

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-12p70 Kit  
**Kit Catalog Numbers:** K152QV-Series  
**Kit Lot Number:** K0081413  
**Expiration Date:** 31 Dec 2021

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-12p70	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%*	< 25%	75% - 125%	< 25%

\* See Additional Comments.

Spot	Name	Value	LLOQ			ULOQ		
			Recovery	CV	Value	Recovery	CV	
1	IL-12p70	179	75% - 125%	< 25%	20600	75% - 125%	< 25%	

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-12p70	Calibrator	31900	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	13500	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	3380	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	806	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
31900	410,000	1.8
7980	83,300	2.7
1990	15,900	3.7
498	3,790	2.5
125	1,010	4.2
31.2	312	4.1
7.79	131	4.7
0.00	75	7.4

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-12p70 Plate	Z0047257	2-8°C	30 Jun 2022
Anti-ms IL-12p70 Antibody	D0081154	2-8°C	30 Apr 2022
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080197	2-8°C	31 Dec 2023
Diluent 41	M0410029	≤ -10°C	31 Dec 2021
Diluent 45	M0450030	≤ -10°C	28 Feb 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 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 3	Not Kit Specific

### Additional Comments:

	Assay	Spot	Specification	Comment
1	IL-12p70	1	Control Conc.	Control 2 % Recovery = 132.9%

### 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		18 FEB 2020