

Certificate of Analysis

VeriPlex™ Mouse Cytokine 9-Plex ELISA Kit

Catalog No: 52500-1

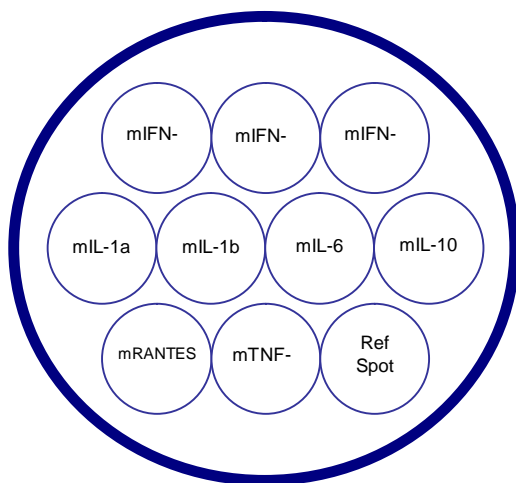
Lot No:

Expiration:

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

Kit Components	Part No.	Lot No.	Quantity	Reconstitution Instructions
Plate(s)	491952MS		1	N/A
Plate Sealers	N/A	N/A	3	N/A
Wash Solution Concentrate	101158GR		1 bottle	ml DI Water
Mouse Cytokine 9-Plex Antigen Standard	491955MS		1 vial	ml Sample Diluent
Mouse Cytokine 9-Plex Detection Mix	491956MS		1 vial	N/A
HRP Concentrate	101173GR		6 ml	N/A
Mouse Cytokine 9-Plex Sample Diluent	491957MS		10 ml	N/A
Substrate A	101159GR		3 ml	Mix equal parts of A and B together and store in the dark
Substrate B+	101120GR		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-		.		
mIFN-		.		
mIFN-		.		
mL-1a		.		
mL-1b		.		
mL-6		.		
mL-10		.		
mRANTES		.		
mTNF-		.		

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.