



Certificate of Analysis

VeriKine™ Human Interferon Alpha Multi-Subtype ELISA Kit

Catalog No: 41105-2

Lot No: 6787

Expiration:

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP047	K4736	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP022-250	K4593	250 ml
Human IFN Alpha Standard 10,000pg/ml	SMP049-2	K4739	1 vial
Dilution Buffer	SMP231-250	K4743	250 ml
Antibody Concentrate	SMP048-2	K4741	1 vial
HRP Conjugate Concentrate	SMP050-1800	K4742	1 vial
Concentrate Diluent	SMP024-60	K4525	60 ml
TMB Substrate Solution	KET-60	150503	60 ml
Stop Solution	SCY-60	44142	60 ml

Lot Specific Preparations

(Refer to VeriKine™ Human Interferon Alpha Multi-Subtype ELISA Kit Protocol for full instructions.)

Antibody Solution Preparation

Micro-plate Strips Used		2	4	6	8	10	12
Antibody Concentrate (µl)	High Sensitivity 12.5-500 pg/ml	20	40	60	80	100	120
	Extended Range 156-5000 pg/ml	10	20	30	40	50	60
Dilution Buffer (ml)		2.0	4.0	6.0	8.0	10.0	12.0

HRP Solution Preparation

Micro-plate Strips Used		2	4	6	8	10	12
HRP Conjugate Concentrate (µl)	High Sensitivity 12.5-500 pg/ml	40	80	120	160	200	240
	Extended Range 156-5000 pg/ml	13	27	40	53	67	80
Concentrate Diluent (ml)		2.0	4.0	6.0	8.0	10.0	12.0

Performance Characteristics

Intra-Assay CV: m8%

Inter-Assay CV: m8%

Authorization

Released by: _____

Date: September 13, 2017

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.