



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® Human TSLP Kit

Catalog Number: K151D3S-Series

Lot Number: K00S0028

Expiration Date: 28 FEB 2023

Precision, Accuracy, and Sensitivity Specifications:

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

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

| LLOD | | |
|------|----------|-------|
| Name | Observed | Units |
| TSLP | 7.77 | 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™ Human TSLP Antibody | D00T0032A | 2-8°C | 30 Nov 2023 |
| Biotin Human TSLP Antibody | B00T0032A | 2-8°C | 31 Mar 2024 |
| Human TSLP Calibrator | A00S0011 | 2-8°C | 31 Jul 2024 |
| Diluent 43 | M0430044 | ≤ -10°C | 30 Sep 2023 |
| Diluent 3 | M0030106 | ≤ -10°C | 31 Aug 2023 |
| S-PLEX Coating Reagent C1 | 329998 | ≤ -70°C | 31 Aug 2023 |
| Blocker S1 | R0140004 | ≤ -10°C | 30 Apr 2023 |
| S-PLEX Enhance E1 | Y0550005 | ≤ -10°C | 31 May 2023 |
| S-PLEX Enhance E2 | Y0560006 | ≤ -10°C | 31 May 2023 |
| S-PLEX Enhance E3 | Y0570005 | ≤ -70°C | 28 Feb 2023 |
| S-PLEX Detect D1 | Y0580006 | ≤ -70°C | 31 Mar 2023 |
| S-PLEX Detect D2 | Y0590005 | ≤ -70°C | 28 Feb 2023 |

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 | | 24 JUN 2021 |