



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human TNF- α Kit
Lot Number: K0082709
Expiration Date: 30 SEP 2027
Catalog Number: K151QW-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu TNF- α Antibody	D0082434	2-8°C	31 AUG 2028
Human TNF- α 96-Well Small Spot SECTOR Plate	Z0048123	2-8°C	31 OCT 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080427	2-8°C	28 FEB 2030
Diluent 2	M0020134	$\leq -10^{\circ}\text{C}$	30 SEP 2027
Diluent 3	M0030140	$\leq -10^{\circ}\text{C}$	30 SEP 2027

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TNF- α	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

		LLOQ			ULOQ		
Spot	Name	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	362	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		09 JUN 2025