



Certificate of Analysis

VeriKine™ Mouse Interferon Alpha ELISA Kit

Catalog No: 42120-1

Lot No: 7247

Expiration: April 30, 2022

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP167	K6162	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K5832	50 ml
Mouse IFN Alpha Standard 10,000 pg/ml	SMP166-1	K6166	1 vial
Sample Buffer	SMP169-60	K6167	50 ml
Antibody Concentrate	SMP170-1	K6168	1 vial
Concentrate Diluent	SMP024-15	K6001	15 ml
HRP Conjugate Concentrate	SMP179-180	K6169	1 vial
Assay Diluent	ASD-15	651172	15 ml
TMB Substrate Solution	KET-15	200210D02	15 ml
Stop Solution	SCY-15	56334	15 ml

Lot Specific Preparations

(Refer to VeriKine™ Mouse Interferon Alpha ELISA Kit Protocol for full instructions.)

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)	16	24	32	40	48	56
Concentrate Diluent (ml)	2	3	4	5	6	7

HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µl)	12	20	24	32	40	48
Assay Diluent (ml)	3	5	6	8	10	12

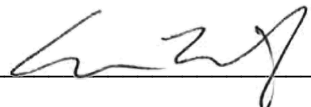
Product Performance Specifications

	Sample Buffer	RPMI (10% FBS)*
Intra-Assay CV	≤ 8%	≤ 10%
Inter-Assay CV	≤ 10%	≤ 10%

Spike Recovery: 115% (range: 82-135%)

* We have noticed variability in the %CV between replicates of individual points on the standard curves prepared in MEM (10% FBS). The range of the variations observed is from 0.3% to 28%. However, these variations do not affect the product's ability to accurately measure Mouse IFN-α independently spiked in MEM (10% FBS).

Authorization

Released by: 

Date: July 28, 2020

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.