



Certificate of Analysis

VeriKine™ Human Interferon Beta ELISA Kit

Only for use with 41410-1B Human Interferon Beta ELISA Standard, packaged separately.

Catalog No: 41410-1A

Lot No: 6279

Expiration: July 31, 2016

Kit Components – Store at 2-8°C	Part No.	Lot No.	Quantity
Plate(s)	SMP055	K3587	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60	K3553	50 ml
Sample Diluent	SMP067-30	K3682	30 ml
Antibody Concentrate	SMP062-1	K3683	1 vial
HRP Conjugate Concentrate	SMP056-180	K3684	1 vial
Concentrate Diluent	SMP024-30	K3339	30 ml
TMB Substrate Solution	KET-15	131026	15 ml
Stop Solution	SCY-15	32954	15 ml

Shipped separately on dry ice, refer to 41410-1B 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	SMP032-1	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

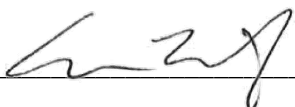
Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Spike Recovery ≥ 90%

Authorization

Released by: 

Date: July 15, 2015

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.



Certificate of Analysis

VeriKine™ Human Interferon Beta ELISA Kit

Only for use with 41410-1A Human Interferon Beta ELISA Kit, packaged separately.

Catalog No: 41410-1B

Lot No: 6279

Expiration: July 31, 2016

Kit Components – Store at -20°C or below	Part No.	Lot No.	Quantity
Human IFN Beta Standard, 100,000 pg/ml	SMP032-1	K3681	3 vials

Shipped separately on wet ice, refer to 41410-1A lot specific CoA for lot and quantity reference information.

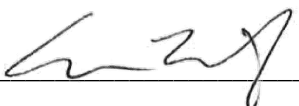
Kit Components – Store at 2-8°C	Part No.	Lot No.	Quantity
Plate(s)	SMP055	N/A	N/A
Plate Sealers	N/A	N/A	N/A
Wash Solution Concentrate	SMP057-60	N/A	N/A
Sample Diluent	SMP067-30	N/A	N/A
Antibody Concentrate	SMP062-1	N/A	N/A
HRP Conjugate Concentrate	SMP056-180	N/A	N/A
Concentrate Diluent	SMP024-30	N/A	N/A
TMB Substrate Solution	KET-15	N/A	N/A
Stop Solution	SCY-15	N/A	N/A

Storage: Human Interferon Beta Standard should be kept at -20°C for one time use only.

Use: Please refer to the VeriKine™ Human Interferon Beta ELISA protocol for full instructions.

Note: This product is provided in two parts, 1A and 1B, due to separate storage conditions. Each part has been validated for use with lot specific counterpart.

Authorization

Released by: _____ 

Date: July 15, 2015

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.