



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-13 Kit
Lot Number: K00S0263
Expiration Date: 30 JUN 2027
Catalog Number: K151Y9S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-13 Antibody	B00T0248	2-8°C	31 AUG 2028
Human IL-13 Calibrator	A00S0124	2-8°C	28 FEB 2030
TURBO-BOOST Human IL-13 Antibody	D00T0248	2-8°C	30 NOV 2028
Blocker S1	R0140016	≤ -10°C	31 DEC 2027
Diluent 2	M0020133	≤ -10°C	30 JUN 2027
S-PLEX Enhance E1	Y0550024	≤ -10°C	31 OCT 2027
S-PLEX Enhance E2	Y0560023	≤ -10°C	31 OCT 2027
S-PLEX Coating Reagent C1	521329	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580033	≤ -70°C	30 NOV 2027
S-PLEX Detect D2	Y0590026	≤ -70°C	31 DEC 2027
S-PLEX Enhance E3	Y0570026	≤ -70°C	31 OCT 2027

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	46.9	fg/mL	80% - 120%	< 20%	7.61	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-13	1X Calibrator	122000	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		31 OCT 2025