

**Extension Memo**

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 β Gen. B Kit
Lot Number: K0082349
New Expiration Date: 31 JUL 2026
Catalog Number(s): K151ANH-Series, K156ANH-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
Human MIP-1 β Gen. B 96-Well Small Spot Plate	Z0047901	28 FEB 2026	31 JUL 2026
Diluent 3	M0030123	31 JAN 2026	31 JAN 2027
Diluent 57	M0570011	31 JUL 2025	31 JUL 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		13 Oct 2025



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 β Gen. B Kit
Lot Number: K0082349
Expiration Date: 31 JUL 2025
Catalog Number: K151ANH-Series, K156ANH-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080377	2-8°C	31 AUG 2028
Human MIP-1 β Antibody	D0082130	2-8°C	28 FEB 2027
Human MIP-1 β Gen. B SECTOR Plate	Z0047901	2-8°C	28 FEB 2026
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, Accuracy, and Sensitivity Specifications:

Spot	Name	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	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	MIP-1 β	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 β	1.20	80% - 120%	< 20%	390	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MIP-1 β	606	pg/mL

Dilution to top of Curve: 1X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		30 MAY 2024