



## CERTIFICATE OF ANALYSIS

**Product Description:** S-PLEX Human IFN- $\alpha$ 2a Kit  
**Lot Number:** K00S0205  
**Expiration Date:** 30 SEP 2026  
**Catalog Number:** K151P3S-Series

**Lot-Specific Components:**

| Description                                 | Lot Number | Storage Temperature        | Expiration Date |
|---------------------------------------------|------------|----------------------------|-----------------|
| Biotin Human IFN- $\alpha$ 2a Antibody      | B00T0218   | 2-8°C                      | 31 OCT 2027     |
| Human IFN- $\alpha$ 2a Calibrator           | A00S0109   | 2-8°C                      | 31 AUG 2029     |
| TURBO-BOOST Human IFN- $\alpha$ 2a Antibody | D00T0218   | 2-8°C                      | 30 NOV 2027     |
| S-PLEX Coating Reagent C1                   | 476750     | $\leq -70^{\circ}\text{C}$ | 31 MAR 2027     |
| S-PLEX Detect D1 with TURBO-TAG™ Reagent    | Y0580028   | $\leq -70^{\circ}\text{C}$ | 31 JAN 2027     |
| S-PLEX Detect D2                            | Y0590021   | $\leq -70^{\circ}\text{C}$ | 31 MAR 2027     |
| S-PLEX Enhance E3                           | Y0570021   | $\leq -70^{\circ}\text{C}$ | 30 NOV 2026     |
| Blocker S1                                  | R0140010   | $\leq -10^{\circ}\text{C}$ | 31 DEC 2026     |
| Diluent 3                                   | M0030129   | $\leq -10^{\circ}\text{C}$ | 30 NOV 2026     |
| Diluent 43                                  | M0430056   | $\leq -10^{\circ}\text{C}$ | 30 SEP 2026     |
| S-PLEX Enhance E1                           | Y0550019   | $\leq -10^{\circ}\text{C}$ | 31 MAR 2027     |
| S-PLEX Enhance E2                           | Y0560020   | $\leq -10^{\circ}\text{C}$ | 31 MAR 2027     |

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

**Precision, Accuracy, and Sensitivity Specifications:**

|             |                  | Calibrator |       |
|-------------|------------------|------------|-------|
| Spot Number | Name             | Recovery   | CV    |
| 1           | IFN- $\alpha$ 2a | 80% - 120% | < 20% |

| LLOQ             |              |          |       |            |       | LLOD     |       |
|------------------|--------------|----------|-------|------------|-------|----------|-------|
| Name             | Range Tested | Observed | Units | Recovery   | CV    | Observed | Units |
| IFN- $\alpha$ 2a | 3.50 - 56.0  | 56.0     | fg/mL | 80% - 120% | < 20% | 7.03     | fg/mL |

**Stock Concentration of Calibrator:**

| Name             | Item            | Concentration | Units |
|------------------|-----------------|---------------|-------|
| IFN- $\alpha$ 2a | 750X Calibrator | 50500000      | fg/mL |

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

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 11 FEB 2025 |