

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse MCP-1 Kit
Kit Catalog Numbers: K152NN-Series
Kit Lot Number: K0080958
Expiration Date: 30 Nov 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms MCP-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms MCP-1	4.42	75% - 125%	< 20%	325	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms MCP-1	Calibrator	482	pg/mL
		Cytokine Panel 1 (mouse) Control 1	489	pg/mL
		Cytokine Panel 1 (mouse) Control 2	149	pg/mL
		Cytokine Panel 1 (mouse) Control 3	64.9	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms MCP-1 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
482	437,000	3.2
121	55,900	4.4
30.1	7,250	5.2
7.53	1,200	7.4
1.88	373	4.0
0.471	154	1.9
0.118	108	4.1
0.00	93	4.3

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse MCP-1 Plate	Z0046760	2-8°C	31 Jul 2020
Anti-ms MCP-1 Antibody	D0080971	2-8°C	30 Nov 2019
Cytokine Panel 1 (mouse) Calibrator Blend	A0080135	2-8°C	30 Jun 2022
Diluent 41	M0410021	≤ -10°C	30 Jun 2020
Diluent 45	M0450020	≤ -10°C	31 May 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Cytokine Panel 1 (mouse) Control 1	Not Kit Specific
Cytokine Panel 1 (mouse) Control 2	Not Kit Specific
Cytokine Panel 1 (mouse) Control 3	Not Kit Specific

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		13 MAR 2018