



### Certificate of Analysis

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

Catalog No: 41110-2

Lot No:

Expiration:

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP047		5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP022-250		250 ml
Human IFN Alpha Standard 10,000 pg/ml	SMP049-2		2.5 ml
Sample Diluent	SMP233-150		150 ml
Antibody Concentrate	SMP048-2		1 vial
HRP Conjugate Concentrate	SMP050-1000		1 vial
Concentrate Diluent	SMP024-60		60 ml
TMB Substrate Solution	KET-60		60 ml
Stop Solution	SCY-60		60 ml
Dilution Buffer	SMP231-150		150 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)						
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)						
Concentrate Diluent (ml)	2	4	6	8	10	12

#### Performance Characteristics

Intra-Assay CV: m8%

Inter-Assay CV: m8%

Spike Recovery: ~ 90%

#### 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.