



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IP-10 Gen B. Kit
Lot Number: K0082677
Expiration Date: 30 JUN 2027
Catalog Number: K151ANL-Series, K156ANL-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (hu) Gen. B Calibrator	A0080431	2-8°C	31 MAR 2030
Human IP-10 Antibody	D0082320	2-8°C	31 JAN 2028
Human IP-10 Gen. B 96-Well Small Spot SECTOR Plate	Z0048052	2-8°C	30 JUN 2027
Diluent 3	M0030139	≤ -10°C	29 SEP 2027
Diluent 57	M0570020	≤ -10°C	31 OCT 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) 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	1980	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		11 JUN 2025