

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IP-10 Kit
Kit Catalog Numbers: K151NV-Series, K156NV-Series
Kit Lot Number: K0081501
Expiration Date: 31 May 2022

Acceptance Data/ Percent Recovery 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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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 Material Tested :

Spot	Name	Item	Conc	Units
1	IP-10	Calibrator	2360	pg/mL
		Chemokine Panel 1 (human) Control 1	589	pg/mL
		Chemokine Panel 1 (human) Control 2	86.8	pg/mL
		Chemokine Panel 1 (human) Control 3	14.3	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IP-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2360	474,000	2.4
590	142,000	3.5
148	36,700	3.9
36.9	9,250	2.9
9.22	2,360	4.0
2.30	638	4.1
0.576	215	2.3
0.00	71	5.4

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0047346	2-8°C	30 Nov 2022
Anti-hu IP-10 Antibody	D0081227	2-8°C	31 May 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080228	2-8°C	31 Aug 2024
Diluent 43	M0430040	≤ -10°C	30 Nov 2022
Diluent 3	M0030100	≤ -10°C	31 Oct 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Chemokine Panel 1 (human) Control 1	Not Kit Specific
Chemokine Panel 1 (human) Control 2	Not Kit Specific
Chemokine Panel 1 (human) Control 3	Not Kit Specific

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		17 AUG 2020