



Certificate of Analysis

VeriKine™ Human Interferon Alpha/Beta Receptor 2 ELISA Kit (custom order)

Catalog No: 41385-1

Lot No: 7402

Expiration: February 28, 2022

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP225	K6600	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K6446	2 x 50 ml
Human IFN Alpha/Beta Receptor 2 (IFNAR2) Standard 10,000 pg/ml	SMP228-1	K6604	1 vial
Sample Diluent	SMP227-60	K6605	60 ml
Antibody Concentrate	SMP229-1	K6606	1 vial
HRP Conjugate Concentrate	SMP056-240	K6607	1 vial
Concentrate Diluent	SMP024-30	K6611	30 ml
Sample Buffer	SMP226-8	K6608	8 ml
TMB Substrate Solution	KET-15	200210D02	15 ml
Stop Solution	SCY-15	59172	15 ml

Lot Specific Preparation

(Refer to VeriKine™ Human IFN- α/β Receptor 2 ELISA Kit Protocol for full instructions.

***For this lot, please refer to updated Standard Curve Preparation Figure 1 on the next page)**

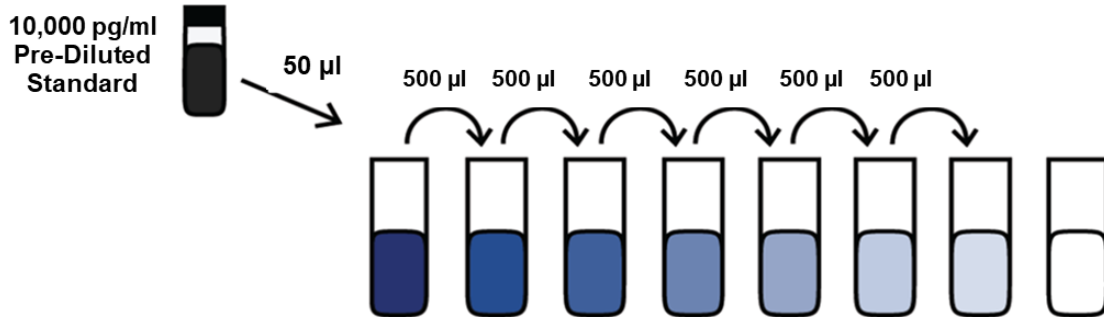
Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (μl)	17	33	50	67	83	100
Concentrate Diluent (ml)	2.0	4.0	6.0	8.0	10.0	12.0

HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
HRP Conjugate Concentrate (μl)	10	20	30	40	50	60
Concentrate Diluent (ml)	2.0	4.0	6.0	8.0	10.0	12.0

***Update to Standard Curve Preparation “Figure 1: 7-Point Standard Curve Prepared in Sample Diluent” below:**



Sample Diluent Label	S7	S6	S5	S4	S3	S2	S1	Blank
Sample Diluent (µl)	950	500	500	500	500	500	500	500
IFNAR2 Conc. (pg/ml)	500	250	125	62.5	31.25	15.63	7.81	0

Authorization

Released by: _____

Date: September 8, 2021

Sold under license from Pestka Biomedical Laboratory, 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.