

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human TNF- $\alpha$  Kit  
**Kit Catalog Numbers:** K151QW-Series  
**Kit Lot Number:** K0081024  
**Expiration Date:** 31 Jan 2020

### 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 TNF- $\alpha$	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu TNF- $\alpha$	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu TNF- $\alpha$	Calibrator	369	pg/mL
		Proinflammatory Panel 1 (human) Control 1	86.7	pg/mL
		Proinflammatory Panel 1 (human) Control 2	16	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.06	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

hu TNF- $\alpha$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
369	159,000	8.1
92.3	40,100	2.9
23.1	10,200	3.0
5.77	2,560	3.1
1.44	699	4.9
0.360	231	3.0
0.0901	110	3.4
0.00	70	5.2

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- $\alpha$ Plate	Z0046845	2-8°C	30 Nov 2020
Anti-hu TNF- $\alpha$ Antibody	D0081020	2-8°C	31 Jan 2020
Proinflammatory Panel 1 (human) Calibrator Blend	A0080169	2-8°C	31 May 2023
Diluent 2	M0020099	$\leq$ -10°C	31 Mar 2020
Diluent 3	M0030082	$\leq$ -10°C	31 Oct 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 (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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		10 JUL 2018