



Certificate of Analysis

VeriKine-HS™ Mouse Interferon Alpha All Subtype ELISA

Catalog No: 42115-1

Lot No:

Expiration:

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP281		1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60		2 x 50 ml
Mouse IFN Alpha 4 Standard, 10,000 pg/ml	SMP282-1		1 vial
Sample Diluent	SMP283-30		30 ml
Antibody Concentrate	SMP284-1		1 vial
HRP Conjugate Concentrate	SMP056-120		1 vial
Antibody Diluent	SMP285-15		12 ml
HRP Diluent	ASDHRP-15		12 ml
TMB Substrate Solution	KET-15		15 ml
Stop Solution	SCY-15		15 ml

Lot Specific Preparations

(Refer to VeriKine-HS™ Mouse Interferon Alpha All Subtype ELISA Protocol for full instructions.)

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)						
Antibody Diluent (ml)	1.25	2.5	3.5	5.0	6.0	7.0

HRP Solution 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: m8%

Inter-Assay CV: m10%

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.