

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Human IL-2 Kit  
**Catalog Number:** K151Z2S-Series; K156Z2S-Series  
**Lot Number:** K00S0181  
**Expiration Date:** 30 NOV 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-2 Antibody	B00T0186	2-8°C	30 JUN 2027
Human IL-2 Calibrator	A00S0091	2-8°C	31 OCT 2028
TURBO-BOOST <sup>™</sup> Human IL-2 Antibody	D00T0186	2-8°C	30 JUN 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 2	M0020127	≤ -10°C	30 NOV 2025
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560015	≤ -10°C	28 FEB 2026
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580019	≤ -70°C	31 DEC 2025
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E3	Y0570015	≤ -70°C	30 NOV 2025

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

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

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

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-2	2.75 - 44.0	22.0	fg/mL	80% - 120%	< 20%	7.76	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		03 APR 2024