



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-2 Kit
Catalog Number: K151Z2S-Series; K156Z2S-Series
Lot Number: K00S0047
Expiration Date: 30 APR 2023

Precision, Accuracy, and Sensitivity Specifications:

| Calibrator | | |
|------------|------------|-------|
| Name | Recovery | CV |
| IL-2 | 80% - 120% | < 20% |

| LLOQ | | | | | |
|------|--------------|----------|-------|------------|-------|
| Name | Range Tested | Observed | Units | Recovery | CV |
| IL-2 | 2.75 - 44.0 | 22.0 | fg/mL | 80% - 120% | < 20% |

| LLOD | | |
|------|----------|-------|
| Name | Observed | Units |
| IL-2 | 5.14 | fg/mL |

Stock Concentration of Calibrator:

| Name | Item | Concentration | Units |
|------|----------------|---------------|-------|
| IL-2 | 30X Calibrator | 1360000 | fg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|----------------------------------|------------|---------------------|-----------------|
| TURBO-BOOST™ Human IL-2 Antibody | D00T0049A | 2-8°C | 30 Sep 2024 |
| Biotin Human IL-2 Antibody | B00T0049A | 2-8°C | 30 Apr 2024 |
| Human IL-2 Calibrator | A00S0053 | 2-8°C | 31 Jul 2026 |
| Diluent 2 | M0020117 | ≤ -10°C | 31 Oct 2023 |
| Diluent 3 | M0030109 | ≤ -10°C | 31 Jan 2024 |
| S-PLEX Coating Reagent C1 | 329998 | ≤ -70°C | 31 Aug 2023 |
| Blocker S1 | R0140004 | ≤ -10°C | 30 Apr 2023 |
| S-PLEX Enhance E1 | R0140006 | ≤ -10°C | 31 Jan 2024 |
| S-PLEX Enhance E2 | Y0560006 | ≤ -10°C | 31 May 2023 |
| S-PLEX Enhance E3 | Y0570007 | ≤ -70°C | 29 Feb 2024 |
| S-PLEX Detect D1 | Y0580008 | ≤ -70°C | 30 Nov 2023 |
| S-PLEX Detect D2 | Y0590006 | ≤ -70°C | 31 Oct 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 | | 05 NOV 2021 |