



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX PLUS Rat TNF- $\alpha$  Kit


**Catalog Number(s) / Size(s):** K153QW-Series

**Lot Number:** K0082345

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082345 has been extended to 30 JUN 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Rat TNF- $\alpha$ Plate	Z0047647	30 JUN 2024	30 JUN 2025
Anti-rat TNF- $\alpha$ Antibody	D0081641	31 AUG 2024	30 JUN 2025
Blocker H	R0100022	31 OCT 2024	30 JUN 2025
Diluent 40	M0400019	31 OCT 2024	30 JUN 2025
Diluent 42	M0420018	31 MAY 2025	30 JUN 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		11 MAR 2024

This memo supersedes all other statements regarding date of expiration for this product.

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX PLUS Rat TNF- $\alpha$  Kit

**Kit Lot Number:** K0082345

**Expiration Date:** 30 Jun 2024

**Kit Catalog Numbers:** K153QW-Series

**Precision and Accuracy 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- $\alpha$	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF- $\alpha$	9.10	75% - 125%	< 25%	793	75% - 125%	< 25%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	TNF- $\alpha$	1283	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat TNF- $\alpha$ Plate	Z0047647	2-8°C	30 Jun 2024
Anti-rat TNF- $\alpha$ Antibody	D0081641	2-8°C	31 Aug 2024
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080376	2-8°C	31 Jul 2028
Blocker H	R0100022	Room Temperature	31 Oct 2024
Diluent 40	M0400019	$\leq$ -10°C	31 Oct 2024
Diluent 42	M0420018	$\leq$ -10°C	31 May 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 SEP 2023