

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: K0080999

Expiration Date: 30 Jun 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%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IP-10	Calibrator	2420	pg/mL
		Chemokine Panel 1 (human) Control 1	606	pg/mL
		Chemokine Panel 1 (human) Control 2	89.2	pg/mL
		Chemokine Panel 1 (human) Control 3	15.8	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
2420	323,000	5.0
605	118,000	6.1
151	31,300	5.0
37.8	8,260	5.8
9.45	2,210	3.3
2.36	615	4.4
0.591	208	7.6
0.00	76	16.0

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0046809	2-8°C	30 Sep 2020
Anti-hu IP-10 Antibody	D0080863	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080131	2-8°C	31 Jul 2022
Diluent 43	M0430024	≤ -10°C	31 Jan 2020
Diluent 3	M0030081	≤ -10°C	30 Sep 2020
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	Jessica Drolet	Quality Assurance		14 May 2019