

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX® Human IL-13 Kit  
**Lot Number:** K00S0236  
**Expiration Date:** 31 MAR 2026  
**Catalog Number:** K151Y9S-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-13 Antibody	B00T0236	2-8°C	30 APR 2028
Human IL-13 Calibrator	A00S0103	2-8°C	31 MAR 2029
TURBO-BOOST Human IL-13 Antibody	D00T0236	2-8°C	31 MAY 2028
S-PLEX Coating Reagent C1	516221	≤ -70°C	31 OCT 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580029	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590022	≤ -70°C	30 APR 2027
S-PLEX Enhance E3	Y0570022	≤ -70°C	30 APR 2027
Blocker S1	R0140012	≤ -10°C	30 APR 2027
Diluent 2	M0020128	≤ -10°C	31 MAR 2026
S-PLEX Enhance E1	Y0550020	≤ -10°C	30 APR 2027
S-PLEX Enhance E2	Y0560020	≤ -10°C	31 MAR 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	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-13	5.86 - 93.8	23.5	fg/mL	80% - 120%	< 20%	9.84	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
<b>Review/Approval</b>	Michelle Likiam	Quality		23 DEC 2024