


**Extension Memo**

Product Description: V-PLEX® and V-PLEX Plus Mouse TNF- α Kit
Lot Number: K0082487
New Expiration Date: 30 NOV 2026
Catalog Number(s): K152QW-Series

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Anti-ms TNF- α Antibody	D0082022	31 AUG 2026	30 NOV 2026
Diluent 41	M0410053	30 NOV 2025	30 NOV 2026
Diluent 45	M0450053	31 MAR 2026	30 NOV 2026

Function	Name	Signature	Date
Quality	Michelle Likiam		02 Dec 2025



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:

		Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TNF- α	80% - 120%	< 20%	100% - 150%	< 25%	100% - 150%	< 25%	75% - 125%	< 25%

		LLOQ			ULOQ		
Spot	Name	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