



### Certificate of Analysis

#### VeriKine™ Human Interferon Beta ELISA Kit

Catalog No: 41410-1  
Lot No: 7193

Expiration: December 31, 2021

| Kit Components – Store at 2-8°C        | Part No.   | Lot No.   | Quantity |
|--|------------|-----------|----------|
| Plate(s)                               | SMP055     | K5954     | 1        |
| Plate Sealers                          | N/A        | N/A       | 4        |
| Wash Solution Concentrate              | SMP057-60  | K5493     | 50 ml    |
| Human IFN Beta Standard, 100,000 pg/ml | SMP146-1   | K5957     | 1 vial   |
| Sample Diluent                         | SMP067-30  | K5958     | 30 ml    |
| Antibody Concentrate                   | SMP062-1   | K5960     | 1 vial   |
| HRP Conjugate Concentrate              | SMP056-180 | K5961     | 1 vial   |
| Concentrate Diluent                    | SMP024-30  | K5962     | 30 ml    |
| TMB Substrate Solution                 | KET-15     | 180803D02 | 15 ml    |
| Stop Solution                          | SCY-15     | 51066     | 15 ml    |

#### Lot Specific Preparations

(Refer to VeriKine™ Human Interferon Beta ELISA Kit Protocol for full instructions.)

#### Antibody Solution Preparation

| Micro-plate Strips Used          | 2   | 4   | 6   | 8   | 10   | 12   |
|----------------------------------|-----|-----|-----|-----|------|------|
| <b>Antibody Concentrate (µl)</b> | 5   | 10  | 15  | 20  | 25   | 30   |
| <b>Concentrate Diluent (ml)</b>  | 2.0 | 4.0 | 6.0 | 8.0 | 10.0 | 12.0 |

#### HRP Solution Preparation

| Micro-plate Strips Used               | 2   | 4   | 6   | 8   | 10   | 12   |
|---------------------------------------|-----|-----|-----|-----|------|------|
| <b>HRP Conjugate Concentrate (µl)</b> | 7   | 13  | 20  | 27  | 33   | 40   |
| <b>Concentrate Diluent (ml)</b>       | 2.0 | 4.0 | 6.0 | 8.0 | 10.0 | 12.0 |

#### Performance Characteristics

Intra-Assay CV: ≤ 8%  
Inter-Assay CV: ≤ 8%

#### Authorization

Released by: 

Date: March 13, 2020

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.