



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-2 Kit
Lot Number: K00S0238
Expiration Date: 30 SEP 2026
Catalog Number: K156Z2S-Series, K151Z2S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-2 Antibody	B00T0238	2-8°C	30 JUN 2028
Human IL-2 Calibrator	A00S0091	2-8°C	31 OCT 2028
TURBO-BOOST Human IL-2 Antibody	D00T0238	2-8°C	31 AUG 2028
Blocker S1	R0140013	≤ -10°C	31 JUL 2027
Diluent 2	M0020130	≤ -10°C	30 SEP 2026
Diluent 3	M0030132	≤ -10°C	31 MAR 2027
S-PLEX Enhance E1	Y0550021	≤ -10°C	30 JUN 2027
S-PLEX Enhance E2	Y0560021	≤ -10°C	30 APR 2027
S-PLEX Coating Reagent C1	516221	≤ -70°C	31 OCT 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580031	≤ -70°C	31 JUL 2027
S-PLEX Detect D2	Y0590023	≤ -70°C	31 JUL 2027
S-PLEX Enhance E3	Y0570024	≤ -70°C	30 JUN 2027

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	44.0	fg/mL	80% - 120%	< 20%	5.64	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
Review/Approval	Karen Hamilla	Quality		09 APR 2025