



Certificate of Analysis

VeriKine-HS™ Human Interferon Alpha All Subtype ELISA Kit

Catalog No: 41115-1

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
Sample Buffer	SMP220-8		8 ml
Standard Diluent	SMP218-60		55 ml
Antibody Concentrate	SMP219-1		1 vial
Antibody Diluent	SMP304-15		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 α All Subtype ELISA Kit Protocol for full instructions.)

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (μl)						
Antibody Diluent (ml)	2	4	6	8	10	12

HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (μl)						
HRP Diluent (ml)	2	4	6	8	10	12

Performance Characteristics

Intra-Assay CV: \leq 8%

Inter-Assay CV: \leq 10%

Average Spike Recovery: 112%

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.