



Certificate of Analysis

VeriKine™ Human Interferon Alpha Multi-Subtype ELISA Kit

Catalog No: 41105-1

Lot No: 7433

Expiration: May 31, 2023

| Kit Components | Part No. | Lot No. | Quantity |
|--------------------------------------|------------|-----------|----------|
| Plate(s) | SMP047 | K6647 | 1 |
| Plate Sealers | N/A | N/A | 4 |
| Wash Solution Concentrate | SMP022-60 | K6446 | 50 ml |
| Human IFN Alpha Standard 10,000pg/ml | SMP049-1 | K6650 | 1 vial |
| Dilution Buffer | SMP231-60 | K6654 | 50 ml |
| Antibody Concentrate | SMP048-1 | K6652 | 1 vial |
| HRP Conjugate Concentrate | SMP050-360 | K6653 | 1 vial |
| Concentrate Diluent | SMP024-15 | K6371 | 15 ml |
| TMB Substrate Solution | KET-15 | 210201D01 | 15 ml |
| Stop Solution | SCY-15 | 59172 | 15 ml |

Lot Specific Preparations

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

Antibody Solution Preparation

| Micro-plate Strips Used | | 2 | 4 | 6 | 8 | 10 | 12 |
|----------------------------------|--------------------------------------|-----|-----|-----|-----|------|------|
| Antibody Concentrate (µl) | High Sensitivity 12.5 - 500 pg/ml | 20 | 40 | 60 | 80 | 100 | 120 |
| | Extended Range 156 - 5000 pg/ml | 10 | 20 | 30 | 40 | 50 | 60 |
| Dilution Buffer (ml) | | 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 |
|---------------------------------------|--------------------------------------|-----|-----|-----|-----|------|------|
| HRP Conjugate Concentrate (µl) | High Sensitivity 12.5 - 500 pg/ml | 27 | 53 | 80 | 107 | 133 | 160 |
| | Extended Range 156 - 5000 pg/ml | 10 | 20 | 30 | 40 | 50 | 60 |
| Concentrate Diluent (ml) | | 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: September 7, 2021

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.