

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: K0081458
Expiration Date: 31 Jan 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	473	pg/mL
		Proinflammatory Panel 1 (human) Control 1	197	pg/mL
		Proinflammatory Panel 1 (human) Control 2	38.8	pg/mL
		Proinflammatory Panel 1 (human) Control 3	7.11	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
473	108,000	4.7
118	29,000	3.8
29.6	7,710	5.6
7.39	1,980	2.4
1.85	539	4.4
0.462	186	3.2
0.115	100	5.7
0.00	71	8.7

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12p70 Plate	Z0047297	2-8°C	31 Aug 2022
Anti-hu IL-12p70 Antibody	D0081127	2-8°C	31 Jan 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080229	2-8°C	30 Sep 2024
Diluent 2	M0020112	≤ -10°C	31 Jul 2022
Diluent 3	M0030097	≤ -10°C	31 Aug 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), 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	Karen Hamilla	Quality		08 MAY 2020