

Certificate of Analysis Rev.05

Product: *Verikine* Mouse IFN-Beta ELISA

Catalog No: 42400-1

Lot No: 5706

Exp: March 31, 2014

Kit Components	Part Number	Lot#	Qty
Plate (s)	SMP028	K2805	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K2160	50 ml
Mouse IFN Beta Standard 500,000pg/ml	SMP210-1	K2808	50
Dilution Buffer	SMP021-30	K2594	30 ml
Antibody Concentrate	SMP211-1	K2809	1 vial
HRP Conjugate Concentrate	SMP179-180	K2810	1vial
Concentrate Diluent	SMP024-30	K2811	30 ml
TMB Substrate Solution	KET-15	120710	15 ml
Stop Solution	SCY-15	24620	15 ml

Lot Specific Preparation

Refer to *Verikine* -Beta ELISA Protocol for full instructions

Antibody Concentrate Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (4	8	12	16	20	24
Concentrate Diluent (ml)	2.0	4.0	6.0	8.0	10.0	12.0

HRP Conjugate Concentrate


Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (20	40	60	80	100	120
Concentrate Diluent (ml)	2.0	4.0	6.0	8.0	10.0	12.0

Performance Characteristics

Intra-Assay CV: ≤ 8 %

Inter-Assay CV: ≤ 8%

Spike Recovery: 94%

Released by: 

Date: March 20, 2013

Sold under license from Pestka Biomedical Laboratory, Inc. d/b/a PBL Interferon Source. 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.