



### Certificate of Analysis

#### VeriKine™ Human Interferon Gamma Receptor 1 ELISA

Catalog No: 41580-1

Lot No:

Expiration:

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP266		1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60		50 ml
Human IFN Gamma R1 Standard, 500,000 pg/ml	SMP267-1		1 vial
Sample Diluent	SMP268-30		25 ml
Antibody Concentrate	SMP269-1		1 vial
Antibody Diluent	SMP270-15		12 ml
HRP Conjugate Concentrate	SMP056-210		1 vial
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™ Human Interferon Gamma Receptor 1 ELISA Protocol for full instructions.)

#### Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)						
Antibody Diluent (ml)	2.0	3.0	4.0	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: ≤ 8%

Inter-Assay CV: ≤ 12%

Average Spike Recovery: 93%

#### 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.





# 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. 00

