

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: K0081251
Expiration Date: 30 Sep 2021

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	2480	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
2480	344,000	2.5
620	133,000	2.4
155	35,200	2.9
38.8	9,020	2.7
9.69	2,380	3.2
2.42	635	4.9
0.605	207	3.2
0.00	67	11.0

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0047099	2-8°C	31 Oct 2021
Anti-hu IP-10 Antibody	D0081164	2-8°C	31 May 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080183	2-8°C	31 Oct 2023
Diluent 43	M0430032	≤ -10°C	30 Sep 2021
Diluent 3	M0030089	≤ -10°C	30 Sep 2021
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		02 JUL 2019