

**Certificate of Analysis** Rev.04

**Product:** VeriKine-HS™ Human IFN-β Serum ELISA Kit

**Catalog No:** 41415-1

**Lot No:**

**Exp:**

Kit Components	Part #	Lot #	Amount
Plate (s)	SMP138		1
Plate Sealers	n/a	n/a	4
Wash Solution Concentrate	SMP057-60		2 x 50 mL
Human IFN Beta Standard 100,000pg/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
TMB Substrate Solution	KET-15		15 mL
Stop Solution	SCY-15		15 mL

**Lot Specific Preparations**

Refer to VeriKine-HS™ Human IFN-β Serum ELISA Kit ELISA Kit Protocol for full instructions

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (μl)						
Assay Diluent (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 %

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

Date:

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL InterferonSource. 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.