



### Certificate of Analysis

#### VeriKine™ Cynomolgus/Rhesus Interferon Alpha Serum ELISA Kit

Catalog No: 46100-1

Lot No: 7185

Expiration: November 30, 2020

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP135	K5934	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K5832	50 ml
Cyno/Rhesus IFN Alpha 2 Standard, 10,000 pg/ml	SMP141-1	K5937	1 vial
Standard Diluent	SMP163-30	K5938	25 ml
Sample Buffer	SMP153-8	K5939	8 ml
Antibody Concentrate	SMP154-1	K5963	1 vial
HRP Conjugate Concentrate	SMP056-330	K5942	1 vial
Concentrate Diluent	SMP024-30	K5962	30 ml
TMB Substrate Solution	KET-15	180803D02	15 ml
Stop Solution	SCY-15	51066	15 ml

#### Lot Specific Preparations

(Refer to VeriKine™ Cynomolgus/Rhesus Interferon Alpha Serum ELISA Kit Protocol for full instructions.)

##### Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
<b>Antibody Concentrate (µl)</b>	20	40	60	80	100	120
<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>	10	20	30	40	50	60
<b>Concentrate Diluent (ml)</b>	2	4	6	8	10	12

##### Product Performance Specifications

	Standard Diluent	Cynomolgus Monkey Serum	Tissue Culture Media (10% FBS)
<b>Intra-Assay CV</b>	≤ 8%	≤ 8%	≤ 8%
<b>Inter-Assay CV</b>	≤ 10%	≤ 15%	≤ 10%

#### Authorization

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

Date: March 4, 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.