

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® AND V-PLEX Plus Mouse TNF- $\alpha$  Kit

**Kit Catalog Numbers:** K152QW-Series

**Kit Lot Number:** K0081383

**Expiration Date:** 28 Feb 2022

### 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	TNF- $\alpha$	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF- $\alpha$	0.980	75% - 125%	< 25%	403	75% - 125%	< 25%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	TNF- $\alpha$	Calibrator	577	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	166	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	29.9	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.93	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

TNF- $\alpha$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
577	104,000	4.9
144	24,400	7.2
36.1	6,090	5.2
9.02	1,550	6.2
2.25	445	3.8
0.563	168	4.3
0.141	100	3.9
0.00	82	6.5

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- $\alpha$ Plate	Z0047234	2-8°C	31 May 2022
Anti-ms TNF- $\alpha$ Antibody	D0081226	2-8°C	31 May 2022
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080197	2-8°C	31 Dec 2023
Diluent 41	M0410030	$\leq$ -10°C	31 Mar 2022
Diluent 45	M0450030	$\leq$ -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

### Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		27 MAR 2020