

Certificate of Analysis

VeriPlex™ Mouse Cytokine 9-Plex ELISA Kit

Catalog No: 52500-1

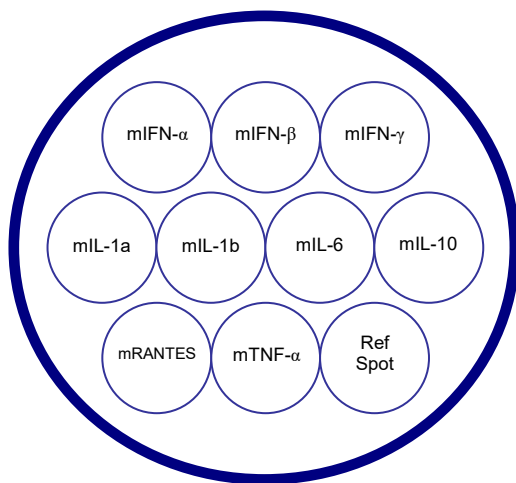
Lot No: 7316

Expiration: November 30, 2021

Software Code (input as product number in Q-View™ software): MPBM170817CY

Kit Components	Part No.	Lot No.	Quantity	Reconstitution Instructions
Plate(s)	491952MS	MPBP201118	1	N/A
Plate Sealers	N/A	N/A	3	N/A
Wash Solution Concentrate	101158GR	PCHW201012	1 bottle	950 ml DI Water
Mouse Cytokine 9-Plex Antigen Standard	491955MS	MPBM170817	1 vial	0.550 ml Sample Diluent
Mouse Cytokine 9-Plex Detection Mix	491956MS	MPBD200306	1 vial	N/A
HRP Concentrate	101173GR	PCHN200909	6 ml	N/A
Mouse Cytokine 9-Plex Sample Diluent	491957MS	MPBS200224	10 ml	N/A
Substrate A	101159GR	PCHA200610	3 ml	Mix equal parts of A and B together and store in the dark
Substrate B+	101120GR	PCHB+200610	3 ml	

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)
mIFN-α	1100	1100 – 1.5	1.0	8.2
mIFN-β	9650	9650 – 13.2	5.3	12.3
mIFN-γ	1000	1000 – 1.4	0.2	4.2
mIL-1a	3500	3500 – 4.8	0.0	4.8
mIL-1b	4500	4500 – 6.2	0.0	6.2
mIL-6	2000	2000 – 2.7	6.6	2.8
mIL-10	1400	1400 – 1.9	8.2	11.0
mRANTES	1450	1450 – 2.0	0.0	2.0
mTNF-α	300	300 – 0.4	0.0	0.4

Authorization

Released by: _____

Date: December 14, 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.