

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX PLUS Human MIP-1 $\beta$  Kit  
**Kit Lot Number:** K0082300  
**Expiration Date:** 31 Oct 2024  
**Kit Catalog Numbers:** K151NR-Series; K156NR-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Recovery	CV
		Recovery	CV	Recovery	CV	Recovery	CV		
1	MIP-1 $\beta$	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 $\beta$	1.88	80% - 120%	< 20%	520	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	MIP-1 $\beta$	1160	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 $\beta$ Plate	Z0047866	2-8°C	31 Oct 2025
Anti-hu MIP-1 $\beta$ Antibody	D0081967	2-8°C	30 Jun 2026
Chemokine Panel 1 (human) Calibrator Blend	A0080347	2-8°C	31 Oct 2027
Diluent 3	M0030121	$\leq$ -10°C	31 Oct 2025
Diluent 43	M0430049	$\leq$ -10°C	31 Oct 2024

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 JUL 2023