

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus mouse IL-6 Kit  
**Kit Catalog Numbers:** K152QX-Series  
**Kit Lot Number:** K0081170  
**Expiration Date:** 31 May 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	ms IL-6	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-6	7.61	75% - 125%	< 25%	3140	75% - 125%	< 25%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-6	Calibrator	4990	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	1050	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	167	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	31.5	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

ms IL-6 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
4990	126,000	3.2
1250	28,100	2.6
312	6,780	2.5
78.0	1,670	3.3
19.5	498	3.2
4.87	196	11.0
1.22	124	8.4
0.00	103	15.0

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-6 Plate	Z0047005	2-8°C	30 Jun 2021
Anti-ms IL-6 Antibody	D0081088	2-8°C	31 Oct 2021
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080152	2-8°C	30 Nov 2022
Diluent 41	M0410026	≤ -10°C	30 Jun 2021
Diluent 45	M0450025	≤ -10°C	31 May 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 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) 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>	Dawn Stollar	Quality Assurance		18 MAR 2019