



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Human IP-10 Gen. B Kit  
**Lot Number:** K0082475  
**Expiration Date:** 31 AUG 2026  
**Catalog Number:** K156ANL-Series, K151ANL-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080377	2-8°C	31 AUG 2028
Human IP-10 Antibody, 1 Plate	D0082219	2-8°C	31 AUG 2027
Human IP-10 Gen. B SECTOR Plate	Z0047984	2-8°C	31 OCT 2026
Diluent 3	M0030128	≤ -10°C	31 OCT 2026
Diluent 57	M0570015	≤ -10°C	31 AUG 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	IP-10	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 JUL 2024