

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Human TSLP Kit  
**Lot Number:** K00S0218  
**Expiration Date:** 30 APR 2026  
**Catalog Number:** K151D3S-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TSLP Antibody	B00T0204	2-8°C	31 AUG 2027
Human TSLP Calibrator	A00S0112	2-8°C	31 JUL 2029
TURBO-BOOST Human TSLP Antibody	D00T0204	2-8°C	31 AUG 2027
Blocker S1	R0140008	≤ -10°C	31 JUL 2026
Diluent 3	M0030128	≤ -10°C	31 OCT 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560018	≤ -10°C	31 AUG 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> Reagent	Y0580025	≤ -70°C	30 SEP 2026
S-PLEX Detect D2	Y0590018	≤ -70°C	31 JUL 2026
S-PLEX Enhance E3	Y0570019	≤ -70°C	31 JUL 2026

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	15.0	fg/mL	80% - 120%	< 20%	8.39	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
TSLP	7X Calibrator	525000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Hamilla, Karen	Quality		16 SEP 2024