

Certificate of Analysis Rev.05

Product: VeriKine™ Cynomolgus/Rhesus IFN-α Serum ELISA Kit

Catalog No: 46100-1

Lot No:

Exp:

Kit Components	Part Number	Lot #	Volume
Plate (s)	SMP135		1
Plate Sealers	n/a	n/a	4
Wash Solution Concentrate	SMP022-60		50 mL
Rhesus/Cyno IFN Alpha 2 Standard 10,000pg/ml	SMP141-1		0.5 mL
Standard Diluent	SMP163-30		25 mL
Sample Buffer	SMP153-8		8 mL
Antibody Concentrate	SMP154-1		1 vial
HRP Conjugate Concentrate	SMP056-330		1 vial
Assay Diluent	SMP024-30		30 mL
TMB Substrate Solution	KET-15		15 mL
Stop Solution	SCY-15		15 mL

Lot Specific Preparations

Refer to VeriKine™ Cynomolgus/Rhesus IFN-α Serum ELISA Kit Protocol for full instructions

Antibody Concentrate Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)						
Assay Diluent (ml)	2	4	6	8	10	12

HRP Conjugate Concentrate Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µl)						
Assay Diluent (ml)	2	4	6	8	10	12

Product Performance Specifications *not lot specific

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

Released by: _____

Date:

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL InterferonSource. 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.