



Extension Memo

Product Description: Diluent 42
Lot Number: M0420021
New Expiration Date: 31 AUG 2027
Catalog Number(s): R50AK-Series

Based on additional testing, the original date of expiration for this diluent has been extended by 5 months.

Function	Name	Signature	Date
Quality	Michelle Likiam		02 Jun 2026



Extension Memo

Product Description: Diluent 40
Lot Number: M0400022
New Expiration Date: 31 AUG 2027
Catalog Number(s): R50AJ-Series

Based on additional testing, the original date of expiration for this diluent has been extended by 9 months.

Function	Name	Signature	Date
Quality	Michelle Likiam		02 Jun 2026



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Rat TNF- α Kit**Lot Number:** K0082968**Expiration Date:** 31 AUG 2027**Catalog Number:** K153QW-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat TNF- α Antibody	D0082393	2-8°C	30 NOV 2028
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080386	2-8°C	28 FEB 2029
Rat TNF- α 96-Well Small Spot SECTOR Plate	Z0048275	2-8°C	30 NOV 2028
Diluent 40	M0400022	$\leq -10^{\circ}\text{C}$	31 AUG 2027
Diluent 42	M0420021	$\leq -10^{\circ}\text{C}$	31 AUG 2027
Blocker H	R0100027	Room Temperature	30 APR 2028

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	9.10	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		08 SEP 2026