



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human TNF- $\alpha$  Kit

**Catalog Number:** K151E3S-Series

**Lot Number:** K00S0054

**Expiration Date:** 30 APR 2023

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human TNF- $\alpha$ Antibody	D00T0063A	2-8°C	31 Jul 2024
Biotin Human TNF- $\alpha$ Antibody	B00T0063	2-8°C	31 Dec 2024
Human TNF- $\alpha$ Calibrator	A00S0030	2-8°C	31 Aug 2025
Diluent 2	M0020117	$\leq -10^{\circ}\text{C}$	31 Oct 2023
Diluent 3	M0030109	$\leq -10^{\circ}\text{C}$	31 Jan 2024
S-PLEX Coating Reagent C1	329998	$\leq -70^{\circ}\text{C}$	31 Aug 2023
Blocker S1	R0140004	$\leq -10^{\circ}\text{C}$	30 Apr 2023
S-PLEX Enhance E1	Y0550006	$\leq -10^{\circ}\text{C}$	31 Oct 2023
S-PLEX Enhance E2	Y0560007	$\leq -10^{\circ}\text{C}$	31 Oct 2023
S-PLEX Enhance E3	Y0570006	$\leq -70^{\circ}\text{C}$	31 Oct 2023
S-PLEX Detect D1	Y0580008	$\leq -70^{\circ}\text{C}$	30 Nov 2023
S-PLEX Detect D2	Y0590006	$\leq -70^{\circ}\text{C}$	31 Oct 2023

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- $\alpha$	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
TNF- $\alpha$	6.25 - 100	12.5	fg/mL	80% - 120%	< 20%	4.28	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TNF- $\alpha$	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		09 DEC 2021