



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151E3S-Series

Lot Number: K00S0006A

Expiration Date: 30 APR 2021

Precision, Accuracy, and Sensitivity Specifications:

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

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

Name	LLOD	
	Observed	Units
TNF- α	6.82	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TNF- α	5X Calibrator	376000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human TNF- α Antibody	D00T0007A	2-8°C	31 OCT 2022
Biotin Human TNF- α Antibody	B00T0007A	2-8°C	31 OCT 2022
Human TNF- α Calibrator	A00S0006	2-8°C	31 JUL 2024
Diluent 2	M0020107	\leq -10°C	31 JAN 2022
Diluent 3	M0030091	\leq -10°C	31 JAN 2022
S-PLEX Coating Reagent C1	265437	\leq -70°C	30 SEP 2021
Blocker S1	R0140001	\leq -10°C	31 MAY 2021
S-PLEX Enhance E1	Y0550001	\leq -10°C	30 APR 2021
S-PLEX Enhance E2	Y0560002	\leq -10°C	31 DEC 2021
S-PLEX Enhance E3	Y0570001	\leq -70°C	31 MAY 2021
S-PLEX Detect D1	Y0580001	\leq -70°C	30 APR 2021
S-PLEX Detect D2	Y0590001	\leq -70°C	30 APR 2021

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		19 MAY 2020