



Certificate of Analysis

VeriKine™ Mouse Interferon Beta ELISA Kit

Catalog No: 42400-1

Lot No: 7546

Expiration: September 30, 2023

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP028	K7042	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K6719	50 ml
Mouse IFN Beta Standard, 500,000 pg/ml	SMP210-1	K7045	1 vial
Dilution Buffer	SMP021-30	K6965	30 ml
Antibody Concentrate	SMP211-1	K7047	1 vial
HRP Conjugate Concentrate	SMP179-240	K7048	1 vial
Concentrate Diluent	SMP024-30	K6929	30 ml
TMB Substrate Solution	KET-15	210201D01	15 ml
Stop Solution	SCY-15	67088	15 ml

Lot Specific Preparations

(Refer to VeriKine™ Mouse Interferon Beta ELISA Kit Protocol for full instructions.)

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)	40	80	120	160	200	240
Concentrate Diluent (ml)	2	4	6	8	10	12

HRP Solution Preparation

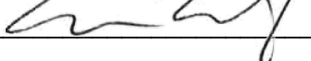
Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µl)	16	32	48	64	80	96
Concentrate Diluent (ml)	2	4	6	8	10	12

Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Authorization

Released by: 

Date: September 15, 2022

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.