



EAGLE
BIOSCIENCES

Rat IFN- γ ELISA KIT

Catalog Number:

IFG21-K01 (1 x 96 wells)

For Research Use Only. Not for use in diagnostic procedures.

v. 1.0

EAGLE BIOSCIENCES, INC.
20A Northwest Blvd., Suite 112, Nashua, NH 03063
Phone: 617-419-2019 Fax: 617-419-1110
WWW.EAGLEBIO.COM



INTENDED USE

The Eagle Biosciences Rat Interferon Gamma (IFN- γ) ELISA Assay Kit (enzyme-linked immunoassay kit) is intended for the quantitative determination of rat γ -interferon (IFN- γ) concentrations in cell culture supernates, serum, and plasma. The Eagle Biosciences Rat Interferon Gamma (IFN- γ) ELISA Assay Kit is for research use only and not to be used in diagnostic procedures.

INTRODUCTION

Interferon gamma (IFN- γ) is a multifunctional protein first observed as an antiviral activity in cultures of Sindbis virus-infected human leukocytes stimulated by PHA (1). Produced by T-lymphocytes and natural killer (NK) cells, interferon gamma (IFN- γ) is now known to be both an inhibitor of viral replication and a regulator of numerous immunological functions. Human interferon gamma (IFN- γ) is reported to be active only on human and non-human primate cells (5). The biochemistry and biological activities of the interferons have been extensively reviewed (2-9).

Rat interferon gamma (IFN- γ) cDNA encodes a 156 amino acid (aa) residue precursor protein with a putative 19 aa residue signal peptide that is cleaved to generate the mature protein which contains two potential N-glycosylation sites (10-12). Rat IFN- γ shares approximately 87% and 39% amino acid sequence identity with mouse interferon gamma (IFN- γ) and human interferon gamma (IFN- γ), respectively. Consistent with their degrees of shared homology, rat

A receptor for interferon gamma (IFN- γ) has been identified and its gene localized to chromosome 6 (13,14). Apparently the product of a single gene, the receptor is a single chain 90 kDa glycoprotein that shows a high degree of species-specific binding of interferon gamma (IFN- γ) (15-18).

Functionally, interferon gamma (IFN- γ) produces a variety of effects. Produced by CD8+, NK, gd, and TH1 T helper cells, interferon gamma (IFN- γ) has documented antiviral, antiprotozoal and immunomodulatory effects on cell proliferation and apoptosis, as well as the stimulation and repression of a variety of genes (19-22) the antiprotozoal activity of interferon gamma (IFN- γ) against *Toxoplasma* and *Chlamydia* is believed to result from indoleamine 2,3-dioxygenase activity, an enzyme induced by interferon gamma (IFN- γ) (23). The immunomodulatory effects of interferon gamma (IFN- γ) are extensive and diverse. In monocyte/macrophages, the activities of interferon gamma (IFN- γ) include: increasing the expression of class I and II MHC antigens; increasing the production of IL-1, platelet-activating factor, H₂O₂, and pterin; protection of monocytes against LAK cell-mediated lysis; down regulation of IL-8 mRNA expression that is up regulated by IL-2; and, with lipopolysaccharide, induction of NO production. Finally, interferon gamma (IFN- γ) has been shown to upregulate ICAM-1, but not E-Selectin or VCAM-1, expression on endothelial cells.

PRINCIPLE OF THE ASSAY

The Eagle Biosciences Rat Interferon Gamma (IFN- γ) ELISA Assay Kit employs the quantitative sandwich enzyme immunoassay technique. A monoclonal antibody specific for interferon gamma (IFN- γ) has been pre-coated onto a microplate. Standards and samples are pipetted into the wells and any interferon gamma (IFN- γ) present is bound by the immobilized antibody. Following incubation unbound samples are removed during a wash step, and then a detection antibody specific for interferon gamma (IFN- γ) is added to the wells and binds to the combination of capture antibody-IFN- γ in sample. Following a wash to remove any unbound



combination, and enzyme conjugate is added to the wells. Following incubation and wash steps a substrate is added. A colored product is formed in proportion to the amount of interferon gamma (IFN- γ) present in the sample. The reaction is terminated by addition of acid and absorbance is measured at 450nm. A standard curve is prepared from seven interferon gamma (IFN- γ) standard dilutions and interferon gamma (IFN- γ) sample concentration determined.

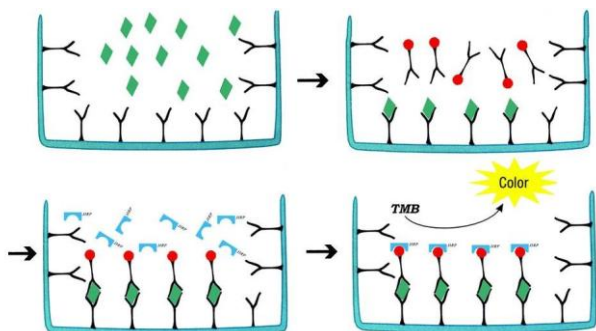


Figure 1:Schematic diagram of the assay

REAGENTS

1. Aluminium pouches with a Microwell Plate coated with monoclonal antibody to rat IFN- γ (8 \times 12)
2. 2 vials human IFN- γ Standard lyophilized, 20000 pg/ml upon reconstitution
3. 2 vials concentrated Biotin-Conjugate anti-human IFN- γ antibody
4. 2 vials Streptavidin-HRP solution,
5. 1 bottle Standard /sample Diluent
6. 1 bottle Biotin-Conjugate antibody Diluent
7. 1 bottle Streptavidin-HRP Diluent
8. 1 bottle Wash Buffer Concentrate 20x (PBS with 1% Tween-20)
9. 1 vial Substrate Solution
10. 1 vial Stop Solution
11. 3 pieces Adhesive Films
12. package insert

NOTE: [96 Tests]



STORAGE

Table 1: Storage of the Rat Interferon Gamma (IFN- γ) ELISA Assay Kit

Unopened Kit	Store at 2 - 8° C. Do not use past kit expiration date.	
Opened/ Reconstituted Reagents	Standard /sample Diluent	May be stored for up to 1 month at 2 - 8° C.**
	Concentrated Biotin-Conjugate	
	Streptavidin-HRP solution	
	Biotin-Conjugate antibody Diluent	
	Streptavidin-HRP Diluent	
	Wash Buffer Concentrate 20x	
	Substrate Solution	
	Stop Solution	
	Standard	Aliquot and store for up to 1 month at $\leq 20^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles. Diluted standard shall not be reused.
	Microplate Wells	Return unused wells to the foil pouch containing the desiccant pack, reseal along entire edge of zip-seal. May be stored for up to 1 month at 2 - 8° C.**

**Provided this is within the expiration date of the kit.

THE REQUIRED ITEMS (not provided):

1. Microplate reader (450nm).
2. Micro-pipette and tips: 0.5-10, 2-20, 20-200, 200-1000ul.
3. 37°C incubator, double-distilled water or deionized water, coordinate paper, graduated cylinder.

PRECAUTIONS FOR USE

1. Store Rat Interferon Gamma (IFN- γ) ELISA Assay Kit reagents between 2° and 8°C. After use all reagents should be immediately returned to cold storage (2° and 8°C).
2. Please perform simple centrifugation to collect the liquid before use.
3. To avoid cross contamination, please use disposable pipette tips.
4. The Stop Solution suggested for use with this Rat Interferon Gamma (IFN- γ) ELISA Assay Kit is an acid solution. Wear eye, hand, face, and clothing protection when using this material. Avoid contact of skin or mucous membranes with Rat Interferon Gamma (IFN- γ) ELISA Assay Kit reagents or specimens. In the case of contact with skin or eyes wash immediately with water.



5. Use clean, dedicated reagent trays for dispensing the washing liquid, conjugate and substrate reagent. Mix all reagents of the Rat Interferon Gamma (IFN- γ) ELISA Assay Kit and samples well before use.
6. After washing microtiter plate should be fully pat dried. Do not use absorbent paper directly into the enzyme reaction wells.
7. Do not mix or substitute reagents of the Rat Interferon Gamma (IFN- γ) ELISA Assay Kit with those from other lots or other sources. Do not use kit reagents beyond expiration date on label.
8. Each sample, standard, blank and optional control samples should be assayed in duplicate or triplicate.
9. Adequate mixing is very important for good result. Use a mini-vortexer at the lowest frequency or Shake by hand at 10min interval when there is no vortexer.
10. Avoid microtiter plates drying during the operation.
11. Dilute samples at the appropriate multiple, and make the sample values fall within the standard curve. If samples generate values higher than the highest standard, dilute the samples and repeat the assay.
12. Any variation in standard diluent, operator, pipetting technique, washing technique, incubation time and temperature, and kit age can cause variation in binding.
13. This Rat Interferon Gamma (IFN- γ) ELISA Assay Kit method can effectively eliminate the interference of the soluble receptors, binding proteins and other factors in biological samples.

SAMPLE COLLECTION AND STORAGE

1. **Cell Culture Supernates** - Remove particulates by centrifugation.
2. **Serum** - Use a serum separator tube (SST) and allow samples to clot for 30 minutes before centrifugation for 15 minutes at approximately 1000 x g. Remove serum, avoid hemolysis and high blood lipid samples.
3. **Plasma** - Recommended EDTA as an anticoagulant in plasma. Centrifuge for 15 minutes at 1000 x g within 30 minutes of collection.
4. Assay immediately or aliquot and store samples at -20°C. Avoid repeated freeze-thaw cycles.
5. Dilute samples at the appropriate multiple (recommended to do pre-test to determine the dilution factor).

Note: Normal rat serum or plasma samples are suggested to make a 1:2 dilution.



REAGENT PREPARATION

1. Bring all reagents of the Rat Interferon Gamma (IFN- γ) ELISA Assay Kit to room temperature before use.
2. **Wash Buffer** - Dilute 10mL of Wash Buffer Concentrate into deionized or distilled water to prepare 200mL of Wash Buffer. If crystals have formed in the concentrate Wash Buffer, warm to room temperature and mix gently until the crystals have completely dissolved.
3. **Standard** - Reconstitute the Standard with 1.0mL of Standard /sample Diluent. This reconstitution produces a stock solution of 20000 pg /mL. Allow the standard to sit for a minimum of 15 minutes with gentle agitation prior to making dilutions.

Pipette 100 μ L of Standard/sample Diluent into the 2000 pg/mL tube and the remaining tubes. Use the stock solution to produce a 2-fold dilution series (below). Mix each tube thoroughly and change pipette tips between each transfer. The 2000 pg/mL standard serves as the high standard. The Standard/ sample Diluent serves as the zero standard (0 pg/mL).

If you do not run out of re-melting standard, store it at -20°C. Diluted standard shall not be reused.

4. Working solution of Biotin-Conjugate anti-rat IFN- γ antibody:
Make a 1:100 dilution of the concentrated Biotin-Conjugate solution with the Biotin-Conjugate antibody Diluent in a clean plastic tube.

The working solution should be used within one day after dilution.

5. Working solution of Streptavidin-HRP: Make a 1:100 dilution of the concentrated Streptavidin-HRP solution with the Streptavidin-HRP Diluent in a clean plastic tube.

The working solution should be used within one day after dilution.

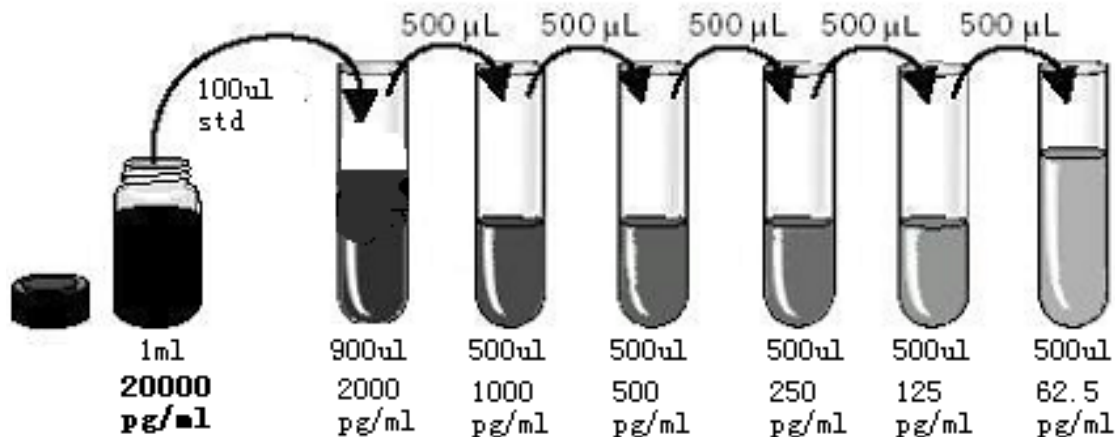


Figure 2: Preparation of IFN- γ standard dilutions

GENERAL ELISA PROTOCOL

1. Prepare all reagents and working standards as directed in the previous sections.
2. Determine the number of microwell strips required to test the desired number of samples plus appropriate number of wells needed for running blanks and standards. Remove extra



microwell strips from holder and store in foil bag with the desiccant provided at 2-8°C sealed tightly.

3. Add 100 μ L of Standard, control, or sample, per well. Cover with the adhesive strip provided. Incubate for 1.5 hours at 37°C.
4. Aspirate each well and wash, repeating the process three times for a total of four washes. Wash by filling each well with Wash Buffer (350 μ L) using a squirt bottle, manifold dispenser or auto-washer. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
5. Add 100 μ L of the working solution of Biotin-Conjugate to each well. Cover with a new adhesive strip and incubate 1 hours at 37°C.
6. Repeat the aspiration/wash as in step 3.
7. Add 100 μ L of the working solution of Streptavidin-HRP to each well. Cover with a new adhesive strip and incubate for 30 minutes at 37°C Avoid placing the plate in direct light.
8. Repeat the aspiration/wash as in step 3.
9. Add 100 μ L of Substrate Solution to each well. Incubate for 10-20 minutes at 37°C. Avoid placing the plate in direct light.
10. Add 100 μ L of Stop Solution to each well. Gently tap the plate to ensure thorough mixing.
11. Determine the optical density of each well immediately, using a microplate reader set to 450 nm.(optionally 630nm as the reference wave length; 610-650nm is acceptable)

TECHNICAL HINTS

1. When mixing or reconstituting protein solutions, always avoid foaming.
2. To avoid cross-contamination, change pipette tips between additions of each standard level, between sample additions, and between reagent additions. Also, use separate reservoirs for each reagent.
3. To ensure accurate results, proper adhesion of plate sealers during incubation steps is necessary.
4. Substrate Solution should remain colorless until added to the plate. Stop Solution should be added to the plate in the same order as the Substrate Solution. Keep Substrate Solution protected from light. Substrate Solution should change from colorless to gradations of blue.
5. A standard curve should be generated for each set of samples assayed. According to the content of tested factors in the sample, appropriate diluted or concentrated samples, it is best to do pre-experiment.



CALCULATION OF RESULTS

1. Average the duplicate readings for each standard, control, and sample and subtract the average zero standard optical density.
2. Create a standard curve by reducing the data using computer software capable of generating a four parameter logistic (4-PL) curve-fit. As an alternative, construct a standard curve by plotting the mean absorbance for each standard on the y-axis against the concentration on the x-axis and draw a best fit curve through the points on the graph.
3. The data may be linearized by plotting the log of the IFN- γ concentrations versus the log of the O.D. and the best fit line can be determined by regression analysis. This procedure will produce an adequate but less precise fit of the data. If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.
4. This standard curve is provided for demonstration only. A standard curve should be generated for each set of samples assayed.

Table 2: Typical data using the Rat Interferon Gamma (IFN- γ) ELISA Assay Kit (Measuring wavelength:450nm, Reference wavelength:630nm)

Standard (pg/ml)	OD.	OD.	Average	Corrected
0	0.048	0.043	0.046	—
31.25	0.115	0.116	0.116	0.115
62.5	0.168	0.160	0.164	0.166
125	0.275	0.278	0.277	0.266
250	0.468	0.459	0.464	0.457
500	0.817	0.824	0.821	0.804
1000	1.345	1.334	1.340	1.358
2000	1.912	1.905	1.909	1.906

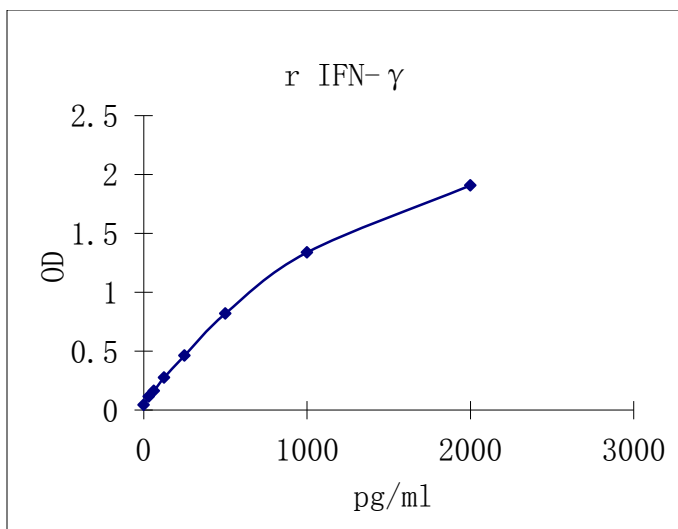


Figure 4: Representative standard curve for Rat Interferon Gamma (IFN- γ) ELISA Assay Kit. IFN- γ was diluted in serial two-fold steps in Sample Diluent.



SENSITIVITY, SPECIFICITY AND REPEATABILITY

1. **REPEATABILITY:** The coefficient of variation of both intra-assay and inter-assay were less than 10%.
2. **SENSITIVITY:** The minimum detectable dose was 7 pg/mL.
3. **SPECIFICITY:** This Rat Interferon Gamma (IFN- γ) ELISA Assay Kit recognizes both natural and recombinant rat IFN- γ . The factors listed below were prepared at 50 ng/ml in Standard /sample Diluent and assayed for cross-reactivity and no significant cross-reactivity or interference was observed.

Table 3: Factors assayed for cross-reactivity

Recombinant human	Recombinant rat
IFN- γ	CINC-1
	GDNF
	β -NGF
	PDGF-BB
	TNF- α

REFERENCES

1. Wheelock, E.F. (1965) Science **149**:310.
2. Ijzermans, J.M. and R.L. Marquet (1989) Immunobiol. **179**:456.
3. Mogensen, S.C. and J.L. Virelizier (1987) Interferon **8**:55.
4. Grossberg, S.E. et al. (1989) Experientia **45**:508.
5. Adolf, G.R. (1985) Oncology (Suppl. 1) **42**:33.
6. Samuel, C.E. (1991) Virology **183**:1.
7. Pellegrini, S. and C. Schindler (1993) Trends Biochem. Sci. **18**:338.
8. Reiter, Z. (1993) J. Interferon Res. **13**:247.
9. Boehm, U. et al. (1997) Annu. Rev. Immunol. **15**:749.
10. Puddu, P. et al. (1997) J. Immunol. **159**:3490.
11. Yoshimoto, T. et al. (1997) Proc. Natl. Acad. Sci. USA **94**:3948.
12. Dijkema, R. et al. (1986) Meth. Enzymol. **119**:453.
13. Rashidbaigi, A. et al. (1986) Proc. Natl. Acad. Sci. USA **83**:384.
14. Pfizenmaier, K. et al. (1988) J. Immunol. **141**:856.
15. Aguet, M. et al. (1988) Cell **55**:273.
16. Fischer, D.G. et al. (1988) J. Biol. Chem. **263**:2632.
17. Calderon, J. et al. (1988) Proc. Natl. Acad. Sci. USA **85**:4837.
18. Paliard, X. et al. (1988) J. Immunol. **141**:849.
19. Christmas, S.E. (1992) Chem. Immunol. **53**:32.
20. Locksley, R.M. and P. Scott (1991) Immunoparasitology Today A58-A61.
21. Billiau, A. and R. Dijkmans (1990) Biochem. Pharmacol. **40**:1433.
22. Sen, G.C. and P. Lengyel (1992) J. Biol. Chem. **267**:5017.
23. Gusella, G.L. et al. (1993) J. Immunol. **151**:2725.



Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.