

**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:** K0081534  
**Expiration Date:** 30 Nov 2022

**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-a	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF-a	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	TNF-a	Calibrator	332	pg/mL
		Proinflammatory Panel 1 (human) Control 1	96.2	pg/mL
		Proinflammatory Panel 1 (human) Control 2	16.9	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.32	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

TNF-a - Spot 1		
Conc. pg/mL	ECL Signal	%CV
332	143,000	1.7
83.0	34,800	2.2
20.8	8,150	2.2
5.19	2,050	2.4
1.30	564	2.5
0.324	193	2.4
0.0811	98	5.8
0.00	73	5.4

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF-a Plate	Z0047375	2-8°C	31 Jan 2023
Anti-hu TNF-a Antibody	D0081324	2-8°C	31 Dec 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080208	2-8°C	30 Apr 2024
Diluent 2	M0020113	≤ -10°C	30 Nov 2022
Diluent 3	M0030101	≤ -10°C	31 Dec 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), 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		09 OCT 2020