



### Certificate of Analysis

#### VeriKine™ Mouse Interferon Alpha ELISA Kit

Catalog No: 42120-1

Lot No:

Expiration:

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP167		1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60		50 ml
Mouse IFN Alpha Standard 10,000 pg/ml	SMP166-1		1 vial
Sample Buffer	SMP169-60		50 ml
Antibody Concentrate	SMP170-1		1 vial
Concentrate Diluent	SMP024-15		15 ml
HRP Conjugate Concentrate	SMP179-180		1 vial
Assay Diluent	ASD-15		15 ml
TMB Substrate Solution	KET-15		15 ml
Stop Solution	SCY-15		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)						
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)						
Assay Diluent (ml)	3	5	6	8	10	12

#### Product Performance Specifications

	Sample Buffer	RPMI (10% FBS)*
Intra-Assay CV	m8%	m10%
Inter-Assay CV	m10%	m10%

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- independently spiked in MEM (10% FBS).

#### Authorization

Released by: \_\_\_\_\_

Date:

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.