

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human TSLP Kit, SECTOR
Lot Number: K00S0297
Expiration Date: 28 FEB 2027
Catalog Number: K151D3S-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human TSLP Antibody	B00T0275	2-8°C	30 APR 2029
Human TSLP Calibrator	A00S0144	2-8°C	30 SEP 2030
TURBO-BOOST Human TSLP Antibody	D00T0275	2-8°C	31 MAY 2029
Blocker S1	R0140018	≤ -10°C	30 APR 2028
Diluent 3	M0030145	≤ -10°C	29 FEB 2028
Diluent 43	M0430058	≤ -10°C	28 FEB 2027
S-PLEX Enhance E1	Y0550026	≤ -10°C	31 JAN 2028
S-PLEX Enhance E2	Y0560026	≤ -10°C	29 FEB 2028
S-PLEX Coating Reagent C1	521330	≤ -70°C	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580036	≤ -70°C	30 APR 2028
S-PLEX Detect D2	Y0590028	≤ -70°C	29 FEB 2028
S-PLEX Enhance E3	Y0570029	≤ -70°C	29 FEB 2028

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	10.9	fg/mL	80% - 120%	< 20%	7.09	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TSLP	7X Calibrator	494000	fg/mL

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 MAR 2026