



Certificate of Analysis

VeriKine-HS™ Mouse Interferon Beta Serum ELISA

Catalog No: 42410-2

Lot No: 6347

Expiration: October 31, 2016

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP199	K3792	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP057-250	K3553	2 x 250 ml
Mouse IFN Beta Serum Standard, 10,000 pg/ml	SMP195-2	K3795	1 vial
Sample Diluent	SMP196-150	K3796	150 ml
Antibody Concentrate	SMP197-2	K3797	1 vial
HRP Conjugate Concentrate	SMP056-750	K3800	1 vial
Antibody Diluent	SMP198-60	K3799	60 ml
HRP Diluent	ASDHRP-60	369809	60 ml
Serum Buffer	SMP213-60	K3798	60 ml
TMB Substrate Solution	KET-60	140602	60 ml
Stop Solution	SCY-60	31345	60 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)	20	30	40	50	60	70
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: October 13, 2015





pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA | T:+1.732.777.9123 | F:+1.732.777.9141 | E: info@pbl assaysci.com | W: pbl assaysci.com

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.

Rev. 01

