



EXTENSION MEMO

Page 1 of 1

Product Description(s): S-PLEX® Human TNF- α Kit


Catalog Number(s) / Size(s): K151E3S-Series

Lot Number: K00S0121

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0121 has been extended to 30 APR 2025.

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

| Description | Lot Number | Original Expiration Date | New Expiration Date |
|---------------------------|------------|--------------------------|---------------------|
| Diluent 2 | M0020122 | 31 OCT 2024 | 30 APR 2025 |
| S-PLEX Coating Reagent C1 | 351987 | 30 APR 2024 | 30 APR 2025 |

| Function | Name | Signature | Date |
|----------|---------------|---|-------------|
| Quality | Karen Hamilla |  | 08 Feb 2024 |

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 TNF- α Kit

Catalog Number: K151E3S-Series

Lot Number: K00S0121

Expiration Date: 30 APR 2024

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------|----------------------------|-----------------|
| TURBO-BOOST [™] Human TNF- α Antibody | D00T0122 | 2-8°C | 30 JUN 2026 |
| Biotin Human TNF- α Antibody | B00T0122 | 2-8°C | 31 MAY 2026 |
| Human TNF- α Calibrator | A00S0079 | 2-8°C | 31 DEC 2027 |
| Diluent 2 | M0020122 | $\leq -10^{\circ}\text{C}$ | 31 OCT 2024 |
| Diluent 3 | M0030119 | $\leq -10^{\circ}\text{C}$ | 30 APR 2025 |
| S-PLEX Coating Reagent C1 | 351987 | $\leq -70^{\circ}\text{C}$ | 30 APR 2024 |
| Blocker S1 | R0010824 | $\leq -10^{\circ}\text{C}$ | 30 JUN 2025 |
| S-PLEX Enhance E1 | Y0550010 | $\leq -10^{\circ}\text{C}$ | 31 MAY 2025 |
| S-PLEX Enhance E2 | Y0560012 | $\leq -10^{\circ}\text{C}$ | 31 MAY 2025 |
| S-PLEX Enhance E3 | Y0570012 | $\leq -70^{\circ}\text{C}$ | 31 MAY 2025 |
| S-PLEX Detect D1 with TURBO-TAG [™] reagent | Y0580016 | $\leq -70^{\circ}\text{C}$ | 31 MAY 2025 |
| S-PLEX Detect D2 | Y0590012 | $\leq -70^{\circ}\text{C}$ | 30 APR 2025 |

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

Precision, Accuracy, and Sensitivity Specifications:

| Name | Calibrator | |
|---------------|------------|-------|
| | Recovery | CV |
| TNF- α | 80% - 120% | < 20% |

| Name | Range Tested | LLOQ | | | | LLOD | |
|---------------|--------------|----------|-------|------------|-------|----------|-------|
| | | Observed | Units | Recovery | CV | Observed | Units |
| TNF- α | 6.25 - 100 | 25.0 | fg/mL | 80% - 120% | < 20% | 6.37 | fg/mL |

Stock Concentration of Calibrator:

| Name | Item | Concentration | Units |
|---------------|---------------|---------------|-------|
| TNF- α | 5X Calibrator | 408000 | fg/mL |

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

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 29 MAR 2023 |