


**Extension Memo**

**Product Description:** V-PLEX® and V-PLEX PLUS Human MIP-1 $\alpha$  Kit  
**Lot Number:** K0082424  
**New Expiration Date:** 31 JUL 2026  
**Catalog Number(s):** K151NQ-Series, K156NQ-Series

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Anti-hu MIP-1 $\alpha$ Antibody	D0081835	30 SEP 2025	31 JUL 2026
Diluent 3	M0030120	31 JUL 2025	31 JUL 2026
Diluent 43	M0430053	31 JAN 2026	31 JUL 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		16 Jan 2026

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX PLUS Human MIP-1 $\alpha$  Kit  
**Kit Lot Number:** K0082424  
**Expiration Date:** 31 Jul 2025  
**Kit Catalog Numbers:** K151NQ-Series; K156NQ-Series

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 $\alpha$ Plate	Z0047951	2-8°C	31 Jul 2026
Anti-hu MIP-1 $\alpha$ Antibody	D0081835	2-8°C	30 Sep 2025
Chemokine Panel 1 (human) Calibrator Blend	A0080364	2-8°C	31 May 2028
Diluent 3	M0030120	$\leq$ -10°C	31 Jul 2025
Diluent 43	M0430053	$\leq$ -10°C	31 Jan 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

### 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 $\alpha$	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 $\alpha$	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%

### Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MIP-1 $\alpha$	1180	pg/mL

Dilution to Top of Curve = 1X

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 Mar 2024