



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human CTLA-4 Kit

Catalog Number: K151AGLS – Series; K156AGLS - Series

Lot Number: K00S0055

Expiration Date: 28 Feb 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human CTLA-4 Detection Antibody	D00T0065A	2-8°C	30 Nov 2024
Biotin Human CTLA-4 Capture Antibody	B00T0065A	2-8°C	30 Nov 2024
Human CTLA-4 Calibrator	A00S0049	2-8°C	31 May 2026
Diluent 58	M0580001	≤ -10°C	28 Feb 2023
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 Aug 2023
Blocker S1	R0140004	≤ -10°C	30 Apr 2023
S-PLEX Enhance E1	Y0550006	≤ -10°C	31 Oct 2023
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 Oct 2023
S-PLEX Enhance E3	Y0570006	≤ -70°C	31 Oct 2023
S-PLEX Detect D1	Y0580008	≤ -70°C	30 Nov 2023
S-PLEX Detect D2	Y0590006	≤ -70°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
CTLA-4	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
CTLA-4	62.4 - 998	125	fg/mL	80% - 120%	< 20%	31.6	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
CTLA-4	1X Calibrator	560000	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		29 NOV 2021