



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Mouse MCP-1 Kit**Lot Number:** K0082883**Expiration Date:** 31 DEC 2027**Catalog Number:** K152NN-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms MCP-1 Antibody	D0082454	2-8°C	31 AUG 2028
Cytokine Panel 1 (mouse) Calibrator Blend	A0080384	2-8°C	28 FEB 2029
Mouse MCP-1, 96-Well Small Spot SECTOR Plate	Z0048222	2-8°C	31 JUL 2028
Diluent 41	M0410059	≤ -10°C	31 DEC 2027
Diluent 45	M0450056	≤ -10°C	29 FEB 2028

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%

*See Additional Comments

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MCP-1	528	pg/mL

Dilution to top of Curve: 1X

Additional Comments:

Spot	Name	Specification	Comment
1	MCP-1	LLOQ Conc.	% LLOQ measured 129% (specification 75-125%) at target 4.42 pg/mL, was fully conforming at 5.17 pg/mL.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 JUL 2026