

Certificate of Analysis Rev.02

Product: VeriKine™ Mouse IFN-α ELISA Kit

Catalog No: 42120-1

Lot No:

Exp:

Kit Components	Part #	Lot #	Amount
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	SMP179-140		1 vial
Assay Diluent	ASD-15		15 mL
TMB Substrate Solution	NGE-15		15 mL
Stop Solution	SCY-15		15 mL

Lot Specific Preparations

Refer to VeriKine™ Mouse IFN-α 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	Tissue Culture Media (10%FBS)
Intra-Assay CV	≤8%	≤10%
Inter-Assay CV	≤10%	≤10%

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

Released by: _____

Date: _____

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL InterferonSource. 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.