

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human TSLP Kit
Lot Number: K00S0231
Expiration Date: 30 SEP 2026
Catalog Number: K151D3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TSLP Antibody	B00T0234	2-8°C	31 MAR 2028
Human TSLP Calibrator	A00S0120	2-8°C	31 OCT 2029
TURBO-BOOST Human TSLP Antibody	D00T0234	2-8°C	31 MAR 2028
Blocker S1	R0140011	≤ -10°C	28 FEB 2027
Diluent 3	M0030130	≤ -10°C	31 JAN 2027
Diluent 43	M0430056	≤ -10°C	30 SEP 2026
S-PLEX Enhance E1	Y0550019	≤ -10°C	31 MAR 2027
S-PLEX Enhance E2	Y0560019	≤ -10°C	30 NOV 2026
S-PLEX Coating Reagent C1	516221	≤ -70°C	31 OCT 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580028	≤ -70°C	31 JAN 2027
S-PLEX Detect D2	Y0590020	≤ -70°C	30 NOV 2026
S-PLEX Enhance E3	Y0570021	≤ -70°C	30 NOV 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	30.0	fg/mL	80% - 120%	< 20%	7.32	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TSLP	7X Calibrator	580000	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		13 DEC 2024