

# VeriKine-HS<sup>™</sup> Human IL-22 ELISA Kit

**Certificate of Analysis & Protocol** 

Assay Range: 0.78 - 50 pg/ml Compatibility: Serum, Plasma, Tissue Culture Media (TCM) Assay Length: 4 hr

#### Catalog No: 41701-1 Lot No:

## Expiration:

Store all components at 2-8°C

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP326		1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60		2 x 50 ml
Human IL-22 Standard, 10,000 pg/ml	SMP328-1		1 vial
Assay Buffer	SMP329-8		8 ml
Standard Diluent	SMP330-60		55 ml
Antibody Concentrate	SMP331-1		1 vial
HRP Conjugate Concentrate	SMP056-240		1 vial
Concentrate Diluent	SMP024-15		15 ml
HRP Diluent	ASDHRP-15		15 ml
TMB Substrate Solution	KET-15		15 ml
Stop Solution	SCY-15		15 ml

## Authorization

Released by:

Date:

Visit the product page on PBL's website (https://pblassaysci.com) to view the technical supplement, including performance characterization and kit specifications.

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

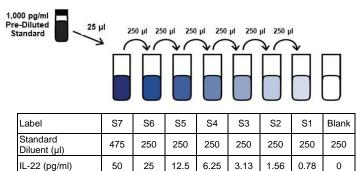
## **PREPARATION OF REAGENTS**

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 IL-22 Standard Curve Preparation:

- a. Prepare a 1:10 working stock of Human IL-22 standard by pipetting 20 µl of IL-22 standard into 180 µl of Standard Diluent. Only use the provided Standard Diluent for this step. Mix thoroughly by gently pipetting up and down 5 times.
- b. Label seven polypropylene tubes (S1 S7).
- c. Add indicated volume of Standard Diluent or Sample Matrix to each tube as indicated in Figure 1.
- d. Using polypropylene tips, add 25 µl of pre-diluted IL-22 Standard working stock to S7 and mix gently. Remove indicated amount from S7 and add to S6. Repeat to complete series to S1.

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



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. Keep on ice (2-8°C) until step 1. Measurements in duplicate are recommended.

Antibody Concentrate: 15 minutes prior to use in step 3, dilute Antibody Concentrate in the volume of Concentrate Diluent as shown below. Keep at RT (22-25°C).

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate						
Concentrate Diluent (ml)	2.0	4.0	6.0	8.0	10.0	12.0

HRP Solution: 15 minutes prior to use in step 4, dilute HRP Conjugate Concentrate in the volume of HRP Diluent as shown below. Keep at RT (22-25°C).

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate						
HRP Diluent (ml)	2.0	4.0	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 2-8ºC
Plate/Plate Sealers	All other components
Standard Diluent	
Assay Buffer	
Concentrate Diluent	
HRP Diluent	
TMB Substrate Solution	
Stop Solution	
Matrices/Samples	

- Incubations: Use plate sealers to cover the plate when directed. All incubations should be conducted in a closed chamber at 22-25°C (RT), keeping the plate away from drafts.
- **Plate Washing**: All wells should be filled with a minimum of 300 µl of Wash Buffer. 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.

#### 2. Total well volume = 100 µl (Step A + Step B)

Step A: Add 50 µl of Assay Buffer to every well.

Step B: Add 50  $\mu l$  of Standard, Test Samples or Blanks (Standard Diluent or appropriate dilution matrix) to each designated well.

Cover with Plate Sealer and shake at 600 rpm at RT for 2 hours.

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

3. Add 100  $\mu$ I of diluted Antibody Solution to each well. Cover with Plate Sealer and shake plate at 600 rpm at RT for 1 hour.

After 1 hour, empty plate contents and wash wells three times.

**4.** Add **100 \muI** of diluted **HRP Solution** to each well. Cover with Plate Sealer and shake plate at 600 rpm at RT for 30 minutes.

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

5. Add 100  $\mu$ I of TMB Substrate Solution to each well. Incubate in the dark at RT for 30 minutes. Do not use a Plate Sealer and DO NOT SHAKE during the incubation.

6. After 30 minutes, DO NOT EMPTY THE WELLS AND DO NOT WASH. Add 100  $\mu l$  of Stop Solution to each well.

**7.** Using a microplate reader, determine the absorbance at 450 nm within 2 minutes after the addition of Stop Solution.

## HUMAN IL-22 ELISA (41701) **ASSAY PROCEDURE – QUICK REFERENCE** Total Time: 4 hr Note: All incubations are at Room Temperature (RT) (22-25°C)\* 1. Add 50 µl Assay Buffer 2. Add 50 µl Standard, Sample or Blank Incubate 2 hr (shake at 600 rpm) at RT\* Aspirate and Wash 3x Add 100 µl diluted Antibody Solution Incubate 1 hr (shake at 600 rpm) at RT\* Aspirate and Wash 3x Add 100 µl diluted HRP Solution Incubate 1 hr (shake at 600 rpm) at RT\* Aspirate and Wash 4x Add 100 µI TMB Substrate Incubate 30 min in the dark at RT\* Do not seal, shake or wash.

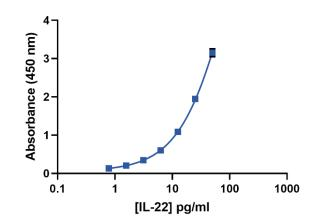


Add **100 µl** Stop Solution Read plate within 2 min (450 nm)

## **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. A 4-parameter logistic plot with  $1/y^2$  weighted analysis is recommended for obtaining optimal fit of standard curve OD values. Blank ODs may be subtracted from the standards and sample ODs to eliminate background.

#### Figure 2: Typical Standard Curve in Standard Diluent



Visit the product page on PBL's website (<u>https://pblassaysci.com</u>) to view the technical supplement, including performance characterization and kit specifications.