

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: K0081402
Expiration Date: 28 Feb 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	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
447	467,000	4.0
112	61,200	7.0
27.9	9,470	3.1
6.98	1,900	3.0
1.75	567	4.3
0.437	221	5.3
0.109	137	3.8
0.00	107	3.7

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse MCP-1 Plate	Z0047246	2-8°C	30 Jun 2022
Anti-ms MCP-1 Antibody	D0081326	2-8°C	31 Mar 2023
Cytokine Panel 1 (mouse) Calibrator Blend	A0080218	2-8°C	31 Oct 2024
Diluent 41	M0410030	≤ -10°C	31 Mar 2022
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
Review/Approval	Karen Hamilla	Quality		26 MAR 2020