

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Human MIP-1 β Gen. B Kit**Lot Number:** K0082606**Expiration Date:** 31 DEC 2026**Catalog Number:** K151ANH-Series; K156ANH-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080389	2-8°C	30 APR 2029
Human MIP-1 β Antibody	D0082110	2-8°C	28 FEB 2027
Human MIP-1 β Gen. B 96-Well Small Spot SECTOR Plate	Z0048051	2-8°C	30 JUN 2027
Diluent 3	M0030131	\leq -10°C	31 JAN 2027
Diluent 57	M0570016	\leq -10°C	31 DEC 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) 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	1.2	80% - 120%	< 20%	390	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MIP-1 β	612	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		31 JAN 2025