



**Certificate of Analysis**

**Protocol B**

**VeriKine-HS™ Human Interferon Beta Serum ELISA Kit**  
*Certificate of Analysis for **PROTOCOL B***

**Catalog No:** 41415-1

**Lot No:** 7292

**Expiration:** October 31, 2021

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP138	K6286	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60	K6285	2 x 50 ml
Human IFN Beta Standard, 100,000 pg/ml	SMP146-1	K6289	1 vial
Standard Diluent	SMP163-30	K6293	25 ml
Sample Buffer	SMP147-15	K6290	15 ml
Antibody Concentrate	SMP148-1	K6291	1 vial
HRP Conjugate Concentrate	SMP056-320	K6292	1 vial
Assay Diluent	ASD-30	795154	25 ml
TMB Substrate Solution	KET-15	200210D02	15 ml
Stop Solution	SCY-15	56334	15 ml

**Lot Specific Preparations**

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

**Antibody Solution Preparation**

Micro-plate Strips Used	2	4	6	8	10	12
<b>Antibody Concentrate (µl)</b>	16	24	32	40	48	56
<b>Assay Diluent (ml)</b>	2	3	4	5	6	7

**HRP Solution Preparation**

Micro-plate Strips Used	2	4	6	8	10	12
<b>HRP Conjugate Concentrate (µl)</b>	17	29	34	46	57	69
<b>Assay Diluent (ml)</b>	3	5	6	8	10	12

**Product Performance Specifications**

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

**Authorization**

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

**Date:** November 16, 2020

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.