



### 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:** 6153

**Expiration:** November 30, 2015

Kit Components – Store at 2-8°C	Part No.	Lot No.	Quantity
Plate(s)	SMP055	K3452	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60	K3147	50 ml
Sample Diluent	SMP067-30	K3456	30 ml
Antibody Concentrate	SMP062-1	K3457	1 vial
HRP Conjugate Concentrate	SMP056-180	K3458	1 vial
Concentrate Diluent	SMP024-30	K3339	30 ml
TMB Substrate Solution	KET-15	131026	15 ml
Stop Solution	SCY-15	30387	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
<b>Antibody Concentrate (µl)</b>	20	40	60	80	100	120
<b>Concentrate Diluent (ml)</b>	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
<b>HRP Conjugate Concentrate (µl)</b>	8	16	24	32	40	48
<b>Concentrate Diluent (ml)</b>	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:** December 4, 2014

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.





# pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA | T:+1.732.777.9123 | F:+1.732.777.9141 | E: [info@pbl assaysci.com](mailto:info@pbl assaysci.com) | W: [pbl assaysci.com](http://pbl assaysci.com)

---

Rev. 07





### 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:** 6153

**Expiration:** November 30, 2015

Kit Components – Store at -20°C or below	Part No.	Lot No.	Quantity
Human IFN Beta Standard, 100,000 pg/ml	SMP032-1	K3455	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:** December 4, 2014

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.





# pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA | T:+1.732.777.9123 | F:+1.732.777.9141 | E: [info@pbl assaysci.com](mailto:info@pbl assaysci.com) | W: [pbl assaysci.com](http://pbl assaysci.com)

---

Rev. 06

