



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX PLUS Human MIP-1 β Kit


Catalog Number(s) / Size(s): K151NR-Series; K156NR-Series

Lot Number: K0082513

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082513 has been extended to 31 OCT 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human MIP-1 β Plate	Z0047697	31 OCT 2024	31 OCT 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		30 JUL 2024

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Human MIP-1β Kit
Lot Number: K0082513
Expiration Date: 31 OCT 2024
Catalog Number: K151NR-Series; K156NR-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu MIP-1β Antibody	D0081967	2-8°C	30 JUN 2026
Chemokine Panel 1 (human) Calibrator Blend	A0080347	2-8°C	31 OCT 2027
Human MIP-1β Plate	Z0047697	2-8°C	31 OCT 2024
Diluent 3	M0030124	≤ -10°C	31 MAR 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026

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) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	MIP-1β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	MIP-1β	1.88	80% - 120%	< 20%	520	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MIP-1β	1160	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		29 JUL 2024