



Certificate of Analysis

VeriKine-HS™ Mouse Interferon Beta Serum ELISA

Catalog No: 42410-2

Lot No:

Expiration:

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP199		5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP057-250		2 x 250 ml
Mouse IFN Beta Serum Standard, 10,000 pg/ml	SMP195-2		1 vial
Sample Diluent	SMP196-150		150 ml
Antibody Concentrate	SMP197-2		1 vial
HRP Conjugate Concentrate	SMP056-750		1 vial
Antibody Diluent	SMP198-60		60 ml
HRP Diluent	ASDHRP-60		60 ml
Serum Buffer	SMP213-60		60 ml
TMB Substrate Solution	KET-60		60 ml
Stop Solution	SCY-60		60 ml

Lot Specific Preparations

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

Antibody Concentrate Preparation

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

HRP Conjugate Concentrate Preparation

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

Product Performance Characteristics

Intra-Assay CV: m6%

Inter-Assay CV: m7%

Average Spike Recovery: 95%

Authorization

Released by: _____

Date:

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.