



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human TNF- α V2 Kit, SECTOR

Lot Number: K0082649

Expiration Date: 31 JUL 2027

Catalog Number: K151AVG-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- α Antibody	D0082464	2-8°C	30 SEP 2028
Human TNF- α V2 96-Well Small Spot SECTOR Plate	Z0048081	2-8°C	31 JUL 2027
Diluent 58	M0580015	\leq -10°C	31 JUL 2027
Cytolytic Panel 1 (human) Calibrator Blend	A0080419	\leq -70°C	31 JUL 2027

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

Precision, Accuracy, and Sensitivity Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF- α	0.898	80% - 120%	< 20%	1360	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TNF- α	40100	pg/mL

Dilution to top of Curve: 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		11 SEP 2025