



Extension Memo


Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Rat TNF- α Kit
Catalog Number(s): K153QW-Series
Lot Number: K0082508
New Expiration Date: 30 Sep 2026

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
Blocker H	R0100024	28 Feb 2026	30 Sep 2026
Diluent 40	M0400020	28 Feb 2026	30 Sep 2026
Diluent 42	M0420019	30 Sep 2025	30 Sep 2026
Anti-rat TNF- α Antibody	D0082001	31 Jul 2026	30 Sep 2026

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		28 Jul 2025



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:

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


		LLOQ			ULOQ		
Spot	Name	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