



RFE MEMO

Page 1 of 1

Product Description(s): S-PLEX® Human IL-6 Kit
Catalog Number(s) / Size(s): K151B3S-Series; K156B3S-Series
Lot Number: K00S0019

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0019 has been extended to 31 DEC 2022.

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

| Description | Lot Number | Original Expiration Date | New Expiration Date |
|-------------------|------------|--------------------------|---------------------|
| Diluent 2 | M0020113 | 30 NOV 2022 | 31 DEC 2022 |
| Blocker S1 | R0140003 | 31 AUG 2022 | 31 DEC 2022 |
| S-PLEX Enhance E1 | Y0550003 | 30 SEP 2022 | 31 DEC 2022 |
| S-PLEX Enhance E2 | Y0560003 | 31 MAR 2022 | 31 DEC 2022 |
| S-PLEX Enhance E3 | Y0570002 | 31 DEC 2021 | 31 DEC 2022 |
| S-PLEX Detect D2 | Y0590003 | 31 AUG 2022 | 31 DEC 2022 |

| Function | Name | Signature | Date |
|----------|---------------|-----------|-------------|
| Quality | Karen Hamilla | | 06 DEC 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[®] Human IL-6 Kit
Catalog Number: K151B3S-Series; K156B3S-Series
Lot Number: K00S0019
Expiration Date: 31 DEC 2021

Precision, Accuracy, and Sensitivity Specifications:

| Calibrator | | |
|------------|------------|-------|
| Name | Recovery | CV |
| IL-6 | 80% - 120% | < 20% |

| LLOQ | | | | | |
|------|--------------|----------|-------|------------|-------|
| Name | Range Tested | Observed | Units | Recovery | CV |
| IL-6 | 0.33 – 5.2 | 1.3 | fg/mL | 80% - 120% | < 20% |

| LLOD | | |
|------|----------|-------|
| Name | Observed | Units |
| IL-6 | 0.35 | fg/mL |

Stock Concentration of Calibrator:

| Name | Item | Concentration | Units |
|------|-----------------|---------------|-------|
| IL-6 | 120X Calibrator | 768000 | fg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|----------------------------------|------------|---------------------|-----------------|
| TURBO-BOOST™ Human IL-6 Antibody | D00T0026A | 2-8°C | 30 SEP 2023 |
| Biotin Human IL-6 Antibody | B00T0026A | 2-8°C | 30 SEP 2023 |
| Human IL-6 Calibrator | A00S0021 | 2-8°C | 30 JUN 2025 |
| Diluent 2 | M0020113 | ≤ -10°C | 30 NOV 2022 |
| Diluent 3 | M0030102 | ≤ -10°C | 31 JAN 2023 |
| S-PLEX Coating Reagent C1 | 314374 | ≤ -70°C | 28 FEB 2023 |
| Blocker S1 | R0140003 | ≤ -10°C | 31 AUG 2022 |
| S-PLEX Enhance E1 | Y0550003 | ≤ -10°C | 30 SEP 2022 |
| S-PLEX Enhance E2 | Y0560003 | ≤ -10°C | 31 MAR 2022 |
| S-PLEX Enhance E3 | Y0570002 | ≤ -70°C | 31 DEC 2021 |
| S-PLEX Detect D1 | Y0580004 | ≤ -70°C | 28 FEB 2023 |
| S-PLEX Detect D2 | Y0590003 | ≤ -70°C | 31 AUG 2022 |

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 OCT 2020 |