

**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:** K0081494  
**Expiration Date:** 31 Oct 2022

**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	484	pg/mL
		Proinflammatory Panel 1 (human) Control 1	158	pg/mL
		Proinflammatory Panel 1 (human) Control 2	31.4	pg/mL
		Proinflammatory Panel 1 (human) Control 3	5.92	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
484	81,800	3.0
121	21,800	4.5
30.3	5,820	3.1
7.56	1,580	3.1
1.89	430	3.3
0.473	156	3.3
0.118	86	2.8
0.00	65	7.4

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12p70 Plate	Z0047337	2-8°C	30 Nov 2022
Anti-hu IL-12p70 Antibody	D0081307	2-8°C	31 Oct 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080254	2-8°C	30 Apr 2025
Diluent 2	M0020113	≤ -10°C	30 Nov 2022
Diluent 3	M0030100	≤ -10°C	31 Oct 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	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>	Karen Hamilla	Quality		22 JUL 2020