



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human TNF- α Kit

Catalog Number: K151E3S-Series

Lot Number: K00S0087

Expiration Date: 30 APR 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human TNF- α Antibody	D00T0090	2-8°C	31 DEC 2025
Biotin Human TNF- α Antibody	B00T0090	2-8°C	31 MAR 2026
Human TNF- α Calibrator	A00S0030	2-8°C	31 AUG 2025
Diluent 2	M0020123	$\leq -10^{\circ}\text{C}$	28 FEB 2025
Diluent 3	M0030117	$\leq -10^{\circ}\text{C}$	31 JAN 2025
S-PLEX Coating Reagent C1	351987	$\leq -70^{\circ}\text{C}$	30 APR 2024
Blocker S1	R0010768	$\leq -10^{\circ}\text{C}$	31 MAY 2024
S-PLEX Enhance E1	Y0550009	$\leq -10^{\circ}\text{C}$	28 FEB 2025
S-PLEX Enhance E2	Y0560011	$\leq -10^{\circ}\text{C}$	31 MAR 2025
S-PLEX Enhance E3	Y0570010	$\leq -70^{\circ}\text{C}$	31 JAN 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580014	$\leq -70^{\circ}\text{C}$	28 FEB 2025
S-PLEX Detect D2	Y0590010	$\leq -70^{\circ}\text{C}$	28 FEB 2025

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

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
TNF- α	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
TNF- α	6.25 – 100	50.0	fg/mL	80% - 120%	< 20%	4.88	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TNF- α	5X Calibrator	353000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 OCT 2022