

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX PLUS Human IP-10 Gen. B Kit  
**Kit Lot Number:** K0082351  
**Expiration Date:** 31 Jul 2025  
**Kit Catalog Numbers:** K151ANL-Series; K156ANL-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Gen. B Plate	Z0047902	2-8°C	28 Feb 2026
Human IP-10 Antibody	D0082131	2-8°C	31 Jan 2027
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080377	2-8°C	31 Aug 2028
Diluent 3	M0030123	≤ -10°C	31 Jan 2026
Diluent 57	M0570011	≤ -10°C	31 Jul 2025

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

**Precision and Accuracy Specifications:**

		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	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IP-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%


		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IP-10	1.29	80% - 120%	< 20%	1240	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IP-10	1920	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 Oct 2023