



Certificate of Analysis

VeriKine-HS™ Mouse Interferon Beta Serum ELISA

Catalog No: 42410-1

Lot No: 6614

Expiration: January 31, 2018

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP199	K4397	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60	K4034	2 x 50 ml
Mouse IFN Beta Serum Standard, 10,000 pg/ml	SMP195-1	K4400	1 vial
Sample Diluent	SMP196-30	K4401	25 ml
Antibody Concentrate	SMP197-1	K4402	1 vial
HRP Conjugate Concentrate	SMP056-150	K4405	1 vial
Antibody Diluent	SMP198-15	K4404	15 ml
HRP Diluent	ASDHRP-15	582486	15 ml
Serum Buffer	SMP213-15	K4403	12 ml
TMB Substrate Solution	KET-15	150502	15 ml
Stop Solution	SCY-15	40651	15 ml

Lot Specific Preparations

(Refer to VeriKine-HS™ Mouse Interferon Beta Serum ELISA Protocol for full instructions.)

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)	33	50	67	83	100	117
Antibody Diluent (ml)	2.0	3.0	4.0	5.0	6.0	7.0

HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µl)	27	40	53	67	80	93
HRP Diluent (ml)	2.0	3.0	4.0	5.0	6.0	7.0

Product Performance Characteristics

Intra-Assay CV: ≤ 6%

Inter-Assay CV: ≤ 7%

Average Spike Recovery: 95%

Authorization

Released by: _____

Date: January 17, 2017

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.