

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse MCP-1 Kit

Kit Lot Number: K0082076

Expiration Date: 31 Jan 2024

Kit Catalog Numbers: K152NN-Series

Precision and Accuracy 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	496	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse MCP-1 Plate	Z0047727	2-8°C	31 Dec 2024
Anti-ms MCP-1 Antibody	D0081583	2-8°C	30 Jun 2024
Cytokine Panel 1 (mouse) Calibrator Blend	A0080268	2-8°C	31 Aug 2025
Diluent 41	M0410044	≤ -10°C	30 Apr 2024
Diluent 45	M0450043	≤ -10°C	31 Jan 2024

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		30 AUG 2022