



### Certificate of Analysis

#### VeriKine™ Mouse Interferon Beta ELISA Kit

Catalog No: 42400-2

Lot No: 7499

Expiration: April 30, 2023

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP028	K6898	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP022-250	K6596	250 ml
Mouse IFN Beta Standard, 500,000 pg/ml	SMP210-2	K6901	1 vial
Dilution Buffer	SMP021-150	K6902	150 ml
Antibody Concentrate	SMP211-2	K6903	1 vial
HRP Conjugate Concentrate	SMP179-1200	K6904	1 vial
Concentrate Diluent	SMP024-150	K6731	150 ml
TMB Substrate Solution	KET-60	210201D02	60 ml
Stop Solution	SCY-60	63243	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
<b>Antibody Concentrate (µl)</b>	40	80	120	160	200	240
<b>Concentrate Diluent (ml)</b>	2	4	6	8	10	12

##### HRP Solution Preparation

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

#### Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

#### Authorization

Released by: \_\_\_\_\_

Date: May 4, 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.