

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX PLUS Human TNF- $\alpha$  Kit  
**Kit Lot Number:** K0082286  
**Expiration Date:** 31 Aug 2025  
**Kit Catalog Numbers:** K151QW-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- $\alpha$ Plate	Z0047860	2-8°C	31 Oct 2025
Anti-hu TNF- $\alpha$ Antibody	D0081825	2-8°C	31 Aug 2025
Proinflammatory Panel 1 (human) Calibrator Blend	A0080363	2-8°C	31 Mar 2028
Diluent 2	M0020126	$\leq$ -10°C	31 Aug 2025
Diluent 3	M0030121	$\leq$ -10°C	31 Oct 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

**Precision and Accuracy 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 Calibrator:**

Spot	Name	Conc.	Units
1	TNF- $\alpha$	368	pg/mL

Dilution to Top of Curve = 1X

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 Sep 2023