

VeriKine-HS[™] Human Interferon Beta TCM ELISA Kit **Certificate of Analysis & Protocol**

Assay Range: 2.34-150 pg/ml Compatibility: Tissue Culture Media Assay Length: 3 hr

Catalog No: 41435-1

Lot No:

Expiration:

Store all components at 2-8°C

Kit Components	Part No.	Lot No.	Quantity	
Plate(s)	SMP138		1	
Plate Sealers	N/A	N/A	4	
Wash Solution Concentrate	SMP057-60		2 x 50 ml	
Human IFN-Beta Standard, 100,000 pg/ml	SMP146-1		1 vial	
Standard Diluent	SMP163-30		25 ml	
Sample Buffer	SMP147-15		15 ml	
Antibody Concentrate	SMP148-1		1 vial	
HRP Conjugate Concentrate	SMP056-320		1 vial	
Assay Diluent	ASD-30		25 ml	
TMB Substrate Solution	KET-15		15 ml	
Stop Solution	SCY-15		15 ml	

Authorization

Released by:

Date:

Note: The 41435-1 ELISA is not intended for use with human serum or plasma samples because of a significantly elevated risk of false positive results.

Visit the product page on PBL's website (https://pblassaysci.com) to view the technical supplement and additional product information.

CAUTION: Components should be handled with appropriate safety precautions and discarded properly. For further information, consult the safety data sheet (SDS).

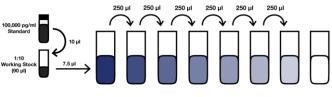
PREPARATION OF REAGENTS

Wash Buffer: Wash Solution Concentrate may contain crystals; place in a warm water bath and gently mix until completely dissolved. Prepare a 1:10 working wash solution (e.g. Add 50 ml Wash Solution Concentrate to 450 ml distilled or deionized water). Mix thoroughly before use. (Note: Prepare fresh Wash Buffer for each assay run.)

Human IFN-Beta Standard Curve Preparation:

- a. Label seven polypropylene tubes (S1 S7).
- b. Add volume of Standard Diluent or Sample Matrix to each tube as indicated in Figure 1.
- c. Using polypropylene tips, add 10 µl of the Human IFN-Beta Standard to 90 µl of Standard Diluent or Sample Matrix.
- d. Add 7.5 µl of prediluted standard to S7 and mix thoroughly.
- e. Transfer 250 µl of S7 to S6 and mix thoroughly. Repeat to complete series to S1. Set aside until use in step 1.

Figure 1: 7-Point Standard Curve Prepared in Standard Diluent



Label	S7	S6	S5	S4	S3	S2	S1	Blank
Standard Diluent (µl)	492.5	250	250	250	250	250	250	250
IFN-Beta (pg/ml)	150.00	75.00	37.50	18.75	9.38	4.69	2.34	0

Sample Preparation: Thaw frozen sample tubes to Room Temperature (RT) (22-25°C) in either tap water or between the fingertips. If samples require dilution, prepare using Standard Diluent or Sample Matrix. Keep at RT until use. Measurements in duplicate are recommended.

Antibody Solution: Prior to use in step 1, dilute Antibody Concentrate in the volume of Assay Diluent as shown below. Keep at RT (22-25°C).

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µI)						
Assay Diluent (ml)	2.0	3.0	4.0	5.0	6.0	7.0

HRP Solution: 15 minutes prior to use in step 2, dilute HRP Conjugate Concentrate in the volume of Assay Diluent as shown below. Keep at RT (22-25°C).

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µI)						
Assay Diluent (ml)	3.0	5.0	6.0	8.0	10.0	12.0

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. © 2020 Pestka Biomedical Laboratories, Inc. All rights reserved. Rev. 00

ASSAY PROCEDURE

Bring to RT (22-25°C)	Keep at 2-8°C		
Plate/Sealers	Human IFN-Beta Standard		
Sample Buffer	Antibody Concentrate		
Standard Diluent	HRP Conjugate Concentrate		
Matrices/Samples			
Assay Diluent			
TMB Substrate Solution			
Stop Solution			

- Incubations: Use plate sealers to cover the plate when directed. All incubations should be conducted in a closed chamber at 22-25°C or at RT, keeping the plate away from drafts.
- Plate Washing: All wells should be filled with a minimum of 300 µl of Wash Solution. Remove plate contents by inverting and blotting the plate on lint-free absorbent paper; tap the plate dry.

1. Determine the number of microplate strips required. We recommend running both the standard and samples at least in duplicate. 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.

Total well volume = 150 µl (Step A + Step B)

Step A: Add 50 μI of Sample Buffer and 50 μI of diluted Antibody Solution to every well.

Step B: Add 50 μ I of Standard, Test Samples or Blanks (Standard Diluent or appropriate dilution matrix) to each designated well.

Cover with Plate Sealer and shake at 450 rpm at RT for 2 hours.

After 2 hours, empty plate contents and wash wells three times.

2. Add 100 µl of diluted HRP Solution to each well.

Cover with Plate Sealer and shake at 450 rpm at RT for 30 minutes.

After 30 minutes, empty plate contents and wash wells four times.

3. Add 100 μl of TMB Substrate Solution to each well. Incubate in the dark at RT for 30 minutes. Do not use a Plate Sealer during the incubation. DO NOT SHAKE.

4. After 30 minutes, DO NOT EMPTY THE WELLS AND DO NOT WASH. Add **100 \muI** of **Stop Solution** to each well.

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

HUMAN IFN-BETA ELISA (41435) ASSAY PROCEDURE – QUICK REFERENCE

Total Time: 3 hr

Note: All incubations are at Room Temperature (RT) (22-25°C)*



 Add 50 µI Sample Buffer
Add 50 µI Diluted Antibody Solution
Add 50 µI Standard, Sample or Blank Incubate 2 hr (shake at 450 rpm) at RT*

Aspirate and Wash 3x



Add **100 µI** diluted HRP Solution Incubate **30 min** (shake at 450 rpm) at RT*

Aspirate and Wash **4x**



Add **100 µI** TMB Substrate Incubate **30 min<u>in the dark</u> Do not seal, shake or wash.**

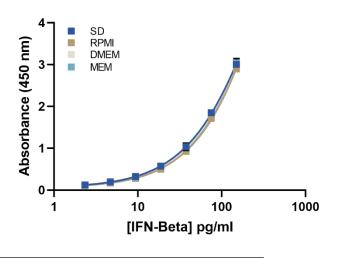


Add **100 µl** Stop Solution Read plate within 5 min (450 nm)

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. Blank ODs may be subtracted from the standards and sample ODs to eliminate background. Use conversion factor of 3 pg/unit to approximate titers in units/ml. A 4-parameter logistic plot with 1/y² weighted analysis is recommended for obtaining optimal fit of standard curve OD values.

Figure 2: Typical Standard Curve



Visit the product page on PBL's website (<u>https://pblassaysci.com</u>) to view technical supplement and additional product information