



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] Human IL-13 Kit
Catalog Number: K151Y9S-Series
Lot Number: K00S0180
Expiration Date: 30 NOV 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-13 Antibody	B00T0185	2-8°C	30 JUN 2027
Human IL-13 Calibrator	A00S0103	2-8°C	31 MAR 2029
TURBO-BOOST™ Human IL-13 Antibody	D00T0185	2-8°C	31 JUL 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 2	M0020127	≤ -10°C	30 NOV 2025
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560017	≤ -10°C	31 JUL 2026
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580023	≤ -70°C	31 JUL 2026
S-PLEX Detect D2	Y0590018	≤ -70°C	31 JUL 2026
S-PLEX Enhance E3	Y0570018	≤ -70°C	30 JUN 2026

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-13	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-13	5.86 - 93.8	11.7	fg/mL	80% - 120%	< 20%	10.2	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-13	1X Calibrator	131000	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		21 MAY 2024