



### Certificate of Analysis

#### VeriKine™ Human Interferon Alpha Multi-Subtype ELISA Kit

Catalog No: 41105-2

Lot No: 7446

Expiration: June 30, 2023

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP047	K6711	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP022-250	K6596	250 ml
Human IFN Alpha Standard 10,000pg/ml	SMP049-2	K6715	1 vial
Dilution Buffer	SMP231-250	K6718	250 ml
Antibody Concentrate	SMP048-2	K6716	1 vial
HRP Conjugate Concentrate	SMP050-1800	K6717	1 vial
Concentrate Diluent	SMP024-60	K6504	60 ml
TMB Substrate Solution	KET-60	210201D02	60 ml
Stop Solution	SCY-60	62731	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
<b>Antibody Concentrate (µl)</b>	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
<b>Dilution Buffer (ml)</b>		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
<b>HRP Conjugate Concentrate (µl)</b>	High Sensitivity 12.5-500 pg/ml	27	53	80	107	133	160
	Extended Range 156-5000 pg/ml	10	20	30	40	50	60
<b>Concentrate Diluent (ml)</b>		2.0	4.0	6.0	8.0	10.0	12.0

#### Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

#### Authorization

Released by: 

Date: October 5, 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.