



Certificate of Analysis

VeriKine-HS™ Human Interferon Alpha All Subtype ELISA Kit

Catalog No: 41115-1

Lot No: 7334

Expiration: January 31, 2022

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP217	K6414	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K6226	2 x 50 ml
Human IFN Alpha Standard 10,000 pg/ml	SMP049-400	K6417	1 vial
Sample Buffer	SMP220-8	K6422	8 ml
Standard Diluent	SMP218-60	K6418	55 ml
Antibody Concentrate	SMP219-1	K6419	1 vial
Antibody Diluent	SMP304-15	K6420	15 ml
HRP Conjugate Concentrate	SMP056-240	K6421	1 vial
HRP Diluent	ASDHRP-15	579381	15 ml
TMB Substrate Solution	KET-15	200210D02	15 ml
Stop Solution	SCY-15	59172	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 (μ l)	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 (μ l)	10	20	30	40	50	60
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: February 12, 2021

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.