

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human IL-13 Kit
Lot Number: K00S0221
Expiration Date: 31 MAR 2026
Catalog Number: K151Y9S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-13 Antibody	B00T0228	2-8°C	29 FEB 2028
Human IL-13 Calibrator	A00S0118	2-8°C	31 AUG 2029
TURBO-BOOST Human IL-13 Antibody	D00T0228	2-8°C	31 DEC 2027
Blocker S1	R0140011	≤ -10°C	28 FEB 2027
Diluent 2	M0020128	≤ -10°C	31 MAR 2026
S-PLEX Enhance E1	Y0550018	≤ -10°C	30 NOV 2026
S-PLEX Enhance E2	Y0560019	≤ -10°C	30 NOV 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580027	≤ -70°C	31 OCT 2026
S-PLEX Detect D2	Y0590020	≤ -70°C	30 NOV 2026
S-PLEX Enhance E3	Y0570020	≤ -70°C	31 OCT 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	23.5	fg/mL	80% - 120%	< 20%	9.19	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-13	1X Calibrator	121000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Hamilla, Karen	Quality		25 OCT 2024