



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX and V-PLEX Plus Human MCP-1 Kit**Lot Number:** K0082816**Expiration Date:** 31 AUG 2027**Catalog Number:** K151NN-Series, K156NN-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu MCP-1 Antibody	D0082215	2-8°C	31 AUG 2027
Chemokine Panel 1 (human) Calibrator Blend	A0080418	2-8°C	28 FEB 2030
Human MCP-1, 96-Well Small Spot SECTOR Plate	Z0048186	2-8°C	31 MAR 2028
Diluent 3	M0030141	≤ -10°C	30 NOV 2027
Diluent 43	M0430059	≤ -10°C	30 SEP 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	Karen Hamilla	Quality		09 DEC 2025