

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-Plex Plus Human MCP-1 Kit

Kit Lot Number: K0082363

Expiration Date: 31 Jul 2025

Kit Catalog Numbers: K151NN-Series; K156NN-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MCP-1 Plate	Z0047908	2-8°C	28 Feb 2026
Anti-hu MCP-1 Antibody	D0081811	2-8°C	31 Jul 2025
Chemokine Panel 1 (human) Calibrator Blend	A0080364	2-8°C	31 May 2028
Diluent 3	M0030123	≤ -10°C	31 Jan 2026
Diluent 43	M0430053	≤ -10°C	31 Jan 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision and Accuracy 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	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	MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MCP-1	553	pg/mL

Dilution to Top of Curve = 1X

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 Nov 2023