

## 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:** K0080909  
**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	388	pg/mL
		Proinflammatory Panel 1 (human) Control 1	89.6	pg/mL
		Proinflammatory Panel 1 (human) Control 2	18	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.72	pg/mL

Controls tested may not be the same lot as shipped.

**CERTIFICATE OF ANALYSIS**

Page 2 of 2

**Standard Curve Data:**

hu TNF- $\alpha$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
388	204,000	1.7
97.0	51,500	1.5
24.3	13,100	0.8
6.06	3,290	2.0
1.52	877	1.9
0.379	271	1.7
0.0947	118	3.7
0.00	69	7.7


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- $\alpha$ Plate	Z0046680	2-8°C	30 Apr 2020
Anti-hu TNF- $\alpha$ Antibody	D0081020	2-8°C	31 Jan 2020
Proinflammatory Panel 1 (human) Calibrator Blend	A0080141	2-8°C	31 Aug 2022
Diluent 2	M0020098	$\leq$ -10°C	28 Feb 2020
Diluent 3	M0030078	$\leq$ -10°C	31 Jan 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		04 JAN 2018