

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

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

**Kit Lot Number:** K0082398

**Expiration Date:** 28 Feb 2026

**Kit Catalog Numbers:** K151NN-Series; K156NN-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human MCP-1 Plate	Z0047932	2-8°C	30 Apr 2026
Anti-hu MCP-1 Antibody	D0081911	2-8°C	28 Feb 2026
Chemokine Panel 1 (human) Calibrator Blend	A0080347	2-8°C	31 Oct 2027
Diluent 3	M0030124	≤ -10°C	31 Mar 2026
Diluent 43	M0430055	≤ -10°C	30 Apr 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	545	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		05 Jan 2024