



### Certificate of Analysis

#### VeriKine-HS™ Human Interferon Alpha All Subtype TCM ELISA Kit

Catalog No: 41135-1

Lot No: 7090

Expiration: May 31, 2020

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP217	K5634	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K5454	2 x 50 ml
Human IFN Alpha Standard 10,000 pg/ml	SMP049-400	K5622	1 vial
Assay Buffer	SMP324-8	K5623	8 ml
Standard Diluent	SMP323-60	K5624	55 ml
Antibody Concentrate	SMP322-1	K5626	1 vial
Antibody Diluent	SMP315-15	K5625	15 ml
HRP Conjugate Concentrate	SMP056-240	K5627	1 vial
HRP Diluent	ASDHRP-15	889225	15 ml
TMB Substrate Solution	KET-15	180803D02	15 ml
Stop Solution	SCY-15	51066	15 ml

#### Lot Specific Preparation

(Refer to VeriKine-HS™ Human IFN  $\alpha$  All Subtype TCM ELISA Kit Protocol for full instructions.)

#### Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate ( $\mu$ l)	20	40	60	80	100	120
Antibody Diluent (ml)	2	4	6	8	10	12

#### HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate ( $\mu$ l)	9	18	27	36	44	53
HRP Diluent (ml)	2	4	6	8	10	12

#### Performance Characteristics

Intra-Assay CV:  $\leq$  8%

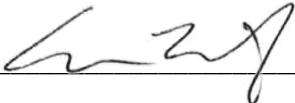
Inter-Assay CV:  $\leq$  10%

Intermediate Precision\* CV:  $\leq$  14%

\*(25 assays performed by 10 operators using 3 lots of components)

Average Spike Recovery: 105%

#### Authorization

Released by: 

Date: September 5, 2019

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.