



### Certificate of Analysis

#### VeriKine™ Human Interferon Alpha Multi-Subtype Serum ELISA Kit

**Catalog No:** 41110-1

**Lot No:** 6377

**Expiration:** March 31, 2017

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP047	K3674	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K3665	50 ml
Human IFN Alpha Standard 10,000 pg/ml	SMP049-1	K3865	1 vial
Sample Diluent	SMP233-30	K3869	30 ml
Antibody Concentrate	SMP048-1	K3867	1 vial
HRP Conjugate Concentrate	SMP050-240	K3868	1 vial
Concentrate Diluent	SMP024-15	K3383	15 ml
TMB Substrate Solution	KET-15	131026	15 ml
Stop Solution	SCY-15	32954	15 ml
Dilution Buffer	SMP231-30	K3866	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
<b>Antibody Concentrate (µl)</b>	20	40	60	80	100	120
<b>Dilution Buffer (ml)</b>	2	4	6	8	10	12

#### HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
<b>HRP Conjugate Concentrate (µl)</b>	13	27	40	53	67	80
<b>Concentrate Diluent (ml)</b>	2	4	6	8	10	12

#### Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Spike Recovery: ≥ 90%

#### Authorization

Released by: \_\_\_\_\_

Date: December 14, 2015

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.