



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®]Human TNF- β Kit
Catalog Number: K151ADPS-Series; K156ADPS - Series
Lot Number: K00S0151
Expiration Date: 30 Apr 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TNF- β Antibody	B00T0044A	2-8°C	30 Apr 2024
TURBO-BOOST [™] Human TNF- β Antibody	D00T0044A	2-8°C	31 May 2024
S-PLEX Coating Reagent C1	417723	$\leq -70^{\circ}\text{C}$	31 Aug 2025
Blocker S1	R0010831	$\leq -10^{\circ}\text{C}$	31 Jul 2025
S-PLEX Enhance E1	Y0550012	$\leq -10^{\circ}\text{C}$	31 Aug 2025
S-PLEX Enhance E2	Y0560014	$\leq -10^{\circ}\text{C}$	31 Oct 2025
S-PLEX Enhance E3	Y0570014	$\leq -70^{\circ}\text{C}$	31 Oct 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580018	$\leq -70^{\circ}\text{C}$	31 Oct 2025
S-PLEX Detect D2	Y0590014	$\leq -70^{\circ}\text{C}$	31 Oct 2025
Human TNF- β Calibrator	A00S0032	2-8°C	31 Aug 2025
Diluent 2	M0020124	$\leq -10^{\circ}\text{C}$	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- β	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
TNF- β	1.39 - 22.2	5.55	fg/mL	80% - 120%	< 20%	2.23	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TNF- β	1X Calibrator	30000	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 JUL 2023