

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Neurology Panel 1 (human) Kit
Lot Number: K00S0198
Expiration Date: 31 JUL 2026
Catalog Number: K15639S-Series; K15640S-Series; K151AKGS-Series; K156AKGS-Series;
 K151AMPS-Series; K156AMPS-Series; K151APSS-Series; K156APSS-Series

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| TURBO-BOOST Hu Neurofilament L Antibody | D00T0207 | 2-8°C | 30 NOV 2027 |
| TURBO-BOOST Human GFAP Antibody | D00T0206 | 2-8°C | 31 AUG 2027 |
| TURBO-BOOST Human Tau Antibody | D00T0208 | 2-8°C | 31 OCT 2027 |
| Blocker S1 | R0140008 | ≤ -10°C | 31 JUL 2026 |
| Blocker S2 | 476747 | ≤ -10°C | 30 APR 2029 |
| S-PLEX Enhance E1 | Y0550017 | ≤ -10°C | 31 AUG 2026 |
| S-PLEX Enhance E2 | Y0560018 | ≤ -10°C | 31 AUG 2026 |
| S-PLEX Detect D1 with TURBO-TAG [™] Reagent | Y0580025 | ≤ -70°C | 30 SEP 2026 |
| S-PLEX Detect D2 | Y0590019 | ≤ -70°C | 30 SEP 2026 |
| S-PLEX Enhance E3 | Y0570019 | ≤ -70°C | 31 JUL 2026 |
| S-PLEX Neurology Panel 1 (human) Calibrator | A00S0106 | ≤ -70°C | 30 SEP 2026 |
| S-PLEX Neurology Panel 1 (human) Coating Solution | 485774 | ≤ -70°C | 31 AUG 2027 |

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

Precision, Accuracy, and Sensitivity Specifications:

| Spot Number | Name | Calibrator | |
|-------------|------------|------------|-------|
| | | Recovery | CV |
| 1 | GFAP | 80% - 120% | < 20% |
| 3 | NF-L | 80% - 120% | < 20% |
| 10 | Tau(total) | 80% - 120% | < 20% |

| Name | Range Tested | Observed | LLOQ | | | LLOD | |
|------------|--------------|----------|-------|------------|-------|----------|-------|
| | | | Units | Recovery | CV | Observed | Units |
| GFAP | 85.0 - 20300 | 464 | fg/mL | 80% - 120% | < 25% | 161 | fg/mL |
| NF-L | 420 - 99800 | 2120 | fg/mL | 80% - 120% | < 25% | 617 | fg/mL |
| Tau(total) | 17.5 - 4160 | 77.5 | fg/mL | 80% - 120% | < 25% | 16.5 | fg/mL |

Stock Concentration of Calibrator:

| Name | Item | Concentration | Units |
|------------|----------------|---------------|-------|
| GFAP | 20X Calibrator | 18540000 | fg/mL |
| NF-L | 20X Calibrator | 84800000 | fg/mL |
| Tau(total) | 20X Calibrator | 3100000 | 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 | | 05 JUL 2024 |