

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Human IL-2 Kit  
**Catalog Number:** K151Z2S-Series  
**Lot Number:** K00S0164  
**Expiration Date:** 31 AUG 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-2 Antibody	B00T0173	2-8°C	31 JAN 2027
Human IL-2 Calibrator	A00S0091	2-8°C	31 OCT 2028
TURBO-BOOST <sup>™</sup> Human IL-2 Antibody	D00T0173	2-8°C	28 FEB 2027
Blocker S1	R0010871	≤ -10°C	31 MAR 2026
Diluent 2	M0020127	≤ -10°C	30 NOV 2025
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
S-PLEX Enhance E1	Y0550014	≤ -10°C	28 FEB 2026
S-PLEX Enhance E2	Y0560015	≤ -10°C	28 FEB 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580020	≤ -70°C	31 MAR 2026
S-PLEX Detect D2	Y0590016	≤ -70°C	30 APR 2026
S-PLEX Enhance E3	Y0570017	≤ -70°C	31 MAR 2026

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

**Precision, Accuracy, and Sensitivity Specifications:**

		Calibrator	
Spot Number	Name	Recovery	CV
1	IL-2	80% - 120%	< 20%

		LLOQ				LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-2	2.75 - 44.0	22.0	fg/mL	80% - 120%	< 20%	8.43	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
IL-2	30X Calibrator	1490000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		14 DEC 2023