

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-Plex® and V-Plex Plus Human IL-12p70 Kit
Kit Catalog Numbers: K151QV-Series
Kit Lot Number: K0080877
Expiration Date: 31 Jul 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1(Human) Control 1		Proinflammatory Panel 1(Human) Control 2		Proinflammatory Panel 1(Human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-12p70	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-12p70	1.22	80% - 120%	< 20%	315	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-12p70	Calibrator	505	pg/mL
		Proinflammatory Panel 1(Human) Control 1	168	pg/mL
		Proinflammatory Panel 1(Human) Control 2	33.2	pg/mL
		Proinflammatory Panel 1(Human) Control 3	6.12	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-12p70 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
505	78,300	3.9
126	22,400	3.5
31.6	5,420	3.3
7.89	1,450	4.1
1.97	406	4.0
0.493	155	4.1
0.123	82	4.6
0.00	62	7.9

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12p70 Plate	Z0046617	2-8°C	31 Jan 2020
Anti-hu IL-12p70 Antibody	D0080921	2-8°C	31 Jul 2019
Proinflammatory Panel 1 (human) Calibrator Blend	A0080139	2-8°C	31 Jul 2022
Diluent 2	M0020096	≤ -10°C	30 Nov 2019
Diluent 3	M0030078	≤ -10°C	31 Jan 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), 100 mL	Not Kit Specific
Proinflammatory Panel 1(human) Control 1	Not Kit Specific
Proinflammatory Panel 1(human) Control 2	Not Kit Specific
Proinflammatory 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		02 OCT 2017