



VeriKine™ Cynomolgus Interferon Beta ELISA Kit

Certificate of Analysis & Protocol

Assay Range: 5.47 - 350 pg/ml
 Compatibility: Serum, Plasma, Tissue Culture Media (TCM)
 Assay Length: 3 hr

Catalog No: 46415-1

Lot No: 7236

Expiration: June 30, 2021

Store all components at 2-8°C

| Kit Components | Part No. | Lot No. | Quantity |
|---------------------------------------|------------|-----------|-----------|
| Plate(s) | SMP138 | K6098 | 1 |
| Plate Sealers | N/A | N/A | 4 |
| Wash Solution Concentrate | SMP057-60 | K6101 | 2 x 50 ml |
| Cyno IFN Beta Standard, 100,000 pg/ml | SMP261-1 | K6102 | 1 vial |
| Standard Diluent | SMP163-30 | K6106 | 25 ml |
| Sample Buffer | SMP147-15 | K6103 | 15 ml |
| Antibody Concentrate | SMP148-100 | K6104 | 1 vial |
| HRP Conjugate Concentrate | SMP056-450 | K6105 | 1 vial |
| Assay Diluent | ASD-30 | 625611 | 25 ml |
| TMB Substrate Solution | KET-15 | 200210D02 | 15 ml |
| Stop Solution | SCY-15 | 56334 | 15 ml |

Authorization

Released by: _____

Date: June 12, 2020

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

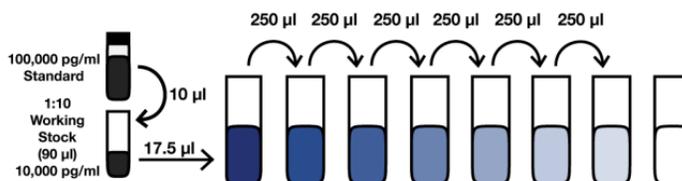
PREPARATION OF REAGENTS

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.

Cyno IFN-Beta Standard Curve Preparation:

- Prepare a 1:10 *working stock* of Cyno IFN-Beta standard by pipetting 10 µl of IFN standard into 90 µl of Standard Diluent or Sample Matrix. Mix thoroughly by gently pipetting up and down 10 times.
- Label seven polypropylene tubes (S1 – S7).
- Add indicated volume of Standard Diluent or Sample Matrix to each tube as indicated in [Figure 1](#).
- Using polypropylene tips, add 17.5 µl of working stock to S7 as indicated and mix thoroughly by pipetting up and down 10 times. Remove indicated amount from S7 and add to S6. Repeat to complete series to S1. *Change tips between each dilution.*

Figure 1: 7-Point Standard Curve Prepared in Sample Matrix



| Label | S7 | S6 | S5 | S4 | S3 | S2 | S1 | Blank |
|----------------------|-------|-----|------|-------|-------|-------|------|-------|
| Dilution Matrix (µl) | 482.5 | 250 | 250 | 250 | 250 | 250 | 250 | 250 |
| IFN-Beta (pg/ml) | 350 | 175 | 87.5 | 43.75 | 21.88 | 10.94 | 5.47 | 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 (µl) | 16 | 24 | 32 | 40 | 48 | 56 |
| 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 (µl) | 24 | 40 | 48 | 64 | 80 | 96 |
| Assay Diluent (ml) | 3.0 | 5.0 | 6.0 | 8.0 | 10.0 | 12.0 |

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

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 | All other components |
| Wash Solution Concentrate | |
| Standard Diluent | |
| Sample Buffer | |
| Assay Diluent | |
| TMB Substrate Solution | |
| Stop Solution | |
| Matrices/Samples | |

- **Incubations:** All incubations should be conducted in a closed chamber at 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.

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

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

Step B: Add 50 µl of diluted **Antibody Solution** to each well.

Step C: Add 50 µl of diluted **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 plate at 450 rpm at RT for 30 minutes.

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

3. Add 100 µl of diluted **TMB Substrate Solution** to each well. Do not use a Plate Sealer and DO NOT SHAKE during the incubation.

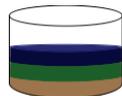
4. After 30 minutes, DO NOT EMPTY THE WELLS AND DO NOT WASH. Add 100 µl of **Stop Solution** to each well.

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

CYNO IFN-BETA ELISA (46415) ASSAY PROCEDURE – QUICK REFERENCE

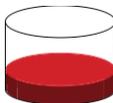
Total Time: 3 hr

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



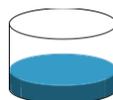
1. Add 50 µl Sample Buffer
 2. Add 50 µl Diluted Antibody Solution
 3. Add 50 µl Standard, Sample or Blank
- Incubate 2 hr (shake at 450 rpm) at RT*

Aspirate and Wash 3x

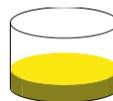


- Add 100 µl diluted HRP Solution
- Incubate 30 min (shake at 450 rpm) at RT*

Aspirate and Wash 4x



- Add 100 µl TMB Substrate
- Incubate 30 min in the dark
- Do not seal, shake or wash.

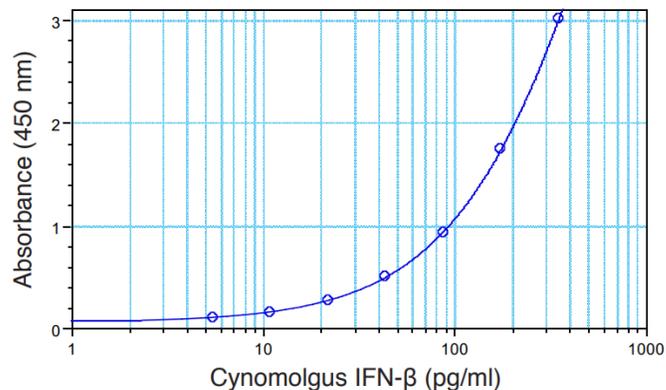


- Add 100 µl Stop Solution
- Read plate within 2 min (450 nm)

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. 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 (<https://pblsaysci.com>) to view the full protocol, including performance characterization and kit specifications.