

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX Human TNF- α Kit
Kit Lot Number: K0082433
Expiration Date: 31 Dec 2025
Kit Catalog Numbers: K151QW-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- α Plate	Z0047957	2-8°C	31 Aug 2026
Anti-hu TNF- α Antibody	D0081878	2-8°C	31 Dec 2025
Proinflammatory Panel 1 (human) Calibrator Blend	A0080385	2-8°C	31 Dec 2028
Diluent 2	M0020129	$\leq -10^{\circ}\text{C}$	30 Jun 2026
Diluent 3	M0030126	$\leq -10^{\circ}\text{C}$	31 Jul 2026

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- 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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 Apr 2024