



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Mouse TNF- α Kit
Lot Number: K0082487
Expiration Date: 30 NOV 2025
Catalog Number: K152QW-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms TNF- α Antibody	D0082022	2-8°C	31 AUG 2026
Mouse TNF- α Plate	Z0047990	2-8°C	30 NOV 2026
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080373	2-8°C	31 AUG 2028
Diluent 41	M0410053	$\leq -10^{\circ}\text{C}$	30 NOV 2025
Diluent 45	M0450053	$\leq -10^{\circ}\text{C}$	31 MAR 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 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TNF- α	80% - 120%	< 20%	100% - 150%	< 25%	100% - 150%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF- α	0.98	75% - 125%	< 25%	403	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TNF- α	610	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		15 JUL 2024