



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human IP-10 Gen. B Kit
Lot Number: K0082572
Expiration Date: 31 AUG 2026
Catalog Number: K151ANL-Series , K156ANL-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	D0082219	2-8°C	31 AUG 2027
Human IP-10 Gen. B 96-Well Small Spot SECTOR Plate	Z0048033	2-8°C	31 MAR 2027
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		26 NOV 2024