



Certificate of Analysis

VeriKine™ Human Interferon Beta ELISA Kit

Catalog No: 41410-1

Lot No:

Expiration:

Kit Components . Store at 2-8°C	Part No.	Lot No.	Quantity
Plate(s)	SMP055		1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP057-60		50 ml
Human IFN Beta Standard, 100,000 pg/ml	SMP146-1		1 vial
Sample Diluent	SMP067-30		30 ml
Antibody Concentrate	SMP062-1		1 vial
HRP Conjugate Concentrate	SMP056-180		1 vial
Concentrate Diluent	SMP024-30		30 ml
TMB Substrate Solution	KET-15		15 ml
Stop Solution	SCY-15		15 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)						
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)						
Concentrate Diluent (ml)	2.0	4.0	6.0	8.0	10.0	12.0

Performance Characteristics

Intra-Assay CV: m8%

Inter-Assay CV: m8%

Authorization

Released by: _____

Date:

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.