

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: K0080736
Expiration Date: 31 Aug 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel (mouse) Control 1		Cytokine Panel (mouse) Control 2		Cytokine Panel (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	491	pg/mL
		Cytokine Panel (mouse) Control 1	479	pg/mL
		Cytokine Panel (mouse) Control 2	153	pg/mL
		Cytokine Panel (mouse) Control 3	55.3	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
491	250,000	4.2
123	28,700	5.1
30.7	4,050	5.7
7.67	789	5.6
1.92	259	5.7
0.479	121	4.9
0.120	89	6.0
0.00	76	4.7



CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse MCP-1 Plate	Z0046397	2-8°C	31 Aug 2019
Anti-ms MCP-1 Antibody	D0080940	2-8°C	31 Aug 2019
Cytokine Panel 1 (mouse) Calibrator Blend	A0080125	2-8°C	31 Mar 2022
Diluent 41	M0410016	≤ -10°C	31 Aug 2019
Diluent 45	M0450016	≤ -10°C	30 Sep 2019
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 (mouse) Control 1	Not Kit Specific
Cytokine Panel (mouse) Control 2	Not Kit Specific
Cytokine Panel (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		23 JUN 2017