



Certificate of Analysis

VeriKine™ Cynomolgus/Rhesus Interferon Alpha Serum ELISA Kit

Catalog No: 46100-1

Lot No: 6549

Expiration: May 31, 2017

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP135	K4256	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K4249	50 ml
Cyno/Rhesus IFN Alpha 2 Standard, 10,000 pg/ml	SMP141-1	K4259	1 vial
Standard Diluent	SMP163-30	K4260	25 ml
Sample Buffer	SMP153-8	K4261	8 ml
Antibody Concentrate	SMP154-1	K4262	1 vial
HRP Conjugate Concentrate	SMP056-330	K4263	1 vial
Concentrate Diluent	SMP024-30	K4264	30 ml
TMB Substrate Solution	KET-15	150502	15 ml
Stop Solution	SCY-15	37395	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
Antibody Concentrate (µl)	20	40	60	80	100	120
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)	7	13	20	27	33	40
Concentrate Diluent (ml)	2	4	6	8	10	12

Product Performance Specifications

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

Authorization

Released by: _____

Date: September 9, 2016

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.