



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse MCP-1 Kit
Catalog Number(s): K152NN-Series
Lot Number: K0082568
New Expiration Date: 31 Dec 2025

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Mouse MCP-1, 96-Well Small Spot SECTOR Plate	Z0047727	31 Dec 2024	31 Dec 2025
Diluent 41	M0410053	30 Nov 2025	31 Dec 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		30 Oct 2024



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Mouse MCP-1 Kit
Lot Number: K0082568
Expiration Date: 31 DEC 2024
Catalog Number: K152NN-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Kit Calibrator	A0080353	2-8°C	29 FEB 2028
Mouse MCP-1, 96-Well Small Spot SECTOR Plate	Z0047727	2-8°C	31 DEC 2024
Mouse/Rat MCP-1 Antibody	D0082024	2-8°C	31 AUG 2026
Diluent 41	M0410053	≤ -10°C	30 NOV 2025
Diluent 45	M0450053	≤ -10°C	31 MAR 2026

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	MCP-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	MCP-1	4.42	75% - 125%	< 20%	325	75% - 125%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MCP-1	479	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 OCT 2024