

**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:** K0080962  
**Expiration Date:** 31 Mar 2019

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

\* See Additional Comments.

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

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	ms TNF- $\alpha$	Calibrator	674	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	152	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	29.1	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	5.1	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

ms TNF- $\alpha$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
674	173,000	1.7
169	39,700	2.2
42.1	9,720	2.9
10.5	2,520	3.8
2.63	703	2.6
0.658	251	1.8
0.165	128	3.4
0.00	93	4.0

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- $\alpha$ Plate	Z0046764	2-8°C	31 Jul 2020
Anti-ms TNF- $\alpha$ Antibody	D0080787	2-8°C	31 Mar 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080138	2-8°C	31 May 2022
Diluent 41	M0410020	$\leq$ -10°C	29 Feb 2020
Diluent 45	M0450020	$\leq$ -10°C	31 May 2020
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

### Additional Comments:

	Assay	Spot	Specification	Comment
1	ms TNF- $\alpha$	1	Control % Recovery	Control-1 recovery 137.4% Control-2 recovery 129.5%

### 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	<i>Dawn Stollar</i>	04 APR 2018