

**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:** K0080987

**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	380	pg/mL
		Proinflammatory Panel 1 (human) Control 1	182	pg/mL
		Proinflammatory Panel 1 (human) Control 2	34	pg/mL
		Proinflammatory Panel 1 (human) Control 3	6.95	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
380	88,100	3.6
95.0	23,800	5.9
23.8	6,330	5.1
5.94	1,620	3.8
1.48	449	2.9
0.371	157	3.6
0.0928	87	4.3
0.00	63	8.3

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12p70 Plate	Z0046797	2-8°C	30 Sep 2020
Anti-hu IL-12p70 Antibody	D0080921	2-8°C	31 Jul 2019
Proinflammatory Panel 1 (human) Calibrator Blend	A0080050	2-8°C	31 Jan 2020
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
<b>Review/Approval</b>	Jessica Drolet	Quality Assurance		17 May 2018