

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human MCP-1 Kit
Kit Catalog Numbers: K151NN-Series, K156NN-Series
Kit Lot Number: K0081197
Expiration Date: 30 Jun 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Human MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Human MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Human MCP-1	Calibrator	537	pg/mL
		Chemokine Panel 1 (human) Control 1	468	pg/mL
		Chemokine Panel 1 (human) Control 2	71.4	pg/mL
		Chemokine Panel 1 (human) Control 3	12.1	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

Human MCP-1 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
537	247,000	2.6
134	60,100	1.8
33.6	14,800	2.4
8.39	3,770	3.4
2.10	1,030	2.3
0.524	302	1.8
0.131	128	2.7
0.00	70	5.9

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MCP-1 Plate	Z0047035	2-8°C	31 Jul 2021
Anti-hu MCP-1 Antibody	D0081124	2-8°C	31 Jan 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080183	2-8°C	31 Oct 2023
Diluent 43	M0430030	≤ -10°C	30 Jun 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 2021
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
Chemokine Panel 1 (human) Control 1	Not Kit Specific
Chemokine Panel 1 (human) Control 2	Not Kit Specific
Chemokine Panel 1 (human) 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		30 APR 2019