



## CERTIFICATE OF ANALYSIS

**Product Description:** S-PLEX Human IL-4 Kit  
**Lot Number:** K00S0265  
**Expiration Date:** 30 JUN 2027  
**Catalog Number:** K151A3S-Series

**Lot-Specific Components:**

| Description                              | Lot Number | Storage Temperature | Expiration Date |
|------------------------------------------|------------|---------------------|-----------------|
| Biotin Human IL-4Antibody                | B00T0250   | 2-8°C               | 31 AUG 2028     |
| TURBO-BOOST Human IL-4 Antibody          | D00T0250   | 2-8°C               | 30 NOV 2028     |
| Human IL-4 Calibrator                    | A00S0111   | 2-8°C               | 31 JAN 2030     |
| Blocker S1                               | R0140015   | ≤ -10°C             | 31 OCT 2027     |
| Diluent 2                                | M0020133   | ≤ -10°C             | 30 JUN 2027     |
| Diluent 3                                | M0030140   | ≤ -10°C             | 30 SEP 2027     |
| S-PLEX Enhance E1                        | Y0550024   | ≤ -10°C             | 31 OCT 2027     |
| S-PLEX Enhance E2                        | Y0560023   | ≤ -10°C             | 31 OCT 2027     |
| S-PLEX Coating Reagent C1                | 521329     | ≤ -70°C             | 31 DEC 2027     |
| S-PLEX Detect D1 with TURBO-TAG™ Reagent | Y0580033   | ≤ -70°C             | 30 NOV 2027     |
| S-PLEX Detect D2                         | Y0590025   | ≤ -70°C             | 30 NOV 2027     |
| S-PLEX Enhance E3                        | Y0570027   | ≤ -70°C             | 31 DEC 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           | IL-4 | 80% - 120% | < 20% |

| LLOQ |              |          |       |            |       | LLOD     |       |
|------|--------------|----------|-------|------------|-------|----------|-------|
| Name | Range Tested | Observed | Units | Recovery   | CV    | Observed | Units |
| IL-4 | 0.306 - 4.90 | 2.45     | fg/mL | 80% - 120% | < 20% | 1.03     | fg/mL |

**Stock Concentration of Calibrator:**

| Name | Item           | Concentration | Units |
|------|----------------|---------------|-------|
| IL-4 | 20X Calibrator | 125000        | 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  |  | 23 SEP 2025 |