



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: 7464

Expiration: November 30, 2022

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP138	K6732	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60	K6742	2 x 50 ml
Human IFN Beta Standard, 100,000 pg/ml	SMP146-1	K6735	1 vial
Standard Diluent	SMP163-30	K6739	25 ml
Sample Buffer	SMP147-15	K6736	15 ml
Antibody Concentrate	SMP148-1	K6737	1 vial
HRP Conjugate Concentrate	SMP056-320	K6738	1 vial
Assay Diluent	ASD-30	157923	25 ml
TMB Substrate Solution	KET-15	210201D01	15 ml
Stop Solution	SCY-15	59172	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
Antibody Concentrate (µl)	7	10	13	17	20	23
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)	15	25	30	40	50	60
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: December 7, 2021

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.