



**RFE MEMO**

Page 1 of 1

**Product Description(s):** S-PLEX® Human IL-13 Kit


**Catalog Number(s) / Size(s):** K151Y9S-Series

**Lot Number:** K00S0033

Based on additional testing, the original date of expiration for this kit has been extended by 10 months. The expiration date of K00S0033 has been extended to 31 MAR 2022.

The expiration date for the following kit component has been extended:

| Description | Lot Number | Original Expiration Date | New Expiration Date |
|-------------|------------|--------------------------|---------------------|
| Blocker S1  | R0140001   | 31 MAY 2021              | 31 MAR 2022         |

| Function | Name          | Signature  | Date        |
|----------|---------------|--|-------------|
| Quality  | Karen Hamilla |  | 07 JUN 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<sup>®</sup> Human IL-13 Kit

**Catalog Number:** K151Y9S-Series

**Lot Number:** K00S0033

**Expiration Date:** 31 MAY 2021

### Precision, Accuracy, and Sensitivity Specifications:

| Calibrator |            |       |
|------------|------------|-------|
| Name       | Recovery   | CV    |
| IL-13      | 80% - 120% | < 20% |

| LLOQ  |              |          |       |            |       |
|-------|--------------|----------|-------|------------|-------|
| Name  | Range Tested | Observed | Units | Recovery   | CV    |
| IL-13 | 5.86 - 93.8  | 46.9     | fg/mL | 80% - 120% | < 20% |

| LLOD  |          |       |
|-------|----------|-------|
| Name  | Observed | Units |
| IL-13 | 9.10     | fg/mL |

### Stock Concentration of Calibrator:

| Name  | Item          | Concentration | Units |
|-------|---------------|---------------|-------|
| IL-13 | 1X Calibrator | 128000        | fg/mL |

### Lot-Specific Components:

| Description                       | Lot Number | Storage Temperature | Expiration Date |
|-----------------------------------|------------|---------------------|-----------------|
| TURBO-BOOST™ Human IL-13 Antibody | D00T0038A  | 2-8°C               | 31 Mar 2024     |
| Biotin Human IL-13 Antibody       | B00T0038A  | 2-8°C               | 29 Feb 2024     |
| Human IL-13 Calibrator            | A00S0027   | 2-8°C               | 31 Aug 2025     |
| Diluent 2                         | M0020112   | ≤ -10°C             | 31 Jul 2022     |
| S-PLEX Coating Reagent C1         | 314374     | ≤ -70°C             | 28 Feb 2023     |
| Blocker S1                        | R0140001   | ≤ -10°C             | 31 May 2021     |
| S-PLEX Enhance E1                 | Y0550003   | ≤ -10°C             | 30 Sep 2022     |
| S-PLEX Enhance E2                 | Y0560003   | ≤ -10°C             | 31 Mar 2022     |
| S-PLEX Enhance E3                 | Y0570003   | ≤ -70°C             | 31 Dec 2022     |
| S-PLEX Detect D1                  | Y0580005   | ≤ -70°C             | 31 Mar 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        |
|------------------------|---------------|----------|-----------|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality  |           | 07 JAN 2021 |