

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IP-10 Kit

Kit Lot Number: K0081689

Expiration Date: 31 Oct 2022

Kit Catalog Numbers: K151NV-Series, K156NV-Series

Precision and Accuracy Specifications:

		Calibrator		Chemokine Panel 1 (human) Control 1		0		0	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IP-10	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IP-10	2300	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0047346	2-8°C	30 Nov 2022
Anti-hu IP-10 Antibody	D0081401	2-8°C	30 Apr 2023
Chemokine Panel 1 (human) Calibrator Blend	A0080274	2-8°C	31 Jan 2026
Diluent 43	M0430040	≤ -10°C	30 Nov 2022
Diluent 3	M0030100	≤ -10°C	31 Oct 2022

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		25 MAY 2021