



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-2 Kit
Catalog Number: K151Z2S-Series; K156Z2S-Series;
Lot Number: K00S0146
Expiration Date: 28 FEB 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL-2 Antibody	D00T0136	2-8°C	30 SEP 2026
Biotin Human IL-2 Antibody	B00T0136	2-8°C	31 JUL 2026
Human IL-2 Calibrator	A00S0069	2-8°C	31 JUL 2027
Diluent 2	M0020123	≤ -10°C	28 FEB 2025
Diluent 3	M0030121	≤ -10°C	31 OCT 2025
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
Blocker S1	R0010824	≤ -10°C	30 JUN 2025
S-PLEX Enhance E1	Y0550011	≤ -10°C	31 JUL 2025
S-PLEX Enhance E2	Y0560013	≤ -10°C	31 AUG 2025
S-PLEX Enhance E3	Y0570013	≤ -70°C	31 JUL 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580017	≤ -70°C	31 AUG 2025
S-PLEX Detect D2	Y0590013	≤ -70°C	31 JUL 2025

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

Precision, Accuracy, and Sensitivity Specifications:

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

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-2	2.75 - 44.0	11.0	fg/mL	80% - 120%	< 20%	4.06	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-2	30X Calibrator	1480000	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		20 JUN 2023