

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IL-2 Kit
Lot Number: K00S0313
Expiration Date: 30 SEP 2027
Catalog Number: K151Z2S-Series, K156Z2S-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-2 Antibody	B00T0284	2-8°C	31 JUL 2029
Human IL-2 Calibrator	A00S0134	2-8°C	30 APR 2030
TURBO-BOOST Human IL-2 Antibody	D00T0284	2-8°C	31 OCT 2029
Blocker S1	R0140022	≤ -10°C	31 OCT 2028
Diluent 2	M0020134	≤ -10°C	30 SEP 2027
Diluent 3	M0030149	≤ -10°C	31 AUG 2028
S-PLEX Enhance E1	Y0550028	≤ -10°C	31 MAR 2028
S-PLEX Enhance E2	Y0560028	≤ -10°C	31 MAY 2028
S-PLEX Coating Reagent C1	521331	≤ -70°C	31 OCT 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580037	≤ -70°C	30 APR 2028
S-PLEX Detect D2	Y0590029	≤ -70°C	31 MAY 2028
S-PLEX Enhance E3	Y0570030	≤ -70°C	31 MAR 2028

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	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-2	2.75 - 44.0	13.4	fg/mL	80% - 120%	< 20%	7.31	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-2	30X Calibrator	1350000	fg/mL

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 JUN 2026