

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Human MIP-1β Kit, SECTOR**Lot Number:** K0082849**Expiration Date:** 31 JUL 2027**Catalog Number:** K151NR-Series, K156NR-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu MIP-1β Antibody	D0082211	2-8°C	31 JUL 2027
Chemokine Panel 1 (human) Calibrator Blend	A0080432	2-8°C	30 APR 2030
Human MIP-1β 96-Well Small Spot SECTOR Plate	Z0048206	2-8°C	31 MAY 2028
Diluent 3	M0030141	≤ -10°C	30 NOV 2027
Diluent 43	M0430059	≤ -10°C	30 SEP 2027

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	
		Recovery	CV	Recovery	CV	Recovery	CV
1	MIP-1β	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%

*See Additional Comments

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	MIP-1β	890	pg/mL

Dilution to top of Curve: 1X

Additional Comments:

Spot	Name	Specification	Comment
1	MIP-1β	LLOQ Conc.	% LLOQ measured 156% (specification 80-120%) at target 1.88pg/mL, was fully conforming at 7.52pg/mL.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		06 MAR 2026