

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human TSLP Kit
Lot Number: K00S0226
Expiration Date: 31 AUG 2025
Catalog Number: K151D3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TSLP Antibody	B00T0179	2-8°C	30 JUN 2027
Human TSLP Antibody	D00T0179	2-8°C	30 JUN 2027
Human TSLP Calibrator	A00S0112	2-8°C	31 JUL 2029
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 [™] 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:

Spot Number	Name	Calibrator	
		Recovery	CV
1	TSLP	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
TSLP	3.75 - 60.0	30.0	fg/mL	80% - 120%	< 20%	6.99	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
Review/Approval	Karen Hamilla	Quality		16 SEP 2024