



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** S-PLEX Human TNF- $\beta$  Kit  
**Catalog Number(s) / Size(s):** K151ADPS-Series; K156ADPS-Series  
**Lot Number:** K00S0217

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0217 has been extended to 30 APR 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Biotin Human TNF- $\beta$ Antibody	B00T0044A	30 APR 2024	30 APR 2025
TURBO-BOOST Human TNF- $\beta$ Antibody	D00T0044A	31 MAY 2024	30 APR 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		23 JUL 2024

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

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX Human TNF- $\beta$  Kit  
**Lot Number:** K00S0217  
**Expiration Date:** 30 APR 2024  
**Catalog Number:** K151ADPS-Series; K156ADPS-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TNF- $\beta$ Antibody	B00T0044A	2-8°C	30 APR 2024
Human TNF- $\beta$ Calibrator	A00S0032	2-8°C	31 AUG 2025
TURBO-BOOST Human TNF- $\beta$ Antibody	D00T0044A	2-8°C	31 MAY 2024
Blocker S1	R0140008	$\leq -10^{\circ}\text{C}$	31 JUL 2026
Diluent 2	M0020128	$\leq -10^{\circ}\text{C}$	31 MAR 2026
S-PLEX Enhance E1	Y0550017	$\leq -10^{\circ}\text{C}$	31 AUG 2026
S-PLEX Enhance E2	Y0560018	$\leq -10^{\circ}\text{C}$	31 AUG 2026
S-PLEX Coating Reagent C1	476750	$\leq -70^{\circ}\text{C}$	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580025	$\leq -70^{\circ}\text{C}$	30 SEP 2026
S-PLEX Detect D2	Y0590019	$\leq -70^{\circ}\text{C}$	30 SEP 2026
S-PLEX Enhance E3	Y0570020	$\leq -70^{\circ}\text{C}$	31 OCT 2026

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

**Precision, Accuracy, and Sensitivity Specifications:**

Spot Number	Name	Calibrator	
		Recovery	CV
1	TNF-beta	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
TNF-beta	1.39 - 22.2	11.0	fg/mL	80% - 120%	< 20%	2.69	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
TNF-beta	1X Calibrator	30000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		23 JUL 2024