



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-3 Kit**Lot Number:** K00S0289**Expiration Date** 31 MAY 2027**Catalog Number:** K151ADQS-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-3 Antibody	B00T0268	2-8°C	31 JAN 2029
Human IL-3 Calibrator	A00S0140	2-8°C	31 JUL 2030
TURBO-BOOST Human IL-3 Antibody	D00T0268	2-8°C	28 FEB 2029
Blocker S1	R0140015	≤ -10°C	31 OCT 2027
Diluent 58	M0580014	≤ -10°C	31 MAY 2027
S-PLEX Enhance E1	Y0550025	≤ -10°C	30 NOV 2027
S-PLEX Enhance E2	Y0560022	≤ -10°C	31 JUL 2027
S-PLEX Coating Reagent C1	521329	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580034	≤ -70°C	31 DEC 2027
S-PLEX Detect D2	Y0590025	≤ -70°C	30 NOV 2027
S-PLEX Enhance E3	Y0570027	≤ -70°C	31 DEC 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-3	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-3	10.6 - 170	10.6	fg/mL	80% - 120%	< 20%	2.92	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-3	1X Calibrator	40200	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		25 SEP 2025