



**RFE MEMO**

Page 1 of 1

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

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

**Lot Number:** K00S0010A

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0010A has been extended to 30 APR 2022.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
S-PLEX Coating Reagent C1	265437	30 SEP 2021	30 APR 2022
Blocker S1	R0140001	31 MAY 2021	30 APR 2022
S-PLEX Enhance E1	Y0550002	31 DEC 2021	30 APR 2022
S-PLEX Enhance E2	Y0560002	31 DEC 2021	30 APR 2022
S-PLEX Enhance E3	Y0570001	31 MAY 2021	30 APR 2022
S-PLEX Detect D1	Y0580001	30 APR 2021	30 APR 2022
S-PLEX Detect D2	Y0590001	30 APR 2021	30 APR 2022
Diluent 43	M0430033	31 JAN 2022	30 APR 2022
Diluent 3	M0030091	31 JAN 2022	30 APR 2022

Function	Name	Signature	Date
Quality	Karen Hamilla		13 MAY 2021

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



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human TSLP Kit

**Catalog Number:** K151D3S-Series

**Lot Number:** K00S0010A

**Expiration Date:** 30 APR 2021

### Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				
	Range Tested	Observed	Units	Recovery	CV
TSLP	3.75 - 60.0	30.0	fg/mL	80% - 120%	< 20%

Name	LLOD	
	Observed	Units
TSLP	9.05	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
TSLP	7X Calibrator	488000	fg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human TSLP Antibody	D00T0006A	2-8°C	31 OCT 2022
Biotin Human TSLP Antibody	B00T0006A	2-8°C	31 OCT 2022
Human TSLP Calibrator	A00S0011	2-8°C	31 JUL 2024
Diluent 43	M0430033	≤ -10°C	31 JAN 2022
Diluent 3	M0030091	≤ -10°C	31 JAN 2022
S-PLEX Coating Reagent C1	265437	≤ -70°C	30 SEP 2021
Blocker S1	R0140001	≤ -10°C	31 MAY 2021
S-PLEX Enhance E1	Y0550002	≤ -10°C	31 DEC 2021
S-PLEX Enhance E2	Y0560002	≤ -10°C	31 DEC 2021
S-PLEX Enhance E3	Y0570001	≤ -70°C	31 MAY 2021
S-PLEX Detect D1	Y0580001	≤ -70°C	30 APR 2021
S-PLEX Detect D2	Y0590001	≤ -70°C	30 APR 2021

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 JUL 2020