



Certificate of Analysis

VeriKine™ Human Interferon Beta ELISA Kit

Only for use with 41410-2B Human Interferon Beta ELISA Standard, packaged separately.

Catalog No: 41410-2A

Lot No: 5973

Expiration: February 28, 2015

Kit Components . Store at 2-8°C	Part No.	Lot No.	Quantity
Plate(s)	SMP055	K3056	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP057-250	K3147	250 ml
Sample Diluent	SMP067-150	K3149	150 ml
Antibody Concentrate	SMP062-2	K3150	1 vial
HRP Conjugate Concentrate	SMP056-900	K3151	1 vial
Concentrate Diluent	SMP024-150	K3152	150 ml
TMB Substrate Solution	KET-60	130606	60 ml
Stop Solution	SCY-60	28629	60 ml

Shipped separately on dry ice, refer to 41410-2B lot specific CoA for lot and quantity reference information.

Kit Components . Store at -20°C or below	Part No.	Lot No.	Quantity
Human IFN Beta Standard, 100,000 pg/ml	SMP032-2	N/A	N/A

Lot Specific Preparations

(Refer to VeriKine™ Human Interferon Beta ELISA Kit Protocol for full instructions.)

Antibody Solution Preparation

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

HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µl)	10	20	30	40	50	60
Concentrate Diluent (ml)	2.0	4.0	6.0	8.0	10.0	12.0

Performance Characteristics

Intra-Assay CV: m8%

Inter-Assay CV: m8%

Spike Recovery ~ 90%

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.