



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human GM-CSF Kit

Catalog Number: K151F3S-Series

Lot Number: K00S0007A

Expiration Date: 30 APR 2021

Precision, Accuracy, and Sensitivity Specifications:

| Name | Calibrator | |
|--------|------------|-------|
| | Recovery | CV |
| GM-CSF | 80% - 120% | < 20% |

| Name | LLOQ | | | | |
|--------|--------------|----------|-------|------------|-------|
| | Range Tested | Observed | Units | Recovery | CV |
| GM-CSF | 0.55 - 8.80 | 4.40 | fg/mL | 80% - 120% | < 20% |

| Name | LLOD | |
|--------|----------|-------|
| | Observed | Units |
| GM-CSF | 1.90 | fg/mL |

Stock Concentration of Calibrator:

| Name | Item | Concentration | Units |
|--------|----------------|---------------|-------|
| GM-CSF | 50X Calibrator | 1300000 | fg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|------------------------------------|------------|---------------------|-----------------|
| TURBO-BOOST™ Human GM-CSF Antibody | D00T0008A | 2-8°C | 31 OCT 2022 |
| Biotin Human GM-CSF Antibody | B00T0008A | 2-8°C | 31 OCT 2022 |
| Human GM-CSF Calibrator | A00S0007 | 2-8°C | 31 JUL 2024 |
| Diluent 43 | M0430032 | ≤ -10°C | 30 SEP 2021 |
| 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 | Y0560002 | ≤ -10°C | 31 DEC 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 FEB 2021 |