



### Certificate of Analysis

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

Catalog No: 41110-1

Lot No: 7523

Expiration: February 28, 2024

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP047	K6916	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K6596	50 ml
Human IFN Alpha Standard 10,000 pg/ml	SMP049-1	K6919	1 vial
Sample Diluent	SMP233-30	K6920	30 ml
Antibody Concentrate	SMP048-1	K6921	1 vial
HRP Conjugate Concentrate	SMP050-240	K6922	1 vial
Concentrate Diluent	SMP024-15	K6731	15 ml
TMB Substrate Solution	KET-15	210201D01	15 ml
Stop Solution	SCY-15	62731	15 ml
Dilution Buffer	SMP231-30	K6923	30 ml

#### Lot Specific Preparations

(Refer to VeriKine™ Human IFN Alpha Multi-Subtype Serum 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
Dilution Buffer (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
Concentrate Diluent (ml)	2	4	6	8	10	12

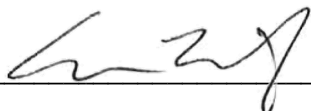
#### Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Spike Recovery: ≥ 90%

#### Authorization

Released by: 

Date: June 2, 2022

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.