

Certificate of Analysis

VeriPlex™ Human Interferon 9-Plex ELISA Kit

Catalog No: 51500-1

Lot No: 7198

Expiration: June 30, 2021

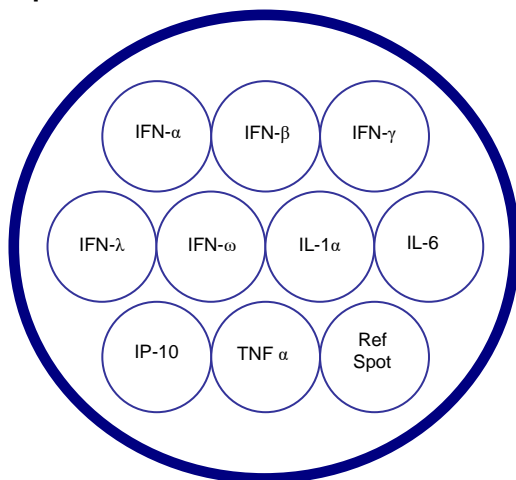
Software Code (input as product number in Q-View™ software): PCHIM180801-PB

Kit Components	Part No.	Lot No.	Quantity	Reconstitution Instructions
Plate(s)	311952HU	PCHIP200220	1	N/A
Plate Sealers	N/A	N/A	3	N/A
Wash Solution Concentrate	101158GR	PCHW200210	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	PCHID200324	1 vial	6 ml DI Water
HRP Concentrate	101173GR	PCHN191231	6 ml	N/A
Sample Diluent	311969HU	PCHIS191210	10 ml	N/A
Substrate A	101159GR	PCHA190911	3 ml	Mix equal parts of A and B together and store in the dark
Substrate B	101160GR	PCHB190911	3 ml	
Diluent Additive II	SMP208-1	DA072	1 vial	N/A
Diluent Additive III	SMP209-3	DA073	1 vial	N/A

Kit Components	Part No.	Lot No.	Quantity	Reconstitution Instructions
Assay Diluent	311983HU	PCHV200304	1 vial	5.0 ml DI Water + 0.480 ml Diluent Additive II + 0.518 ml Diluent Additive III to reconstitute Assay Diluent

***Please note that Assay Diluent contents may not fully dissolve after reconstitution and precipitate may be observed. Performance should not be affected. Continue to use component as directed within the protocol.**

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	3700	3700 – 5.1	1.4	13.8
IFN-β	20700	20700 – 28.4	14.1	28.2
IFN-γ	5800	5800 – 8.0	1.4	8.5
IFN-λ	16600	16600 – 22.8	14.0	119.0
IFN-ω	4000	4000 – 5.5	5.5	10.2
IL-1α	1100	1100 – 1.5	0.4	4.1
IL-6	20600	20600 – 28.3	13.4	80.6
IP-10	1600	1600 – 2.2	0.7	6.5
TNF-α	2800	2800 – 3.8	3.7	4.1

Authorization

Released by: _____

Date: April 29, 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.

Rev. 10