



Certificate of Analysis

VeriKine™ Human Interferon Alpha Multi-Subtype Serum ELISA Kit

Catalog No: 41110-1

Lot No: 7232

Expiration: February 28, 2022

| Kit Components | Part No. | Lot No. | Quantity |
|---------------------------------------|------------|-----------|----------|
| Plate(s) | SMP047 | K6076 | 1 |
| Plate Sealers | N/A | N/A | 4 |
| Wash Solution Concentrate | SMP022-60 | K5832 | 50 ml |
| Human IFN Alpha Standard 10,000 pg/ml | SMP049-1 | K6079 | 1 vial |
| Sample Diluent | SMP233-30 | K6080 | 30 ml |
| Antibody Concentrate | SMP048-1 | K6081 | 1 vial |
| HRP Conjugate Concentrate | SMP050-240 | K6082 | 1 vial |
| Concentrate Diluent | SMP024-15 | K6001 | 15 ml |
| TMB Substrate Solution | KET-15 | 200210D02 | 15 ml |
| Stop Solution | SCY-15 | 56334 | 15 ml |
| Dilution Buffer | SMP231-30 | K6083 | 30 ml |

Lot Specific Preparations

(Refer to VeriKine™ Human IFN Alpha Multi-Subtype Serum ELISA Kit Protocol for full instructions.)

Antibody Solution Preparation

| Micro-plate Strips Used | 2 | 4 | 6 | 8 | 10 | 12 |
|---------------------------|----|----|----|----|-----|-----|
| Antibody Concentrate (µl) | 20 | 40 | 60 | 80 | 100 | 120 |
| Dilution Buffer (ml) | 2 | 4 | 6 | 8 | 10 | 12 |

HRP Solution Preparation

| Micro-plate Strips Used | 2 | 4 | 6 | 8 | 10 | 12 |
|--------------------------------|----|----|----|----|----|----|
| HRP Conjugate Concentrate (µl) | 10 | 20 | 30 | 40 | 50 | 60 |
| Concentrate Diluent (ml) | 2 | 4 | 6 | 8 | 10 | 12 |

Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Spike Recovery: ≥ 90%

Authorization

Released by: _____

Date: June 2, 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.