

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 α Kit

Kit Lot Number: K0081780

Expiration Date: 30 Sep 2023

Kit Catalog Numbers: K151NQ-Series, K156NQ-Series

Precision and Accuracy Specifications:

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

		LLOQ			ULOQ		
Spot	Name	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	1090	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 α Plate	Z0047559	2-8°C	31 Dec 2023
Anti-hu MIP-1 α Antibody	D0081508	2-8°C	30 Sep 2023
Chemokine Panel 1 (human) Calibrator Blend	A0080279	2-8°C	28 Feb 2026
Diluent 43	M0430045	$\leq$ -10°C	30 Nov 2023
Diluent 3	M0030108	$\leq$ -10°C	30 Nov 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		09 SEP 2021