

**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:** K0081115  
**Expiration Date:** 31 Jan 2020

**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	496	pg/mL
		Proinflammatory Panel 1 (human) Control 1	186	pg/mL
		Proinflammatory Panel 1 (human) Control 2	35.6	pg/mL
		Proinflammatory Panel 1 (human) Control 3	6.99	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
496	101,000	3.0
124	27,900	3.7
31.0	7,150	5.1
7.75	1,800	3.6
1.94	480	2.2
0.484	164	5.5
0.121	89	3.6
0.00	61	9.4

## 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	D0081019	2-8°C	31 Jan 2020
Proinflammatory Panel 1 (human) Calibrator Blend	A0080102	2-8°C	30 Sep 2021
Diluent 2	M0020098	≤ -10°C	28 Feb 2020
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), 100mL	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 NOV 2018