



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Human TNF- α Kit QuickPlex**Lot Number:** K0082645**Expiration Date:** 31 MAR 2027**Catalog Number:** K151QW-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu TNF- α Antibody	D0082218	2-8°C	31 AUG 2027
Human TNF- α 96-Well Small Spot QuickPlex Ultra Plate	Z0048078	2-8°C	31 JUL 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080385	2-8°C	31 DEC 2028
Diluent 2	M0020132	$\leq -10^{\circ}\text{C}$	31 MAR 2027
Diluent 3	M0030136	$\leq -10^{\circ}\text{C}$	31 MAY 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- α	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%


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

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TNF- α	354	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
Review/Approval	Karen Hamilla	Quality		04 APR 2025