



Certificate of Analysis

VeriKine™ Human Interferon Alpha ELISA Kit

Catalog No: 41100-2

Lot No: 7765

Expiration: December 31, 2027

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP064	K7864	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP022-250	K7693	250 ml
Human IFN Alpha Standard 10,000pg/ml	SMP049-2	K7868	1 vial
Dilution Buffer	SMP021-250	K7694	250 ml
Antibody Concentrate	SMP066-2	K7869	1 vial
HRP Conjugate Concentrate	SMP050-1200	K7870	1 vial
Concentrate Diluent	SMP024-60	K7769	60 ml
TMB Substrate Solution	KET-60	241003D03	60 ml
Stop Solution	SCY-60	80958	60 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
Antibody Concentrate (µl)	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
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	13	27	40	53	67	80
	Extended Range 156-5000 pg/ml	8	16	24	32	40	48
Concentrate Diluent (ml)		2.0	4.0	6.0	8.0	10.0	12.0

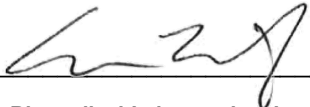
Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Spike Recovery: 80-120%

Authorization

Released by: 

Date: October 20, 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.