



Extension Memo

Page 1 of 1

Product Description: S-PLEX Human TNF- β Kit
Catalog Number(s): K151ADPS-Series; K156ADPS-Series
Lot Number: K00S0230
New Expiration Date: 31 Aug 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
Human TNF- β Calibrator	A00S0032	31 Aug 2025	31 Aug 2026
S-PLEX Coating Reagent C1	417723	31 Aug 2025	31 Aug 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		26 Mar 2025



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human TNF- β Kit
Lot Number: K00S0230
Expiration Date: 31 AUG 2025
Catalog Number: K151ADPS-Series; K156ADPS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TNF- β Antibody	B00T0233	2-8°C	31 MAR 2028
Human TNF- β Calibrator	A00S0032	2-8°C	31 AUG 2025
TURBO-BOOST Human TNF- β Antibody	D00T0233	2-8°C	31 MAR 2028
Diluent 2	M0020133	$\leq -10^{\circ}\text{C}$	30 JUN 2027
S-PLEX Enhance E1	Y0550015	$\leq -10^{\circ}\text{C}$	31 MAY 2026
S-PLEX Enhance E2	Y0560016	$\leq -10^{\circ}\text{C}$	31 MAY 2026
S-PLEX Coating Reagent C1	417723	$\leq -70^{\circ}\text{C}$	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580022	$\leq -70^{\circ}\text{C}$	31 MAY 2026
S-PLEX Detect D2	Y0590016	$\leq -70^{\circ}\text{C}$	30 APR 2026
S-PLEX Enhance E3	Y0570017	$\leq -70^{\circ}\text{C}$	31 MAR 2026
Blocker S1	R0140013	-10°C	31 JUL 2027

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- β	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
TNF- β	1.39 - 22.2	11.1	fg/mL	80% - 120%	< 20%	3.53	fg/mL

Stock Concentration of Calibrator:

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 MAR 2025