



Certificate of Analysis

Protocol A

VeriKine-HS™ Human Interferon Beta Serum ELISA Kit

Certificate of Analysis for **PROTOCOL A** (recommended protocol for improved serum performance)

Catalog No: 41415-1

Lot No:

Expiration:

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP138		1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60		2 x 50 ml
Human IFN Beta Standard, 100,000 pg/ml	SMP146-1		1 vial
Standard Diluent	SMP163-30		25 ml
Sample Buffer	SMP147-15		15 ml
Antibody Concentrate	SMP148-1		1 vial
HRP Conjugate Concentrate	SMP056-320		1 vial
Assay Diluent	ASD-30		25 ml
Diluent Additive III	SMP209-2		ml
TMB Substrate Solution	KET-15		15 ml
Stop Solution	SCY-15		15 ml

Lot Specific Preparations

(Refer to VeriKine-HS™ Human Interferon Beta Serum ELISA Kit **Protocol A** for full instructions.)

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)						
Assay Diluent (ml)						
Diluent Additive (III)						
Total Volume (Assay Diluent + Diluent Additive III) (ml)	2	3	4	5	6	7

HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µl)						
Assay Diluent (ml)	3	5	6	8	10	12

Product Performance Specifications

	Standard Diluent	Human Serum	Tissue Culture Media (10% FBS)
Intra-Assay CV	≤ 8%	≤ 6%	≤ 8%
Inter-Assay CV	≤ 10%	≤ 10%	≤ 10%

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.