

**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:** K0081245  
**Expiration Date:** 31 Mar 2021

**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	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	IL-12p70	1.22	80% - 120%	< 20%	315	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	IL-12p70	Calibrator	533	pg/mL
		Proinflammatory Panel 1 (human) Control 1	195	pg/mL
		Proinflammatory Panel 1 (human) Control 2	38.3	pg/mL
		Proinflammatory Panel 1 (human) Control 3	7.56	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

IL-12p70 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
533	119,000	2.7
133	33,700	3.8
33.3	8,780	3.5
8.33	2,290	2.3
2.08	604	3.9
0.521	186	3.5
0.130	88	2.9
0.00	56	8.7

## CERTIFICATE OF ANALYSIS

Page 2 of 2

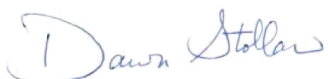
### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12 p70 Plate	Z0047090	2-8°C	31 Oct 2021
Anti-hu IL-12p70 Antibody	D0081161	2-8°C	31 Mar 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080158	2-8°C	31 May 2023
Diluent 2	M0020103	≤ -10°C	31 Mar 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 2021
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>	Dawn Stollar	Quality		11 JUN 2019