



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human CTLA-4 Kit

**Catalog Number:** K151AGLS - Series

**Lot Number:** K00S0154

**Expiration Date:** 30 NOV 2024

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human CTLA-4 Antibody	D00T0065A	2-8°C	30 NOV 2024
Biotin Human CTLA-4 Antibody	B00T0065A	2-8°C	30 NOV 2024
Human CTLA-4 Calibrator	A00S0049	2-8°C	31 MAY 2026
Diluent 58	M0580009	≤ -10°C	30 NOV 2025
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
Blocker S1	R0010831	≤ -10°C	31 JUL 2025
S-PLEX Enhance E1	Y0550012	≤ -10°C	31 AUG 2025
S-PLEX Enhance E2	Y0560014	≤ -10°C	31 OCT 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 OCT 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 OCT 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
CTLA-4	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
CTLA-4	62.4 - 998	125	fg/mL	80% - 120%	< 20%	33.1	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
<b>Review/Approval</b>	Karen Hamilla	Quality		09 AUG 2023