



Certificate of Analysis

VeriKine™ Cynomolgus Interferon Beta ELISA Kit

Catalog No: 46415-1

Lot No: 7102

Expiration: July 31, 2020

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP138	K5696	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60	K5493	2 x 50 ml
Cyno IFN Beta Standard, 100,000 pg/ml	SMP261-1	K5775	1 vial
Standard Diluent	SMP163-30	K5704	25 ml
Sample Buffer	SMP147-15	K5701	15 ml
Antibody Concentrate	SMP148-100	K5702	1 vial
HRP Conjugate Concentrate	SMP056-450	K5703	1 vial
Assay Diluent	ASD-30	314264	25 ml
TMB Substrate Solution	KET-15	180803D02	15 ml
Stop Solution	SCY-15	51066	15 ml

Lot Specific Preparations

(Refer to VeriKine™ Cynomolgus Interferon Beta ELISA Kit Protocol for full instructions.)

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)	16	24	32	40	48	56
Assay Diluent (ml)	2	3	4	5	6	7

HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (µl)	20	33	40	53	67	80
Assay Diluent (ml)	3	5	6	8	10	12

Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 10%

Average Spike Recovery: 94%

Authorization

Released by: _____

Date: October 9, 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.