

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Mouse MCP-1 Kit  
**Kit Lot Number:** K0082045  
**Expiration Date:** 30 Jun 2023  
**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	Z0047443	2-8°C	30 Jun 2023
Anti-ms MCP-1 Antibody	D0081498	2-8°C	30 Nov 2023
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		21 JUN 2022