
3. Repeater pipet for dispensing 50 and 100  $\mu$ L.
4. Disposable beakers for diluting buffer concentrates.
5. Graduated cylinders.
6. Adsorbent paper for blotting.
7. Microcentrifuge to prepare Streptavidin-HRP Solution.
8. Microplate reader capable of reading at 450 nm, preferably with correction between 570 nm and 590 nm.
9. Log-log graph paper for plotting the standard curve.

## **Sample Handling**

The IFN- $\gamma$  (human), ELISA kit is compatible with human IFN- $\gamma$  samples in a wide range of matrices.

Culture fluids, serum, EDTA, heparin and sodium citrate plasma and urine are suitable for use in the assay. Samples containing a visible precipitate must be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic specimens. Samples in the majority of tissue culture media can also be read in the assay, provided the standards have been diluted into the tissue culture media instead of Standard Diluent. There will be a small change in binding associated with running the standards and samples in media. Users should only use standard curves generated in media, or Standard Diluent to calculate concentrations of human IFN- $\gamma$  in the appropriate matrix.

Samples must be stored frozen to avoid loss of bioactive human IFN- $\gamma$ . If samples are to be run within 24 hours, they may be stored at 4°C. Otherwise, samples must be stored frozen at or below -70°C to avoid loss of bioactive human IFN- $\gamma$ . Up to three freeze/thaw cycles of serum has been shown to have no effect on human IFN- $\gamma$  levels. Nonetheless, excessive freeze/thaw cycles should be avoided. Prior to assay, frozen sera should be brought to room temperature slowly and gently mixed by hand. Do not thaw samples in a 37°C incubator. Do not vortex or sharply agitate samples.

## **Procedural Notes**

1. Do not mix components from different kit lots or use reagents beyond the kit expiration date.
2. Allow all reagents to warm to room temperature for at least 30 minutes before opening.
3. Standards must be prepared in polypropylene or polyethylene tubes. Do not use polystyrene, polycarbonate or glass tubes.
4. Pre-rinse the pipet tip with reagent, use fresh pipet tips for each sample, standard and reagent.
5. Pipet standards and samples to the bottom of the wells.
6. Add the reagents to the side of the well to avoid contamination.
7. This kit uses plates with removable strips. Unused strips must be kept desiccated at 4°C in the sealed bag provided. The wells should be used in the frame provided.
8. **Prior to addition of conjugate and substrate, ensure that there is no residual wash buffer in the wells. Any remaining wash buffer may cause variation in assay results.**
9. **It is important that the matrix for the standards and samples be as similar as possible. Human IFN- $\gamma$  samples diluted with Standard Diluent should be run with a standard curve diluted in the same buffer. Serum, plasma and urine samples should be evaluated against a standard curve run in Standard Diluent while Tissue Culture samples should be read against a standard curve diluted in the same complete but non-conditioned media. See Reagent Preparation, step #2.**

## **Reagent Preparation**

### **1. Wash Buffer**

Prepare the Wash Buffer by diluting 50 mL of the supplied concentrate with 1,450 mL of deionized water. This can be stored at room temperature until the kit expiration, or for 3 months, whichever is earlier.

### **2. human IFN- $\gamma$ Standards**

Reconstitute standard with deionized water. Reconstitution volume is stated on the standard vial label. Let it sit at room temperature for 5 minutes. Mix gently. This solution contains 2,500 pg/mL human IFN- $\gamma$ . When testing serum, plasma or urine samples, use the Standard Diluent provided to prepare standard curve serial dilutions. When using cell culture supernatants use tissue culture media to prepare the standard curve serial dilutions.

Label five 12x75 mm test tubes #1 through #5. Pipet 240  $\mu$ L of Standard Diluent or tissue culture media into tubes #1 through #5. Add 160  $\mu$ L of the 2,500 pg/mL Standard to tube #1. Vortex thoroughly. Add 160  $\mu$ L of tube #1 to tube #2 and vortex thoroughly. Add 160  $\mu$ L of tube #2 to #3 and vortex thoroughly. Continue this for tubes #4 and #5.

**The concentration of human IFN- $\gamma$  in tubes #1 through #5 will be 1,000, 400, 160, 64 and 25.6 pg/mL respectively. See human IFN- $\gamma$  Assay Layout Sheet for dilution details.**

**Diluted standards should be used within 60 minutes of preparation. Do not store reconstituted standards.**

### **3. Streptavidin-HRP Solution**

Prepare Streptavidin-HRP solution **immediately before use**. Do not store prepared Streptavidin-HRP solution. Use a plastic tube to prepare Streptavidin-HRP solution. Briefly centrifuge the Streptavidin-HRP Concentrate to force entire vial contents to the bottom. For each strip used, mix 2.5  $\mu$ L of Streptavidin-HRP Concentrate with 1 mL of Streptavidin-HRP Dilution Buffer.

## **Assay Procedure**

**Bring all reagents to room temperature for at least 30 minutes prior to opening.**

**All standards, controls and samples should be run in duplicate.**

1. Refer to the Assay Layout Sheet to determine the number of strips to be used and put any remaining strips with the desiccant back into the pouch and seal the ziploc. Store unused wells at 4°C.
2. Pipet 50  $\mu\text{L}$  of the Biotinylated Antibody into each well, except the Blank.
3. Pipet 50  $\mu\text{L}$  of standard diluent or Tissue Culture Media into the S0 (0 pg/mL standard) wells.
4. Pipet 50  $\mu\text{L}$  of Standards #1 through #5 into the appropriate wells.
5. Pipet 50  $\mu\text{L}$  of the Samples into the appropriate wells.
6. Tap the plate gently to mix the contents, and seal with the plate sealer provided.
7. Incubate the plate at room temperature for 2 hours.
8. Empty the contents of the wells and wash by adding 400  $\mu\text{L}$  of wash solution to every well. Repeat the wash 2 more times for a total of **3 washes**. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
9. Add 100  $\mu\text{L}$  of the freshly prepared Streptavidin-HRP Conjugate to each well, except the Blank.
10. Seal the plate and incubate at room temperature for 30 minutes.
11. Empty the contents of the wells and wash by adding 400  $\mu\text{L}$  of wash solution to every well. Repeat the wash 2 more times for a total of **3 washes**. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
12. Pipet 100  $\mu\text{L}$  of Substrate Solution into each well.
13. Incubate for 30 minutes at room temperature in the dark.
14. Pipet 100  $\mu\text{L}$  Stop Solution to each well.
15. Blank the plate reader against the Blank wells, read the optical density at 450 nm, preferably with correction between 570 and 590 nm. If the plate reader is not able to be blanked against the Blank wells, manually subtract the mean optical density of the Blank wells from all the readings.

## Calculation of Results

Several options are available for the calculation of the concentration of human IFN- $\gamma$  in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic curve fitting program. If data reduction software is not readily available, the concentration of human IFN- $\gamma$  can be calculated as follows:

1. Calculate the average net Optical Density (OD) bound for each standard and sample by subtracting the average Blank OD from the average OD for each standard and sample.

$$\text{Average Net OD} = \text{Average OD} - \text{Average Blank OD}$$

2. Plot the Average Net OD for each standard versus concentration of human IFN- $\gamma$  in each standard. Approximate a straight line through the points. The concentration of h IFN- $\gamma$  in the unknowns can be determined by interpolation.

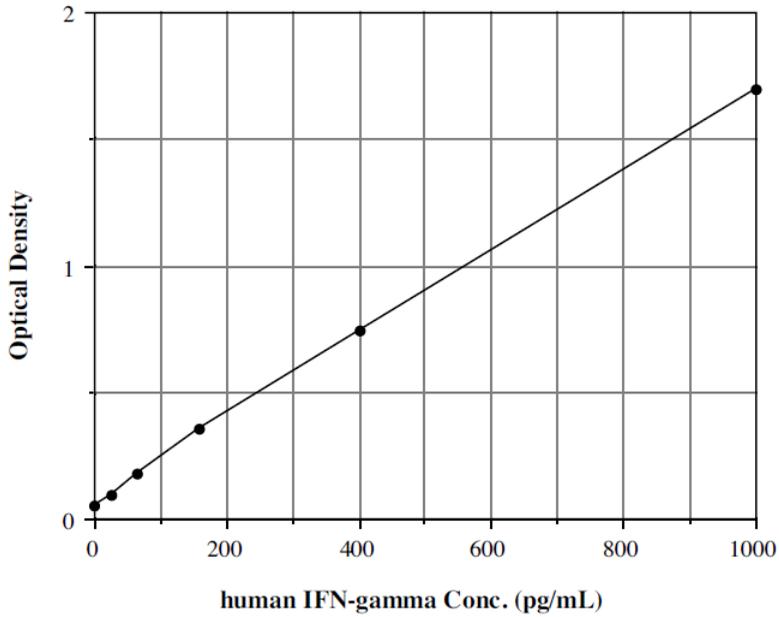
## Typical Results

The results shown below are for illustration only and **should not** be used to calculate results from another assay.

<u>Sample</u>	<u>Average OD</u>	<u>Net OD</u>	<b>h IFN-<math>\gamma</math> (pg/mL)</b>
Blank	(0.043)		
S0	0.097	0.054	<b>0</b>
S1	1.742	1.699	<b>1,000</b>
S2	0.788	0.745	<b>400</b>
S3	0.397	0.354	<b>160</b>
S4	0.221	0.178	<b>64</b>
S5	0.135	0.092	<b>25.6</b>
Unknown 1	1.405	1.362	<b>783.7</b>
Unknown 2	0.307	0.264	<b>113.3</b>

### Typical Standard Curve

A typical standard curve is shown below. This curve **must not** be used to calculate human IFN- $\gamma$  concentrations; each user must run a standard curve for each assay.



### Calibration

The standards in this kit have been calibrated to the NIAID recombinant IFN-g standard lot Gxg23-902-535. One (1) pg of Standard = 0.03 NIAID units.

## **Performance Characteristics**

### **Sensitivity: < 2 pg/mL**

The sensitivity or Lower Limit of Detection (LLD) is determined by assaying replicates of zero and the standard curve. The mean signal of zero + 2 standard deviations read in dose from the standard curve is the LLD. This value is the smallest dose that is not zero with 95% confidence.

### **Linearity**

A sample containing 686.9 pg/mL human IFN- $\gamma$  was serially diluted 4 times 1:2 in the standard diluent supplied in the kit and measured in the assay. The data was plotted graphically as actual human IFN- $\gamma$  concentration versus measured human IFN- $\gamma$  concentration.

The line obtained had a slope of 1.046 with a correlation coefficient of 0.998.

### **Precision**

Intra-assay CV: < 10%

Inter-assay CV: < 10%

### **Cross Reactivities**

The IFN- $\gamma$  (human) ELISA Kit is specific for natural and recombinant human IFN- $\gamma$ . It is unaffected by the presence of human IL-1 $\alpha$ , IL-1 $\beta$ , IL-2, IL-3, IL-6, IL-7, IL-8, TNF- $\alpha$ , or GM-CSF.

### **Sample Recoveries**

Cytokine recovery is determined by spiking low and high levels of recombinant human IFN- $\gamma$  in to normal human serum, plasma, and urine samples, and a Standard Diluent control buffer.

Mean recoveries are as follows:

<b><u>Spike Level</u></b>	<b><u>900 pg/mL</u></b>	<b><u>150 pg/mL</u></b>
Mean Serum Recovery	106%	84%
Mean Plasma Recovery	95%	65%
Mean Urine Recovery	105%	80%

## **References**

1. T. Chard, "An Introduction to Radioimmunoassay & Related Techniques 4th Ed.", (1990) Amsterdam: Elsevier.
2. P. Tijssen, "Practice & Theory of Enzyme Immunoassays", (1985) Amsterdam: Elsevier.
3. Alberts, et al., "The Cell", (2002) New York, Garland Science: 1412-1415.
4. P. Chomarat, et al., "The Cytokine Handbook", (1998) Academic Press: 491-508.
5. E Rinderknecht, et al., J. Biol. Chem., (1984) 259(11): 6790-7.
6. R Devos, et al., Nucleic Acids Res., (1982) 10(8): 2487-501.
7. YK Yip, et al., Proc. Natl. Acad. Sci. USA, (1982) 79(6): 1820-4.
8. Cytokines Online Pathfinder Encyclopaedia, "IFN-gamma".
9. <http://www.copewithcytokines.de/cope.cgi?3172> Accessed 23 July, 2002.
10. P.W. Gray and D.V. Goeddel, Proc. Natl. Acad. Sci. USA, (1983) 80(19): 5842-6.

## Notes



## **USE FOR RESEARCH PURPOSES ONLY**

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products

**[www.enzolifesciences.com](http://www.enzolifesciences.com)**  
***Enabling Discovery in Life Science®***

## **LIMITED WARRANTY; DISCLAIMER OF WARRANTIES**

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

## **TRADEMARKS AND PATENTS**

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

### **Global Headquarters**

#### **Enzo Life Sciences Inc.**

10 Executive Blvd  
Farmingdale, NY 11735  
(p) 1-800-942-0430  
(f) 1-631-694-7501  
(e) [info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

#### **Enzo Life Sciences (ELS) AG**

Industriestrasse 17, Postfach  
CH-4415 Lause / Switzerland  
(p) +41/0 61 926 89 89  
(f) +41/0 61 926 89 79  
(e) [info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)

Please visit our website at [www.enzolifesciences.com](http://www.enzolifesciences.com) for additional contact information.

