



**pbl** assay science

---

# **VeriKine™ Human IFN Beta ELISA Kit**

Catalog No. 41410

Assay Range: 50 – 4000 pg/ml

Store IFN- $\beta$  standard at -20°C  
All other components at 2 - 8°C

*Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a  
PBL Assay Science. For research use only.*

*Not for diagnostic or clinical use in, or administration to, humans. Not for resale in  
original or any modified form, including inclusion in a kit, for any purpose.*

*Not for use in the preparation of any commercial product.*

*© Copyright 2013 Pestka Biomedical Laboratories, Inc. All rights reserved.*

## INTRODUCTION

Interferon beta (IFN- $\beta$ ) is synthesized and secreted by fibroblasts and many other cell types in response to pathogens. These pathogens include viruses and bacteria and signaling can occur through Toll-like receptor dependent or independent pathways. Following secretion, IFN- $\beta$  binds to type I interferon receptors on proximal or distal cells activating the JAK1-STAT signaling pathway. Activation of this signal transduction pathway leads to the expression of 2'-5' oligoadenylate synthetases (2'5' OAS), protein kinase R (PKR), MxA proteins, and interferon regulatory factor 7 (IRF-7). The upregulation of IRF-7 expression can exert a positive feedback on IFN- $\beta$  production, whereas the induction of 2'5' OAS activates a latent endonuclease known as RNase L. RNase L cleaves both viral and cellular single stranded mRNA, thereby limiting viral replication and dissemination.

The PBL Assay Science VeriKine™ Human IFN- $\beta$  ELISA kit uses the sandwich immunoassay technique for the quantitative measurement of IFN- $\beta$  in media. It is developed for superior performance with intra-assay and inter-assay CVs of  $\leq 8\%$ .

## **MATERIALS PROVIDED**

- Pre-coated microtiter plate(s)
- Plate sealers
- Wash Solution Concentrate
- Human Interferon Beta Standard, 100,000 pg/ml
- Sample Diluent
- Antibody Concentrate
- HRP Conjugate Concentrate
- Concentrate Diluent
- TMB Substrate Solution
- Stop Solution

## **ADDITIONAL MATERIALS REQUIRED (NOT PROVIDED)**

- Microplate reader capable of reading a wavelength of 450 nm
- Variable volume microtiter pipettes
- Adjustable multichannel pipette (50-200  $\mu$ l)
- Reagent reservoirs
- Wash bottle or plate washing system
- Distilled or deionized water
- Serological pipettes (1, 5, 10 or 25 ml)
- Disposable pipette tips (polypropylene)

**Specifications:** This VeriKine™ kit quantitates human interferon beta in buffers or tissue culture media (TCM) using a sandwich immunoassay.<sup>1,2</sup> The lowest concentration of Hu-IFN-β that can be detected in a test sample is 50 pg/ml. This kit is not designed to measure Hu-IFN-β in human serum. The kit is based on an ELISA with streptavidin conjugated to horseradish peroxidase (HRP). Tetramethyl-benzidine (TMB) is the substrate. The assay is based on the international reference standard for human interferon beta (Hu-IFN-β) provided by the National Institutes of Health.<sup>3</sup>

**Speed:** Incubation time, 3 hr 15 min

**Specificity:** Human IFN-β. No cross reactivity detected with human IFN-α, human IFN-γ, mouse IFN-β or rat IFN-β.

**Storage Conditions/Comments:** For retention of full activity, all reagents except the IFN standard should be kept at 2-8°C in the dark. The standard should be kept at -20°C for one time use only.

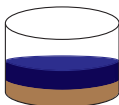
Please note that the dilutions of the Antibody Concentrate and HRP Conjugate Concentrate differ from lot to lot as a result of calibrating each kit for optimal sensitivity. Please refer to the lot specific Certificate of Analysis (COA) for their preparation.

**CAUTION:** The Wash Solution Concentrate, Sample Diluent and Concentrate Diluent contain 0.1% Kathon CG/ICP as a preservative; they should be handled with appropriate safety precautions and discarded properly. For further information, consult the material safety data sheet for Kathon CG/ICP.

For laboratory research use only. Not for use in human diagnostic or therapeutic procedures.

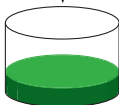
## ASSAY PROCEDURE - QUICK REFERENCE

**Total Time:**  
3 hr, 15 min



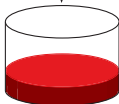
Add **50  $\mu$ l** Sample Diluent  
Add **50  $\mu$ l** Standard, Blank  
or Sample

Incubate **1 hr**  
Aspirate and Wash **3x**



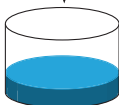
Add **100  $\mu$ l** Diluted  
Ab Solution

Incubate **1 hr**  
Aspirate and Wash **3x**



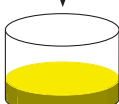
Add **100  $\mu$ l** Diluted  
HRP Solution

Incubate **1 hr**  
Aspirate and Wash **3x**



Add **100  $\mu$ l** TMB  
Substrate

Incubate **15 min** in the dark  
**Do not seal or wash.**



**Note:** ALL  
incubations are at  
room temperature  
(22-25°C)

Add **100  $\mu$ l** Stop Solution  
Read plate within 5 min  
(450 nm)

## PREPARATION OF REAGENTS

Before starting the assay, the plate(s), Wash Solution Concentrate, applicable dilution matrices (e.g. tissue culture media), Sample Diluent, Concentrate Diluent, Stop Solution and samples should be equilibrated to room temperature (RT), 22-25°C. The TMB Substrate Solution should be equilibrated to RT (22-25°C) during step 3 of the Assay Procedure. Supplied Human IFN Beta Standard, Antibody Concentrate and HRP Conjugate Concentrate should be kept on ice (4°C).

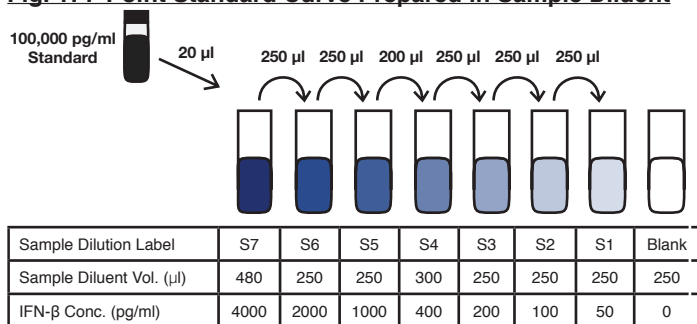
**Wash Buffer:** The Wash Solution Concentrate may contain crystals. Place the bottle in a warm water bath and gently mix until completely dissolved. Prepare a 1:10 working Wash Buffer by adding 50 ml of Wash Solution Concentrate to 450 ml of distilled or deionized water. Mix thoroughly before use. The diluted Wash Buffer can be stored at RT (22-25°C).

**Human Interferon Beta Solution:** Using the Human IFN Beta Standard, provided at 100,000 pg/ml, construct a standard curve (50 - 4000 pg/ml), as shown in Figure 1, in Sample Diluent. In certain situations, test samples may contain substances that can interfere with assay results. Therefore, it is recommended to run the Human IFN Beta standard curve in your test sample matrix.

### **Standard Curve Preparation:**

- a) Label seven polypropylene tubes (S7-S1).
- b) Fill tubes with Sample Diluent as indicated in Figure 1.
- c) Using polypropylene tips, add the indicated amount of Human IFN- $\beta$  Standard to S7 and mix gently. Change tips between each dilution.
- d) Remove indicated amount from S7 and add to S6. Repeat to complete series to S1.
- e) Set aside on ice (4°C) until use in step 1 of the Assay Procedure.

**Fig. 1: 7-Point Standard Curve Prepared in Sample Diluent**



**Sample Preparation:** Prepare test samples of unknown IFN concentration to be tested using Sample Diluent (or applicable matrices) as required. Measurements in duplicate are recommended. Set aside on ice (4°C) until use in step 1 of the Assay Procedure.

**Antibody Solution:** Dilute Antibody Concentrate in Concentrate Diluent. Refer to the lot specific Certificate of Analysis (COA) for the correct volumes to use. Prepare 15 minutes prior to use in step 2 of the Assay Procedure and keep at RT (22-25°C).

**HRP Solution:** Dilute HRP Conjugate Concentrate in Concentrate Diluent. Refer to the lot specific Certificate of Analysis (COA) for the correct volumes to use. Prepare 15 minutes prior to use in step 3 of the Assay Procedure and keep at RT (22-25°C).

## ASSAY PROCEDURE

All incubations should be performed at room temperature (RT), 22-25°C, keeping the plate away from drafts and other temperature fluctuations. Use plate sealers to cover the plate as directed. During all wash steps remove contents of plate by inverting and shaking over a sink and blotting the plate on lint-free absorbent paper; tap the plate. All wells should be filled with a minimum of 250 µl of diluted Wash Buffer at each wash step. Refer to Preparation of Reagents for dilution of concentrated solutions.

**Figure 2: Example of a Typical Plate Setup**

	1	2	3	4	5	6	7	8	9	10	11	12
A	B	B	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa
B	S1	S1	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa
C	S2	S2	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa
D	S3	S3	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa
E	S4	S4	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa
F	S5	S5	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa
G	S6	S6	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa
H	S7	S7	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa	Sa

B = Blanks

S1-S7 = Standard Curve

Sa = Samples

1. **Standards and Test Samples:** Determine the number of microplate strips required to test the desired number of samples plus the appropriate number of wells needed to run blanks and standards. We recommend running the Human IFN Beta Standard, blanks and samples in duplicate or triplicate (see Figure 2 for example plate setup). A standard curve is required for each assay. Remove extra microtiter strips from the frame, seal in the foil bag provided and store at 2-8°C. Unused strips can be used in later assays.



Add 50  $\mu$ l Sample Diluent to the wells. Add 50  $\mu$ l of the diluted Standard Curve, blanks or test samples. Cover with plate sealer and incubate for 1 hour.

After 1 hour, empty the contents of the plate and wash the wells three times with diluted Wash Buffer (refer to Preparation of Reagents).

2. **Antibody Solution:** Add 100  $\mu$ l of diluted Antibody Solution (refer to Preparation of Reagents) to each well. Cover with plate sealer and incubate for 1 hour.

After 1 hour, empty the contents of the plate and wash the wells three times with diluted Wash Buffer.

3. **HRP:** Add 100  $\mu$ l of diluted HRP Solution (refer to Preparation of Reagents) to each well. Cover with plate sealer and incubate for 1 hour. During this incubation period, warm the TMB Substrate Solution to RT (22-25°C).

After 1 hour, empty the contents of the plate and wash the wells three times with diluted Wash Buffer.

4. **TMB Substrate:** Add 100  $\mu$ l of the TMB Substrate Solution to each well. Incubate, in the dark, at RT (22-25°C), for 15 minutes. Do not use a plate sealer during the incubation.

5. **Stop Solution:** After the 15 minute incubation of TMB, DO NOT EMPTY THE WELLS AND DO NOT WASH. Add 100  $\mu$ l of Stop Solution to each well.

6. **Read:** Using a microplate reader, determine the absorbance at 450 nm within 5 minutes after the addition of the Stop Solution.

## CALCULATION OF RESULTS

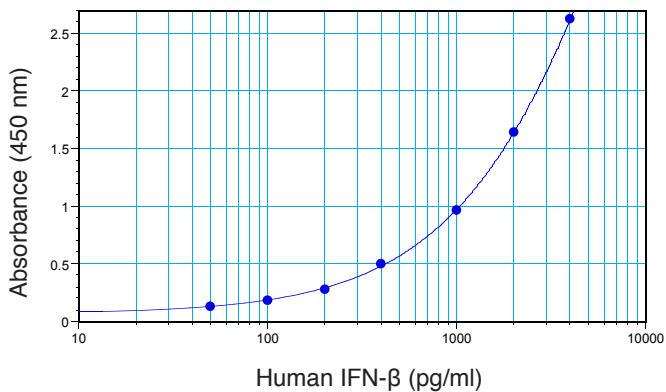
By plotting the optical densities (OD) using a 4-parameter fit for the standard curve, the interferon titer in the samples can be determined. Based on user preference, blank ODs may be subtracted from the standards and sample ODs to eliminate background.

Because the interferon samples are titrated against the international standard, the values from the curves can be determined in units/ml as well as pg/ml. The conversion factor of about 3 – 10 pg/unit of Hu-IFN- $\beta$ , mammalian, is applicable for human interferon beta.<sup>4</sup> Nevertheless, this conversion factor is only an approximation.

A shift in optical densities is typical between users and kit lots. The back fit concentration extrapolated from the standard curve is a more accurate determination of the sample titer and performance of the kit. Variations from the typical curve provided can be a result of operator technique, altered incubation time, fluctuations in temperature, and kit age.

Results of a typical standard curves using a 4-parameter fit are provided in Figure 3 for demonstration only and should not be used to obtain test results. A standard curve must be run for each set of samples assayed.

**Figure 3: Typical Standard Curve**



## REFERENCES

1. Staehelin, T., Stähli, C., Hobbs, D.S., and Pestka, S. (1981) "A Rapid Quantitative Assay of High Sensitivity for Human Leukocyte Interferon with Monoclonal Antibodies," in *Methods in Enzymology*, Vol. 79 (S. Pestka, ed.), Academic Press, New York, 589-595.
2. Kelder, B., Rashidbaigi, A., and Pestka, S. (1986) "A Sandwich Radioimmunoassay for Human IFN $\beta$ ," in *Methods in Enzymology*, Vol. 119 (S. Pestka, ed.), Academic Press, New York, 582-587.
3. Human IFN- $\beta$  international reference standard provided by the NIH, reference no. Gb23-902-531. Pestka, S. (1986) "Interferon Standards and General Abbreviations," in *Methods in Enzymology*, Vol. 119 (S. Pestka, ed.), Academic Press, New York, 14-23.
4. Moschera, J., Woehle, D., Tsai, K., Chen, C. and Tarnowski, J. (1981) "Purification of Recombinant Human Fibroblast Interferon Produced in *Escherichia coli*," in *Methods of Enzymology*, Vol.119 (S. Pestka ed.), Academic Press, New York, 177 – 183.

## PLATE LAYOUT

Use this plate layout as a record of standards and samples assayed.

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H

## NOTES

## NOTES

**PBL Assay Science**

131 Ethel Road West, Suite 6, Piscataway, NJ 08854 USA

Tel: +1 732-777-9123, Fax: +1 732-777-9141

Email: [info@pblassaysci.com](mailto:info@pblassaysci.com), Website: [www.pblassaysci.com](http://www.pblassaysci.com)

---

© 2013 Pestka Biomedical Laboratories, Inc. All rights reserved.

