

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Human IP-10 Kit
Kit Lot Number: K0082259
Expiration Date: 31 May 2024
Kit Catalog Numbers: K151NV-Series; K156NV-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) 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	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IP-10	2510	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0047678	2-8°C	31 Aug 2024
Anti-hu IP-10 Antibody	D0081608	2-8°C	31 May 2024
Chemokine Panel 1 (human) Calibrator Blend	A0080320	2-8°C	31 Mar 2027
Diluent 3	M0030116	≤ -10°C	30 Nov 2024
Diluent 43	M0430049	≤ -10°C	31 Oct 2024

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		04 MAY 2023