



Certificate of Analysis

VeriKine-HS™ Mouse Interferon Beta Serum ELISA

Catalog No: 42410-1

Lot No:

Expiration:

| Kit Components | Part No. | Lot No. | Quantity |
|---|------------|---------|-----------|
| Plate(s) | SMP199 | | 1 |
| Plate Sealers | N/A | N/A | 4 |
| Wash Solution Concentrate | SMP057-60 | | 2 x 50 ml |
| Mouse IFN Beta Serum Standard, 10,000 pg/ml | SMP195-1 | | 1 vial |
| Sample Diluent | SMP196-30 | | 25 ml |
| Antibody Concentrate | SMP197-1 | | 1 vial |
| HRP Conjugate Concentrate | SMP056-150 | | 1 vial |
| Antibody Diluent | SMP198-15 | | 15 ml |
| HRP Diluent | ASDHRP-15 | | 15 ml |
| Serum Buffer | SMP213-15 | | 12 ml |
| TMB Substrate Solution | KET-15 | | 15 ml |
| Stop Solution | SCY-15 | | 15 ml |

Lot Specific Preparations

(Refer to VeriKine-HS™ Mouse Interferon Beta Serum ELISA Protocol for full instructions.)

Antibody Solution Preparation

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

HRP Solution Preparation

| Micro-plate Strips Used | 2 | 4 | 6 | 8 | 10 | 12 |
|--------------------------------|-----|-----|-----|-----|-----|-----|
| HRP Conjugate Concentrate (µl) | | | | | | |
| HRP Diluent (ml) | 2.0 | 3.0 | 4.0 | 5.0 | 6.0 | 7.0 |

Product Performance Characteristics

Intra-Assay CV: m6%

Inter-Assay CV: m7%

Average Spike Recovery: 95%

Authorization

Released by: _____

Date:

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.