

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Rat TNF- α Kit
Lot Number: K0082508
Expiration Date: 30 SEP 2025
Catalog Number: K153QW-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat TNF- α Antibody	D0082001	2-8°C	31 JUL 2026
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080386	2-8°C	28 FEB 2029
Rat TNF- α Plate	Z0047997	2-8°C	31 DEC 2026
Diluent 40	M0400020	$\leq -10^{\circ}\text{C}$	28 FEB 2026
Diluent 42	M0420019	$\leq -10^{\circ}\text{C}$	30 SEP 2025
Blocker H	R0100024	Room Temperature	28 FEB 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 (rat) Control 1		Proinflammatory Panel 1 (rat) Control 2		Proinflammatory Panel 1 (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TNF- α	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF- α	9.1	75% - 125%	< 25%	793	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TNF- α	1200	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		27 AUG 2024