

Certificate of Analysis

VeriPlex™ Human Interferon 9-Plex ELISA Kit

Catalog No: 51500-1

Lot No: 7404

Expiration: July 5, 2022

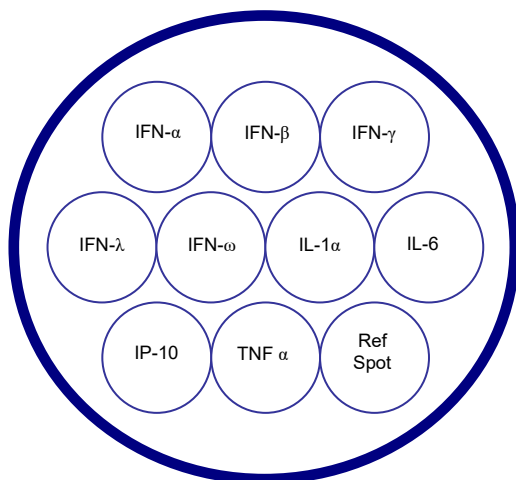
Software Code (input as product number in Q-View™ software): PCHIM180801-PB

Kit Components	Part No.	Lot No.	Quantity	Reconstitution Instructions
Plate(s)	311952HU	PCHIP210519	1	N/A
Plate Sealers	N/A	N/A	3	N/A
Wash Solution Concentrate	101158GR	PCHW210428	1 bottle	950 ml DI Water
Human IFN Multiplex Antigen Standard	311955HU	PCHIM180801	1 vial	0.600 ml Sample Diluent
Human IFN Multiplex Detection Mix	311956HU	PCHID210527	1 vial	6 ml DI Water
HRP Concentrate	101173GR	PCHN210105	6 ml	N/A
Sample Diluent	311969HU	PCHIS210407	10 ml	N/A
Substrate A	101159GR	PCHA210407	3 ml	Mix equal parts of A and B together and store in the dark
Substrate B	101160GR	PCHB210407	3 ml	
Diluent Additive II	SMP208-1	DA078	1 vial	N/A
Diluent Additive III	SMP209-3	DA079	1 vial	N/A

Kit Components	Part No.	Lot No.	Quantity	Reconstitution Instructions
Assay Diluent	311983HU	PCHV210517	1 vial	5.0 ml DI Water + 0.480 ml Diluent Additive II + 0.518 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	3400	3400 – 4.7	0.0	4.4
IFN-β	20100	20100 – 27.6	8.6	31.4
IFN-γ	5600	5600 – 7.7	0.1	8.6
IFN-λ	12900	12900 – 17.7	7.9	94.9
IFN-ω	3800	3800 – 5.2	0.3	5.1
IL-1α	1100	1100 – 1.5	0.3	2.6
IL-6	22000	22000 – 30.2	0.0	65.5
IP-10	1400	1400 – 1.9	0.3	6.1
TNF-α	2900	2900 – 4.0	0.0	4.1

Authorization

Released by: _____

Date: June 29, 2021

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.