



Certificate of Analysis

VeriKine-HS™ Human Interferon Alpha All Subtype ELISA Kit

Catalog No: 41115-1

Lot No: 7146

Expiration: December 31, 2020

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP217	K5822	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K5454	2 x 50 ml
Human IFN Alpha Standard 10,000 pg/ml	SMP049-400	K5825	1 vial
Sample Buffer	SMP220-8	K5830	8 ml
Standard Diluent	SMP218-60	K5826	55 ml
Antibody Concentrate	SMP219-1	K5827	1 vial
Antibody Diluent	SMP304-15	K5828	15 ml
HRP Conjugate Concentrate	SMP056-240	K5829	1 vial
HRP Diluent	ASDHRP-15	935569	15 ml
TMB Substrate Solution	KET-15	180803D02	15 ml
Stop Solution	SCY-15	51066	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)	20	40	60	80	100	120
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)	10	20	30	40	50	60
HRP Diluent (ml)	2	4	6	8	10	12

Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 10%

Average Spike Recovery: 112%

Please note for this lot:

A 4-parameter logistic plot with 1/y weighted analysis may be recommended for obtaining optimal fit of standard curve OD values.

Authorization

Released by: _____

Date: December 17, 2019

Sold under license from Pestka Biomedical Laboratory, 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.

Rev. 03