

Certificate of Analysis Rev.03

Product: *VeriKine*<sup>TM</sup> Human IFN- $\alpha$  Serum Sample ELISA Kit

Catalog No: 41110-1

Lot No: 5775

Exp: August 31, 2014

Kit Components	Part #	Lot #	Amount
Plate (s)	SMP047	K2885	1
Plate Sealers	n/a	n/a	4
Wash Solution Concentrate	SMP022-60	K2845	50 mL
Human IFN Alpha Standard 10,000 pg/ml	SMP049-1	K2882	1 vial
Sample Diluent	SMP069-30	K2891	30 mL
Antibody Concentrate	SMP048-1	K2883	1 vial
HRP Conjugate Concentrate	SMP050-240	K2884	1 vial
Concentrate Diluent	SMP024-15	K2876	15 mL
TMB Substrate Solution	KET-15	130514	15 mL
Stop Solution	SCY-15	27059	15 mL
Dilution Buffer	SMP021-30	K2594	30 mL

Lot Specific Preparations

Refer to *VeriKine*<sup>TM</sup> Human IFN Alpha Serum Sample ELISA Kit Protocol for full instructions

Antibody Concentrate Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate ( $\mu$ l)	20	40	60	80	100	120
Dilution Buffer (ml)	2	4	6	8	10	12

HRP Conjugate Concentrate Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate ( $\mu$ l)	16	32	48	64	80	96
Concentrate Diluent (ml)	2	4	6	8	10	12

*Performance Characteristics*

Intra-Assay CV: 8 %  
 Inter-Assay CV: 8 %  
 Spike Recovery: 90%

Released by: 

Date: July 3, 2013

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.