

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP IP-10 Kit
Kit Catalog Numbers: K156NV-Series
Kit Lot Number: K0080925
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	NHP 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	NHP IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IP-10	Calibrator	2380	pg/mL
		Chemokine Panel 1 (human) Control 1	590	pg/mL
		Chemokine Panel 1 (human) Control 2	90	pg/mL
		Chemokine Panel 1 (human) Control 3	16	pg/mL

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

Standard Curve Data:

NHP IP-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2380	371,000	3.6
595	133,000	3.2
149	34,600	2.2
37.2	9,280	2.8
9.30	2,330	1.8
2.32	641	1.9
0.581	203	2.2
0.00	74	6.6

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0046698	2-8°C	31 May 2020
Anti-hu IP-10 Antibody	D0080863	2-8°C	30 Jun 2019
Chemokine Panel 1 (Human) Calibrator Blend	A0080159	2-8°C	31 Dec 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

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		14 FEB 2018