

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: K0081196
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	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	550	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:

hu IL-12p70 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
550	65,800	1.9
138	18,600	3.5
34.4	4,910	1.6
8.59	1,240	2.5
2.15	343	1.7
0.537	122	3.8
0.134	70	7.4
0.00	50	9.6

CERTIFICATE OF ANALYSIS

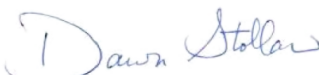
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12p70 Plate	Z0047036	2-8°C	31 Jul 2021
Anti-hu IL-12p70 Antibody	D0081161	2-8°C	31 Mar 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080172	2-8°C	30 Jun 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), 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		17 APR 2019