



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** S-PLEX® Human TSLP Kit


**Catalog Number(s) / Size(s):** K151D3S - Series

**Lot Number:** K00S0119

Based on additional testing, the original date of expiration for this kit has been extended by 5 months. The expiration date of K00S0119 has been extended to 31 JAN 2024.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
S-PLEX Coating Reagent C1	329998	31 AUG 2023	31 JAN 2024
S-PLEX Enhance E2	Y0560007	31 OCT 2023	31 JAN 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		21 JUL 2023

This memo supersedes all other statements regarding date of expiration for this product.



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX® Human TSLP Kit

**Catalog Number:** K151D3S-Series

**Lot Number:** K00S0119

**Expiration Date:** 31 AUG 2023

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human TSLP Antibody	D00T0120	2-8°C	31 MAY 2026
Biotin Human TSLP Antibody	B00T0120	2-8°C	30 APR 2026
Human TSLP Calibrator	A00S0047	2-8°C	30 APR 2026
Diluent 3	M0030116	≤ -10°C	30 NOV 2024
Diluent 43	M0430047	≤ -10°C	31 MAY 2024
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 AUG 2023
Blocker S1	R0010768	≤ -10°C	31 MAY 2024
S-PLEX Enhance E1	Y0550009	≤ -10°C	28 FEB 2025
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 OCT 2023
S-PLEX Enhance E3	Y0570011	≤ -70°C	31 MAR 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580015	≤ -70°C	30 APR 2025
S-PLEX Detect D2	Y0590010	≤ -70°C	28 FEB 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	30.0	fg/mL	80% - 120%	< 20%	13.2	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TSLP	7X Calibrator	429000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		12 DEC 2022