

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: K0081390
Expiration Date: 31 Oct 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	159	pg/mL
		Proinflammatory Panel 1 (human) Control 2	32.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	6.4	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	101,000	3.7
133	28,500	3.0
33.3	7,660	2.7
8.33	1,950	3.1
2.08	541	3.2
0.521	179	4.4
0.130	85	3.1
0.00	57	5.3

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12p70 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	M0020109	≤ -10°C	28 Feb 2022
Diluent 3	M0030094	≤ -10°C	31 May 2022
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	Karen Hamilla	Quality		24 JAN 2020