

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: K0081493
Expiration Date: 31 Mar 2022

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	447	pg/mL
		Cytokine Panel 1 (mouse) Control 1	439	pg/mL
		Cytokine Panel 1 (mouse) Control 2	133	pg/mL
		Cytokine Panel 1 (mouse) Control 3	55.3	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
447	283,000	3.4
112	38,600	4.9
27.9	6,550	3.6
6.98	1,420	4.2
1.75	444	3.7
0.437	185	2.7
0.109	126	6.3
0.00	104	4.0



CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse MCP-1 Plate	Z0047336	2-8°C	30 Nov 2022
Anti-ms MCP-1 Antibody	D0081498	2-8°C	30 Nov 2023
Cytokine Panel 1 (mouse) Calibrator Blend	A0080218	2-8°C	31 Oct 2024
Diluent 41	M0410030	≤ -10°C	31 Mar 2022
Diluent 45	M0450031	≤ -10°C	31 Mar 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
Review/Approval	Karen Hamilla	Quality		14 JUL 2020