



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX Human TNF- α Kit
Lot Number: K0082538
Expiration Date: 30 NOV 2026
Catalog Number: K151QW-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu TNF- α Antibody	D0082120	2-8°C	31 JAN 2027
Human TNF- α Plate	Z0048014	2-8°C	31 JAN 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080395	2-8°C	31 MAR 2029
Diluent 2	M0020131	$\leq -10^\circ\text{C}$	31 DEC 2026
Diluent 3	M0030129	$\leq -10^\circ\text{C}$	30 NOV 2026

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	0.69	80% - 120%	< 20%	248	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TNF- α	363	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		11 OCT 2024