



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: 7101

Expiration: July 15, 2020

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP138	K5696	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60	K5493	2 x 50 ml
Human IFN Beta Standard, 100,000 pg/ml	SMP146-1	K5699	1 vial
Standard Diluent	SMP163-30	K5704	25 ml
Sample Buffer	SMP147-15	K5701	15 ml
Antibody Concentrate	SMP148-1	K5702	1 vial
HRP Conjugate Concentrate	SMP056-320	K5703	1 vial
Assay Diluent	ASD-30	314264	25 ml
TMB Substrate Solution	KET-15	180803D02	15 ml
Stop Solution	SCY-15	51066	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)	16	24	32	40	48	56
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)	17	29	34	46	57	69
Assay Diluent (ml)	3	5	6	8	10	12

Product Performance Specifications

	Standard Diluent	Human Serum	Tissue Culture Media (10% FBS)
Intra-Assay CV	m8%	m6%	m8%
Inter-Assay CV	m10%	m10%	m10%

Authorization

Released by: _____

Date: August 13, 2019

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.