

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Human MCP-1 Kit
Lot Number: K0082735
Expiration Date: 31 MAY 2026
Catalog Number: K151NN-Series; K156NN-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu MCP-1 Antibody	D0081963	2-8°C	31 MAY 2026
Chemokine Panel 1 (human) Calibrator Blend	A0080418	2-8°C	28 FEB 2030
Human MCP-1, 96-Well Small Spot SECTOR Plate	Z0048141	2-8°C	30 NOV 2027
Diluent 3	M0030141	≤ -10°C	30 NOV 2027
Diluent 43	M0430058	≤ -10°C	28 FEB 2027

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

Precision, Accuracy, and Sensitivity 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	546	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		25 AUG 2025