



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX and V-PLEX Plus Human TNF- α Kit


Catalog Number(s) / Size(s): K151QW-Series

Lot Number: K0082444

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082444 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
Human TNF- α Plate	Z0047638	30 JUN 2024	30 JUN 2025
Diluent 2	M0020123	28 FEB 2025	30 JUN 2025
Diluent 3	M0030118	31 MAR 2025	30 JUN 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		05 APR 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 Human TNF- α Kit

Kit Lot Number: K0082444

Expiration Date: 30 Jun 2024

Kit Catalog Numbers: K151QW-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- α Plate	Z0047638	2-8°C	30 Jun 2024
Anti-hu TNF- α Antibody	D0081777	2-8°C	31 Aug 2025
Proinflammatory Panel 1 (human) Calibrator Blend	A0080355	2-8°C	31 Jan 2028
Diluent 2	M0020123	$\leq -10^\circ\text{C}$	28 Feb 2025
Diluent 3	M0030118	$\leq -10^\circ\text{C}$	31 Mar 2025

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

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TNF- α	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%


Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF- α	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TNF- α	290	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 Apr 2024