

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IP-10 Kit
Kit Catalog Numbers: K151NV-Series
Kit Lot Number: K0080825
Expiration Date: 31 Jul 2019

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	hu 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	hu IP-10	1.37	80% - 120%	< 20%	500	80% - 120%*	< 20%

* See Additional Comments.

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IP-10	Calibrator	2660	pg/mL
		Chemokine Panel 1 (human) Control 1	473	pg/mL
		Chemokine Panel 1 (human) Control 2	74.9	pg/mL
		Chemokine Panel 1 (human) Control 3	11.7	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IP-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2660	95,700	6.2
665	82,100	5.0
166	29,800	2.5
41.6	8,100	2.9
10.4	2,130	1.9
2.60	564	6.9
0.649	189	3.2
0.00	64	3.7

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0046531	2-8°C	30 Nov 2019
Anti-hu IP-10 Antibody	D0080901	2-8°C	31 Jul 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080122	2-8°C	28 Feb 2022
Diluent 43	M0430021	≤ -10°C	30 Sep 2019
Diluent 3	M0030076	≤ -10°C	30 Sep 2019
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

Additional Comments:

	Assay	Spot	Specification	Comment
1	hu IP-10	1	ULOQ Conc.	ULOQ recovery 124.8%

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		08 AUG 2017