

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 α Kit
Kit Lot Number: K0081950
Expiration Date: 30 Sep 2023
Kit Catalog Numbers: K151NQ-Series; K156NQ-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	MIP-1 α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	MIP-1 α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	MIP-1 α	1110	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 α Plate	Z0047653	2-8°C	31 Jul 2024
Anti-hu MIP-1 α Antibody	D0081508	2-8°C	30 Sep 2023
Chemokine Panel 1 (human) Calibrator Blend	A0080306	2-8°C	30 Sep 2026
Diluent 43	M0430045	\leq -10°C	30 Nov 2023
Diluent 3	M0030109	\leq -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		06 MAY 2022