

| | |
|-----------------|-------------|
| Product Number: | 103520 |
| Lot Number: | 503840 |
| Expiration: | 11-Nov-2023 |
| Platform(s): | HD-X & SR-X |



Data below represents results generated on the Simoa® HD-X Analyzer.

| Component | Part Number | Lot Number |
|---------------------|-------------|------------------|
| Bead Reagent | 103510 | 231507 |
| Detector Reagent | 103512 | 307606 |
| SBG Reagent | 103514 | 232007 |
| Sample Diluent | 103516 | 232008 |
| Calibrators | 103517 | 305807 |
| Control 1 | 103518 | 305808 |
| Control 2 | 103519 | 305808 |
| SBG Diluent (SR-X)† | 103158 | N/A ¹ |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP / SBG Diluent (SR-X) are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa® Neuro 2-Plex B SR-X Kit Instructions (KI-0062).

Release Materials

NF-light®

| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|--------|----------------|---------------------|
| Control 1 | 305808 | 6.56 | 4.45 - 8.21 |
| Control 2 | 305808 | 457 | 403 - 604 |
| Panel 1 | 310007 | 6.68 | 4.21 - 9.64 |
| Panel 2 | 310007 | 4.07 | 2.12 - 6.42 |
| Panel 3 | 306904 | 87.3 | 73.9 - 111 |
| Panel 4 | 306904 | 430 | 362 - 594 |

GFAP®

| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|--------|----------------|---------------------|
| Control 1 | 305808 | 119 | 79.2 - 144 |
| Control 2 | 305808 | 10909 | 9066 - 13598 |
| Panel 1 | 310007 | 69.7 | 43.2 - 85.0 |
| Panel 2 | 310007 | 57.5 | 41.1 - 69.8 |
| Panel 3 | 306904 | 1228 | 742 - 1457 |
| Panel 4 | 306904 | 7616 | 6356 - 9535 |

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Sonia Khillan

Name

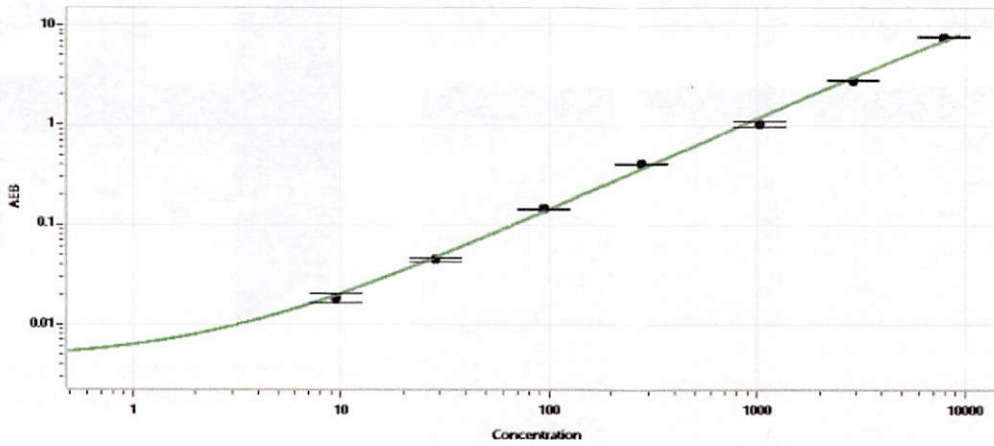
QC

Title

Sonia Khillan 01/11/2023

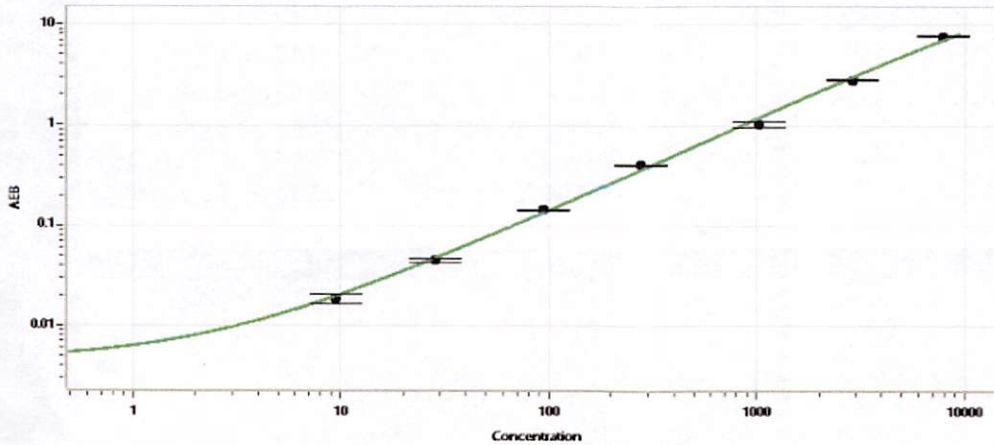
Signature/Date

NF-light® Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.552 |
| C | 1.53 |
| D | 4.81 |
| E | 12.2 |
| F | 42.7 |
| G | 122 |
| H | 387 |

GFAP® Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 9.54 |
| C | 28.5 |
| D | 93.9 |
| E | 279 |
| F | 1029 |
| G | 2898 |
| H | 7883 |