



Certificate of Analysis

VeriKine™ Mouse Interferon Beta ELISA Kit

Catalog No: 42400-1
Lot No: 7321

Expiration: January 31, 2022

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP028	K6362	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K6226	50 ml
Mouse IFN Beta Standard, 500,000 pg/ml	SMP210-1	K6367	1 vial
Dilution Buffer	SMP021-30	K6368	30 ml
Antibody Concentrate	SMP211-1	K6369	1 vial
HRP Conjugate Concentrate	SMP179-240	K6370	1 vial
Concentrate Diluent	SMP024-30	K6371	30 ml
TMB Substrate Solution	KET-15	200210D02	15 ml
Stop Solution	SCY-15	56334	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)	10	20	30	40	50	60
Concentrate Diluent (ml)	2	4	6	8	10	12

Performance Characteristics

Intra-Assay CV: ≤ 8%
Inter-Assay CV: ≤ 8%

Authorization

Released by: _____ 

Date: January 8, 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.