



### Certificate of Analysis

#### VeriKine-HS™ Human Interferon Alpha All Subtype TCM ELISA Kit

Catalog No: 41135-1

Lot No:  

Expiration:  

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP217		1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60		2 x 50 ml
Human IFN Alpha Standard 10,000 pg/ml	SMP049-400		1 vial
Assay Buffer	SMP324-8		8 ml
Standard Diluent	SMP323-60		55 ml
Antibody Concentrate	SMP322-1		1 vial
Antibody Diluent	SMP315-15		15 ml
HRP Conjugate Concentrate	SMP056-240		1 vial
HRP Diluent	ASDHRP-15		15 ml
TMB Substrate Solution	KET-15		15 ml
Stop Solution	SCY-15		15 ml

#### Lot Specific Preparation

(Refer to VeriKine-HS™ Human IFN  $\alpha$  All Subtype TCM ELISA Kit Protocol for full instructions.)

#### Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
<b>Antibody Concentrate (<math>\mu</math>l)</b>						
<b>Antibody Diluent (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 (<math>\mu</math>l)</b>						
<b>HRP Diluent (ml)</b>	2	4	6	8	10	12

#### Performance Characteristics

Intra-Assay CV: m8%

Inter-Assay CV: m10%

Intermediate Precision\* CV: m14%

\*(25 assays performed by 10 operators using 3 lots of components)

Average Spike Recovery: 105%

#### 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.