



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Human MIP-1 α Gen. B Kit
Catalog Number(s): K151ANM-Series, K156ANM-Series
Lot Number: K0082354
New Expiration Date: 31 Jul 2026

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
Human MIP-1 α Gen. B SECTOR Plate	Z0047905	28 Feb 2026	31 Jul 2026
Diluent 3	M0030123	31 Jan 2026	31 Jul 2026
Diluent 57	M0570011	31 Jul 2025	31 Jul 2026

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 Jun 2025

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 α Gen. B Kit
Kit Lot Number: K0082354
Expiration Date: 31 Jul 2025
Kit Catalog Numbers: K151ANM-Series; K156ANM-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 α Gen. B SECTOR Plate	Z0047905	2-8°C	28 Feb 2026
Human MIP-1 α Antibody	D0082134	2-8°C	31 Jan 2027
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080377	2-8°C	31 Aug 2028
Diluent 3	M0030123	$\leq$ -10°C	31 Jan 2026
Diluent 57	M0570011	$\leq$ -10°C	31 Jul 2025

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

Precision and Accuracy Specifications:

		Calibrator		Chemokine Panel 1 (human) Gen. B Control 1		Chemokine Panel 1 (human) Gen. B Control 2		Chemokine Panel 1 (human) Gen. B 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	0.357	80% - 120%	< 20%	403	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MIP-1 α	614	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		17 Oct 2023