



### 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: 6456

Expiration: April 30, 2017

Kit Components – Store at 2-8°C	Part No.	Lot No.	Quantity
Plate(s)	SMP055	K3996	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP057-250	K3822	250 ml
Sample Diluent	SMP067-150	K4027	150 ml
Antibody Concentrate	SMP062-2	K4028	1 vial
HRP Conjugate Concentrate	SMP056-900	K4029	1 vial
Concentrate Diluent	SMP024-150	K3720	150 ml
TMB Substrate Solution	KET-60	140602	60 ml
Stop Solution	SCY-60	31345	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

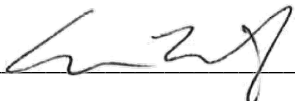
#### Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Spike Recovery ≥ 90%

#### Authorization

Released by: 

Date: April 18, 2016

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 41410-2A Human Interferon Beta ELISA Kit, packaged separately.*

**Catalog No:** 41410-2B

**Lot No:** 6456

**Expiration:** April 30, 2017

Kit Components – Store at -20°C or below	Part No.	Lot No.	Quantity
Human IFN Beta Standard, 100,000 pg/ml	SMP032-2	K4026	15 vials

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

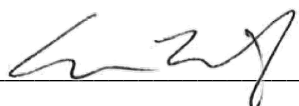
Kit Components – Store at 2-8°C	Part No.	Lot No.	Quantity
Plate(s)	SMP055	N/A	N/A
Plate Sealers	N/A	N/A	N/A
Wash Solution Concentrate	SMP057-250	N/A	N/A
Sample Diluent	SMP067-150	N/A	N/A
Antibody Concentrate	SMP062-2	N/A	N/A
HRP Conjugate Concentrate	SMP056-900	N/A	N/A
Concentrate Diluent	SMP024-150	N/A	N/A
TMB Substrate Solution	KET-60	N/A	N/A
Stop Solution	SCY-60	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, 2A and 2B, due to separate storage conditions. Each part has been validated for use with lot specific counterpart.

**Authorization**

**Released by:** \_\_\_\_\_ 

**Date:** April 18, 2016

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.