



### Certificate of Analysis

#### VeriKine™ Human Interferon Alpha ELISA Kit

Catalog No: 41100-1

Lot No: 7762

Expiration: September 30, 2027

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP064	K7687	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K7693	50 ml
Human IFN Alpha Standard, 10,000pg/ml	SMP049-1	K7690	1 vial
Dilution Buffer	SMP021-60	K7694	50 ml
Antibody Concentrate	SMP066-1	K7691	1 vial
HRP Conjugate Concentrate	SMP050-240	K7692	1 vial
Concentrate Diluent	SMP024-15	K7592	15 ml
TMB Substrate Solution	KET-15	241003D02	15 ml
Stop Solution	SCY-15	80958	15 ml

#### Lot Specific Preparations

(Refer to VeriKine™ Human Interferon Alpha 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	13	27	40	53	67	80
	Extended Range 156-5000 pg/ml	5	10	15	20	25	30
<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	20	40	60	80	100	120
	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

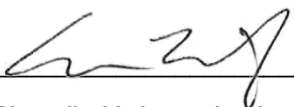
#### Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Spike Recovery: 80-120%

#### Authorization

Released by: 

Date: June 27, 2025

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.