



### Certificate of Analysis

#### VeriKine™ Mouse Interferon Beta ELISA Kit

Catalog No: 42400-2

Lot No: 7250

Expiration: July 31, 2021

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP028	K6191	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP022-250	K5510	250 ml
Mouse IFN Beta Standard, 500,000 pg/ml	SMP210-2	K6194	1 vial
Dilution Buffer	SMP021-150	K5891	150 ml
Antibody Concentrate	SMP211-2	K6195	1 vial
HRP Conjugate Concentrate	SMP179-1200	K6196	1 vial
Concentrate Diluent	SMP024-150	K6001	150 ml
TMB Substrate Solution	KET-60	180803D03	60 ml
Stop Solution	SCY-60	51066	60 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: July 14, 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.