

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Human TSLP Kit  
**Catalog Number:** K151D3S-Series  
**Lot Number:** K00S0172  
**Expiration Date:** 31 AUG 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TSLP Antibody	B00T0179	2-8°C	30 JUN 2027
Human TSLP Calibrator	A00S0087	2-8°C	31 JUL 2028
TURBO-BOOST <sup>™</sup> Human TSLP Antibody	D00T0179	2-8°C	30 JUN 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 3	M0030125	≤ -10°C	31 MAY 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
S-PLEX Enhance E1	Y0550014	≤ -10°C	28 FEB 2026
S-PLEX Enhance E2	Y0560016	≤ -10°C	31 MAY 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580022	≤ -70°C	31 MAY 2026
S-PLEX Detect D2	Y0590015	≤ -70°C	31 JAN 2026
S-PLEX Enhance E3	Y0570015	≤ -70°C	30 NOV 2025

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	TSLP	80% - 120%	< 20%

		LLOQ				LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
TSLP	3.75 - 60.0	30.0	fg/mL	80% - 120%	< 20%	4.49	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
TSLP	7X Calibrator	462000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		28 FEB 2024