

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human MIP-1 $\alpha$  Kit  
**Kit Lot Number:** K0082269  
**Expiration Date:** 31 Mar 2025  
**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 $\alpha$	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 $\alpha$	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	MIP-1 $\alpha$	1160	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 $\alpha$ Plate	Z0047847	2-8°C	30 Sep 2025
Anti-hu MIP-1 $\alpha$ Antibody	D0081835	2-8°C	30 Sep 2025
Chemokine Panel 1 (human) Calibrator Blend	A0080347	2-8°C	31 Oct 2027
Diluent 3	M0030120	$\leq$ -10°C	31 Jul 2025
Diluent 43	M0430050	$\leq$ -10°C	31 Mar 2025

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		12 MAY 2023