

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse MCP-1 Kit
Kit Lot Number: K0081560
Expiration Date: 31 Mar 2022
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	491	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse MCP-1 Plate	Z0047400	2-8°C	31 Mar 2023
Anti-ms MCP-1 Antibody	D0081498	2-8°C	30 Nov 2023
Cytokine Panel 1 (mouse) Calibrator Blend	A0080125	2-8°C	31 Mar 2022
Diluent 41	M0410036	≤ -10°C	31 Mar 2023
Diluent 45	M0450037	≤ -10°C	31 Mar 2023

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		07 DEC 2020