

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Human TNF- $\alpha$  Kit**Lot Number:** K0082603**Expiration Date:** 31 MAR 2027**Catalog Number:** K151QW-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu TNF- $\alpha$ Antibody	D0082164	2-8°C	30 APR 2027
Human TNF- $\alpha$ 96-Well Small Spot SECTOR Plate	Z0048050	2-8°C	31 MAY 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080411	2-8°C	31 OCT 2029
Diluent 2	M0020132	$\leq -10^\circ\text{C}$	31 MAR 2027
Diluent 3	M0030132	$\leq -10^\circ\text{C}$	31 MAR 2027

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

**Precision, Accuracy, and Sensitivity 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- $\alpha$	369	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		24 MAR 2025