

## 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:** K0081369  
**Expiration Date:** 31 Jul 2020

**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	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	MCP-1	4.42	75% - 125%	< 20%	325	75% - 125%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	MCP-1	Calibrator	482	pg/mL
		Cytokine Panel 1 (mouse) Control 1	443	pg/mL
		Cytokine Panel 1 (mouse) Control 2	135	pg/mL
		Cytokine Panel 1 (mouse) Control 3	56	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

MCP-1 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
482	387,000	3.0
121	46,800	6.1
30.1	6,010	5.4
7.53	1,140	5.1
1.88	342	3.0
0.471	152	4.2
0.118	106	8.9
0.00	91	8.4

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse MCP-1 Plate	Z0046760	2-8°C	31 Jul 2020
Anti-ms MCP-1 Antibody	D0081214	2-8°C	31 Jul 2022
Cytokine Panel 1 (mouse) Calibrator Blend	A0080135	2-8°C	30 Jun 2022
Diluent 41	M0410029	≤ -10°C	31 Dec 2021
Diluent 45	M0450030	≤ -10°C	28 Feb 2022
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
<b>Review/Approval</b>	Karen Hamilla	Quality		05 DEC 2019