



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human MCP-1 Gen. B Kit**Lot Number:** K0082681**Expiration Date:** 31 MAY 2027**Catalog Number:** K151ANPD-Series, K151ANPG-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080416	2-8°C	31 DEC 2029
Human MCP-1 Antibody	D0082401	2-8°C	30 JUN 2028
Human MCP-1 Gen. B 96-Well Small Spot SECTOR Plate	Z0048102	2-8°C	31 AUG 2027
Diluent 3	M0030137	≤ -10°C	31 JUL 2027
Diluent 57	M0570018	≤ -10°C	31 MAY 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) Gen. B Control 1		Chemokine Panel 1 (human) Gen. B Control 2		Chemokine Panel 1 (human) Gen. B 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	1.45	80% - 120%	< 20%	247	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MCP-1	379	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		17 APR 2025