



Certificate of Analysis

VeriPlex™ Human Interferon 9-Plex ELISA Kit

Catalog No: 51500-1

Lot No:

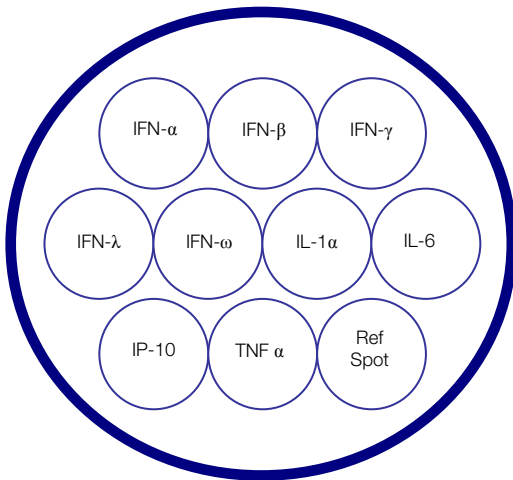
Expiration:

Software Code (input as product number in Q-View™ software):

Kit Components	Part No.	Lot No.	Quantity	Reconstitution Instructions
Plate(s)	311952HU		1	N/A
Plate Sealers	N/A	N/A	3	N/A
Wash Solution Concentrate	101158GR		1 bottle	950 ml DI Water
Human IFN Multiplex Antigen Standard	311955HU		1 vial	ml Sample Diluent
Human IFN Multiplex Detection Mix	311956HU		1 vial	ml DI Water
HRP Concentrate	101173GR		6 ml	N/A
Sample Diluent	311969HU		10 ml	N/A
Substrate A	101159GR		3 ml	Mix equal parts of A and B together and store in the dark
Substrate B	101160GR		3 ml	
Diluent Additive II	SMP208-1		1 vial	N/A
Diluent Additive III	SMP209-3		1 vial	N/A

Kit Components	Part No.	Lot No.	Quantity	Reconstitution Instructions
Assay Diluent	311983HU		1 vial	ml DI Water + ml Diluent Additive II + ml Diluent Additive III to reconstitute Assay Diluent

Spot Location in each well:



Assay Range for Analytes:

Analyte	Assigned High Value (pg/ml)	Assay Range (pg/ml)	LLOD (pg/ml)	LLOQ (pg/ml)
IFN-α		-		
IFN-β		-		
IFN-γ		-		
IFN-λ		-		
IFN-ω		-		
IFN-1α		-		
IL-6		-		
IP-10		-		
TNF-α		-		

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.