

| | |
|-----------------|-------------|
| Product Number: | 102153 |
| Lot Number: | 503063 |
| Expiration: | 22-Apr-2022 |
| Platform(s): | HD-X & SR-X |



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

| Component | Part Number | Lot Number |
|---------------------|-------------|----------------------------|
| Bead Reagent | 102141 | 120008 |
| Detector Reagent | 102143 | 120017 |
| SBG Reagent | 102145 | 120016 |
| Sample Diluent | 102147 | 120019 |
| Calibrators | 102150 | A-E: 113203 F-H: 114804 |
| Control 1 | 102151 | 113204 |
| Control 2 | 102152 | 113204 |
| SBG Diluent (SR-X)† | 103158 | N/A ¹ |
| RGP Reagent | 103159 | N/A ¹ |

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

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex A SR-X Kit Instructions (KI-0031).

Release Materials

| Tau | | | |
|-----------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 113204 | 2.70 | 1.96 - 2.94 |
| Control 2 | 113204 | 92.7 | 77.0 - 120 |
| Panel 1 | 111709 | 3.11 | 159 - 3.43 |
| Panel 2 | 111709 | 3.47 | 1.73 - 3.76 |
| Panel 3 | 111709 | 30.7 | 23.8 - 35.8 |

| NF-light® | | | |
|-----------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 113204 | 8.37 | 6.23 - 9.35 |
| Control 2 | 113204 | 747 | 596 - 894 |
| Panel 1 | 111709 | 5.64 | 4.28 - 6.41 |
| Panel 2 | 111709 | 11.0 | 8.26 - 12.4 |
| Panel 3 | 111709 | 211 | 154 - 230 |

| GFAP® | | | |
|-----------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 113204 | 33.9 | 25.6 - 38.4 |
| Control 2 | 113204 | 1177 | 975 - 1473 |
| Panel 1 | 111709 | 32.8 | 25.5 - 38.2 |
| Panel 2 | 111709 | 109 | 80.9 - 121 |
| Panel 3 | 111709 | 1716 | 1236 - 1957 |

| UCHL1® | | | |
|-----------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 113204 | 114 | 83.1 - 139 |
| Control 2 | 113204 | 10060 | 7613 - 13364 |
| Panel 1 | 111709 | 102 | 65.4 - 115 |
| Panel 2 | 111709 | 123 | 78.1 - 155 |
| Panel 3 | 111709 | 2072 | 1460 - 2190 |

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

Review/Approval

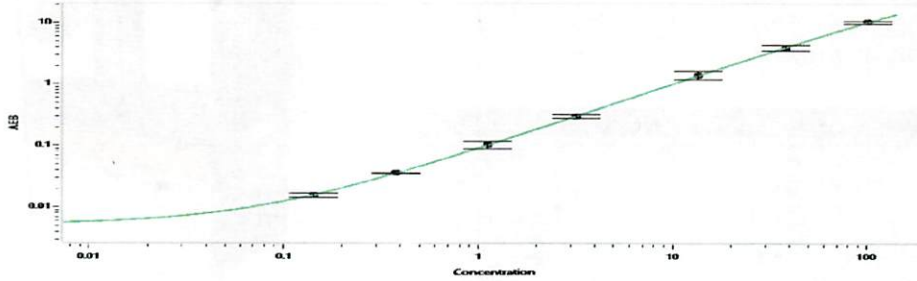
Junpil Park
Name

QC Supervisor
Title


Signature/Date

02 Aug 2021

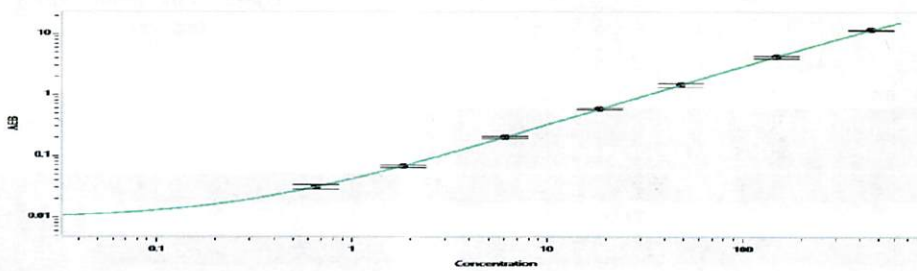
Tau Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.144 |
| C | 0.381 |
| D | 1.12 |
| E | 3.18 |
| F | 13.5 |
| G | 38.5 |
| H | 102 |

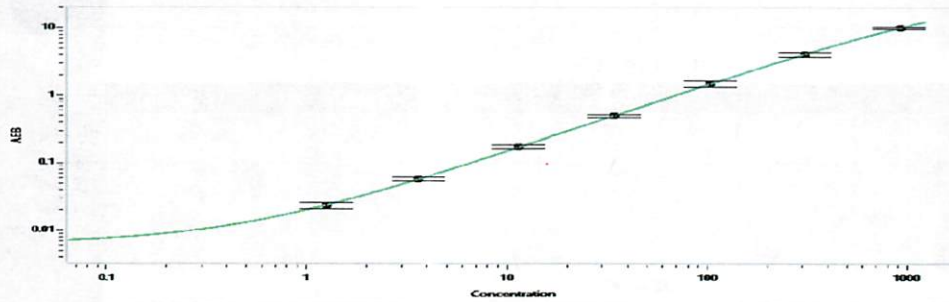
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.650 |
| C | 1.85 |
| D | 6.04 |
| E | 18.5 |
| F | 47.9 |
| G | 148 |
| H | 455 |

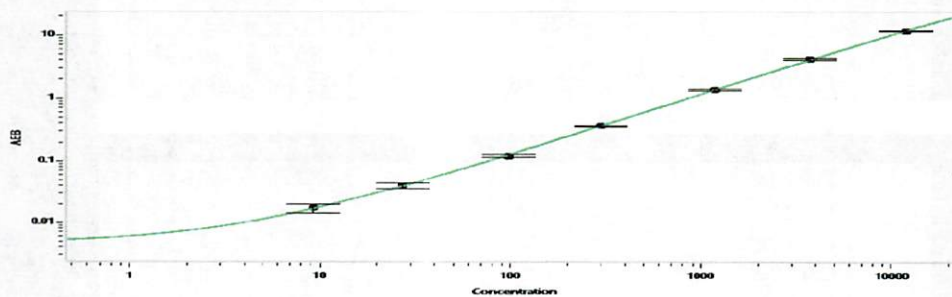
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 1.25 |
| C | 3.60 |
| D | 11.4 |
| E | 34.6 |
| F | 104 |
| G | 305 |
| H | 918 |

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 9.20 |
| C | 27.4 |
| D | 97.0 |
| E | 299 |
| F | 1186 |
| G | 3696 |
| H | 11863 |