



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IFN- α 2a Kit

Catalog Number: K151P3S-Series

Lot Number: K00S0103

Expiration Date: 31 AUG 2023

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|----------------------------|-----------------|
| TURBO-BOOST [™] Human IFN- α 2a Antibody | D00T0099 | 2-8°C | 31 MAR 2026 |
| Biotin Human IFN- α 2a Antibody | B00T0099 | 2-8°C | 31 MAR 2026 |
| Human IFN- α 2a Calibrator | A00S0058 | 2-8°C | 30 NOV 2026 |
| Diluent 43 | M0430049 | $\leq -10^{\circ}\text{C}$ | 31 OCT 2024 |
| Diluent 3 | M0030117 | $\leq -10^{\circ}\text{C}$ | 31 JAN 2025 |
| S-PLEX Coating Reagent C1 | 329998 | $\leq -70^{\circ}\text{C}$ | 31 AUG 2023 |
| Blocker S1 | R0010811 | $\leq -10^{\circ}\text{C}$ | 28 FEB 2025 |
| S-PLEX Enhance E1 | Y0550010 | $\leq -10^{\circ}\text{C}$ | 31 MAY 2025 |
| S-PLEX Enhance E2 | Y0560011 | $\leq -10^{\circ}\text{C}$ | 31 MAR 2025 |
| S-PLEX Enhance E3 | Y0570011 | $\leq -70^{\circ}\text{C}$ | 31 MAR 2025 |
| S-PLEX Detect D1 with TURBO-TAG [™] reagent | Y0580015 | $\leq -70^{\circ}\text{C}$ | 30 APR 2025 |
| S-PLEX Detect D2 | Y0590011 | $\leq -70^{\circ}\text{C}$ | 31 MAR 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 |
| IFN- α 2a | 80% - 120% | < 20% |

| Name | LLOQ | | | | | LLOD | |
|------------------|--------------|----------|-------|------------|-------|----------|-------|
| | Range Tested | Observed | Units | Recovery | CV | Observed | Units |
| IFN- α 2a | 3.50 - 56.0 | 28.0 | fg/mL | 80% - 120% | < 20% | 4.45 | fg/mL |

Stock Concentration of Calibrator:

| Name | Item | Concentration | Units |
|------------------|-----------------|---------------|-------|
| IFN- α 2a | 750X Calibrator | 51000000 | 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 | | 17 JAN 2023 |