



### Certificate of Analysis

#### VeriKine-HS™ Mouse Interferon Alpha All Subtype ELISA

Catalog No: 42115-1

Lot No: 7209

Expiration: April 30, 2021

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP281	K5982	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K5832	2 x 50 ml
Mouse IFN Alpha 4 Standard, 10,000 pg/ml	SMP282-1	K5985	1 vial
Sample Diluent	SMP283-30	K6137	30 ml
Antibody Concentrate	SMP284-1	K5987	1 vial
HRP Conjugate Concentrate	SMP056-120	K5988	1 vial
Antibody Diluent	SMP285-15	K5989	12 ml
HRP Diluent	ASDHRP-15	782465	12 ml
TMB Substrate Solution	KET-15	200210D02	15 ml
Stop Solution	SCY-15	56334	15 ml

#### Lot Specific Preparations

(Refer to VeriKine-HS™ Mouse Interferon Alpha All Subtype ELISA Protocol for full instructions.)

#### Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)	100	200	280	400	480	560
Antibody Diluent (ml)	1.25	2.5	3.5	5.0	6.0	7.0

#### HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µl)	10	15	20	25	30	35
HRP Diluent (ml)	2.0	3.0	4.0	5.0	6.0	7.0

#### Product Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 10%

#### Authorization

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

Date: June 12, 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.