

**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:** K0081388  
**Expiration Date:** 30 Sep 2021

**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	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	TNF- $\alpha$	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	TNF- $\alpha$	Calibrator	347	pg/mL
		Proinflammatory Panel 1 (human) Control 1	87.8	pg/mL
		Proinflammatory Panel 1 (human) Control 2	15.7	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.02	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
347	113,000	1.9
86.8	26,900	3.4
21.7	6,800	4.0
5.42	1,770	3.8
1.36	484	2.6
0.339	158	2.9
0.0847	81	4.3
0.00	52	7.4

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- $\alpha$ Plate	Z0047066	2-8°C	30 Sep 2021
Anti-hu TNF- $\alpha$ Antibody	D0081099	2-8°C	30 Nov 2021
Proinflammatory Panel 1 (human) Calibrator Blend	A0080187	2-8°C	31 Dec 2023
Diluent 2	M0020108	$\leq$ -10°C	28 Feb 2022
Diluent 3	M0030090	$\leq$ -10°C	31 Dec 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 (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>	Karen Hamilla	Quality		06 DEC 2019