



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human TSLP Kit

Catalog Number: K151D3S-Series

Lot Number: K00S0150

Expiration Date: 31 JUL 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human TSLP Antibody	D00T0155	2-8°C	31 DEC 2026
Biotin Human TSLP Antibody	B00T0155	2-8°C	31 DEC 2026
Human TSLP Calibrator	A00S0087	2-8°C	31 JUL 2028
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
Diluent 43	M0430053	≤ -10°C	31 JAN 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
Blocker S1	R0010831	≤ -10°C	31 JUL 2025
S-PLEX Enhance E1	Y0550012	≤ -10°C	31 AUG 2025
S-PLEX Enhance E2	Y0560014	≤ -10°C	31 OCT 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 OCT 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 OCT 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
TSLP	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
TSLP	3.75 – 60.0	15.0	fg/mL	80% - 120%	< 20%	4.87	fg/mL

Stock Concentration of Calibrator:

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 OCT 2023