



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human TNF- β Kit
Catalog Number: K151ADPS-Series; K156ADPS-Series
Lot Number: K00S0036
Expiration Date: 31 JUL 2022

Precision, Accuracy, and Sensitivity Specifications:

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

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

LLOD		
Name	Observed	Units
TNF- β	3.34	fg/mL

Stock Concentration of Calibrator:

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

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human TNF- β Antibody	D00T0044A	2-8°C	31 May 2024
Biotin Human TNF- β Antibody	B00T0044A	2-8°C	30 Apr 2024
Human TNF- β Calibrator	A00S0032	2-8°C	31 Aug 2025
Diluent 2	M0020112	\leq -10°C	31 Jul 2022
S-PLEX Coating Reagent C1	329998	\leq -70°C	31 Aug 2023
Blocker S1	R0140004	\leq -10°C	30 Apr 2023
S-PLEX Enhance E1	Y0550004	\leq -10°C	28 Feb 2023
S-PLEX Enhance E2	Y0560004	\leq -10°C	31 Jan 2023
S-PLEX Enhance E3	Y0570004	\leq -70°C	31 Jan 2023
S-PLEX Detect D1	Y0580006	\leq -70°C	31 Mar 2023
S-PLEX Detect D2	Y0590004	\leq -70°C	31 Jan 2023

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 APR 2023