



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human MCP-1 Gen. B Kit
Catalog Number(s): K151ANP-Series; K156ANP-Series
Lot Number: K0082353
New Expiration Date: 31 Jan 2026

Based on additional testing, the original date of expiration for this kit has been extended by 6 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 57	M0570011	31 Jul 2025	31 Jan 2026

Function	Name	Signature	Date
Quality	Michelle Likiam		19 May 2025



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human MCP-1 Gen. B Kit
Lot Number: K0082353
Expiration Date: 31 JUL 2025
Catalog Number: K151ANP-Series; K156ANP-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080377	2-8°C	31 AUG 2028
Human MCP-1 Antibody	D0082133	2-8°C	31 JAN 2027
Human MCP-1 Gen. B SECTOR Plate	Z0047904	2-8°C	28 FEB 2026
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
Diluent 57	M0570011	≤ -10°C	31 JUL 2025

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	395	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		30 MAY 2024