



Certificate of Analysis

VeriKine™ Human Interferon Beta ELISA Kit

Catalog No: 41410-2

Lot No: 7274

Expiration: May 31, 2022

Kit Components – Store at 2-8°C	Part No.	Lot No.	Quantity
Plate(s)	SMP055	K6243	5
Plate Sealers	N/A	N/A	20
Wash Solution Concentrate	SMP057-250	K6246	250 ml
Human IFN Beta Standard, 100,000 pg/ml	SMP146-2	K6247	1 vial
Sample Diluent	SMP067-150	K6248	150 ml
Antibody Concentrate	SMP062-2	K6250	1 vial
HRP Conjugate Concentrate	SMP056-900	K6251	1 vial
Concentrate Diluent	SMP024-150	K6001	150 ml
TMB Substrate Solution	KET-60	180803D03	60 ml
Stop Solution	SCY-60	56334	60 ml

Lot Specific Preparations

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

Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
Antibody Concentrate (µl)	5	10	15	20	25	30
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)	5	10	15	20	25	30
Concentrate Diluent (ml)	2.0	4.0	6.0	8.0	10.0	12.0

Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 8%

Authorization

Released by: 

Date: September 1, 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.