

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human TNF- α Kit
Lot Number: K00S0256
Expiration Date: 30 NOV 2025
Catalog Number: K151E3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TNF- α Antibody	B00T0203	2-8°C	31 AUG 2027
Human TNF- α Calibrator	A00S0079	2-8°C	31 DEC 2027
TURBO-BOOST Human TNF- α Antibody	D00T0203	2-8°C	31 AUG 2027
Blocker S1	R0140007	$\leq -10^{\circ}\text{C}$	30 JUN 2026
Diluent 2	M0020127	$\leq -10^{\circ}\text{C}$	30 NOV 2025
Diluent 3	M0030128	$\leq -10^{\circ}\text{C}$	31 OCT 2026
S-PLEX Enhance E1	Y0550016	$\leq -10^{\circ}\text{C}$	31 JUL 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 TM Reagent	Y0580030	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Detect D2	Y0590022	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Enhance E3	Y0570023	$\leq -70^{\circ}\text{C}$	30 APR 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- α	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
TNF- α	6.25 - 100	12.5	fg/mL	80% - 120%	< 20%	3.72	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TNF- α	5X Calibrator	408000	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		05 MAR 2025