

Certificate of Analysis Rev.04

Product: VeriKine™ Mouse IFN-α ELISA Kit

Catalog No: 42120-2

Lot No: 5496

Exp: April 30, 2013

Kit Components	Part #	Lot #	Amount
Plate (s)	SMP167	K2435	5
Plate Sealers	n/a	n/a	20
Wash Solution Concentrate	SMP022-250	K2446	250 mL
Mouse IFN Alpha Standard 10,000 pg/ml	SMP166-2	K2548	1 vial
Sample Buffer	SMP169-250	K2549	250 mL
Antibody Concentrate	SMP170-2	K2550	1 vial
Concentrate Diluent	SMP024-60	K2365	60 mL
HRP Conjugate Concentrate	SMP179-700	K2551	1 vial
Assay Diluent	ASD-60	883946	60 mL
TMB Substrate Solution	KET-60	110602	60 mL
Stop Solution	SCY-60	23860	60 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)	20	30	40	50	60	70
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)	6	10	12	16	20	24
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%)

Released by: _____

Date: May 16, 2012

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.

*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).