



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-6 Kit  
**Catalog Number:** K151B3S-Series, K156B3S-Series  
**Lot Number:** K00S0003A  
**Expiration Date:** 30 APR 2021

#### Precision, Accuracy, and Sensitivity Specifications:

| Name | Calibrator |       |
|------|------------|-------|
|      | Recovery   | CV    |
| IL-6 | 80% - 120% | < 20% |

| Name | LLOQ         |          |       |            |       |
|------|--------------|----------|-------|------------|-------|
|      | Range Tested | Observed | Units | Recovery   | CV    |
| IL-6 | 0.32 - 5.2   | 1.3      | fg/mL | 80% - 120% | < 20% |

| Name | LLOD     |       |
|------|----------|-------|
|      | Observed | Units |
| IL-6 | 1.07     | fg/mL |

#### Stock Concentration of Calibrator:

| Name | Item            | Concentration | Units |
|------|-----------------|---------------|-------|
| IL-6 | 120X Calibrator | 720000        | fg/mL |

#### Lot-Specific Components:

| Description                                  | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| TURBO-BOOST <sup>™</sup> Human IL-6 Antibody | D00T0004A  | 2-8°C               | 31 OCT 2022     |
| Biotin Human IL-6 Antibody                   | B00T0004A  | 2-8°C               | 31 OCT 2022     |
| Human IL-6 Calibrator                        | A00S0003   | 2-8°C               | 31 JUL 2024     |
| Diluent 2                                    | M0020107   | ≤ -10°C             | 31 JAN 2022     |
| Diluent 3                                    | M0030091   | ≤ -10°C             | 31 JAN 2022     |
| S-PLEX Coating Reagent C1                    | 265437     | ≤ -70°C             | 30 SEP 2021     |
| Blocker S1                                   | R0140001   | ≤ -10°C             | 31 MAY 2021     |
| S-PLEX Enhance E1                            | Y0550001   | ≤ -10°C             | 30 APR 2021     |
| S-PLEX Enhance E2                            | Y0560001   | ≤ -10°C             | 30 APR 2021     |
| S-PLEX Enhance E3                            | Y0570001   | ≤ -70°C             | 31 MAY 2021     |
| S-PLEX Detect D1                             | Y0580001   | ≤ -70°C             | 30 APR 2021     |
| S-PLEX Detect D2                             | Y0590001   | ≤ -70°C             | 30 APR 2021     |

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        |
|-----------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality  |           | 22 JUN 2020 |