

| | |
|-----------------|------------------|
| Product Number: | 102153 |
| Lot Number: | 502001 |
| Expiration: | 19-Feb-2020 |
| Platform(s): | HD-1/HD-X & SR-X |



Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component | Part Number | Lot Number |
|---------------------|---------------|------------------|
| Bead Reagent | 102141 | 922801 |
| Detector Reagent | 102143 | 922803 |
| SBG Reagent | 102145 | 908107 |
| Sample Diluent | 102147 | 905009 |
| Calibrators | 102150 | 920401 |
| Controls | 102151-102152 | 920402 |
| SBG Diluent (SR-X)† | 103158 | N/A ¹ |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is 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 | 920402 | 1.66 | 1.10-1.71 |
| Control 2 | 920402 | 70.6 | 57.1-85.6 |
| Panel 1 | MF-101617-01 | 1.36 | 0.605-1.92 |
| Panel 2 | MF-101617-02 | 30.5 | 22.7-44.5 |
| Panel 3 | MF-101617-03 | 128 | 96.8-194 |

NF-light®

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 920402 | 5.43 | 4.23-6.34 |
| Control 2 | 920402 | 402 | 371-556 |
| Panel 1 | MF-101617-01 | 5.10 | 3.08-6.38 |
| Panel 2 | MF-101617-02 | 36.5 | 16.7-50.4 |
| Panel 3 | MF-101617-03 | 364 | 234-541 |

GFAP®

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 920402 | 19.7 | 16.5-24.7 |
| Control 2 | 920402 | 822 | 809-1213 |
| Panel 1 | MF-101617-01 | 45.4 | 40.9-62.1 |
| Panel 2 | MF-101617-02 | 185 | 111-227 |
| Panel 3 | MF-101617-03 | 2165 | 2085-3303 |

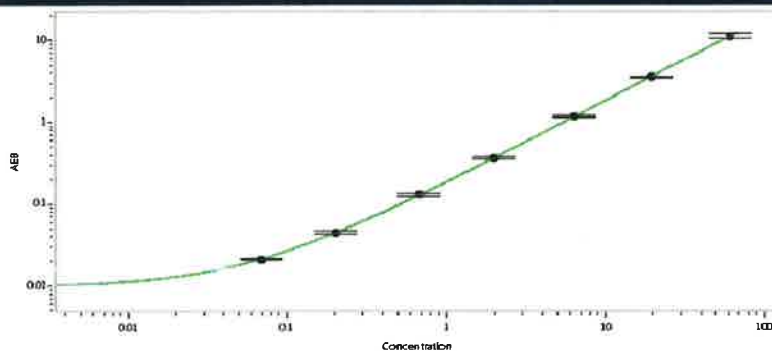
UCHL1®

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 920402 | 89.3 | 78.3-138 |
| Control 2 | 920402 | 8632 | 7808-11713 |
| Panel 1 | MF-101617-01 | 59.4 | 42.9-87.8 |
| Panel 2 | MF-101617-02 | 408 | 249-503 |
| Panel 3 | MF-101617-03 | 4673 | 4232-6348 |

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

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 31 Oct 2019
 Name Title Signature/Date

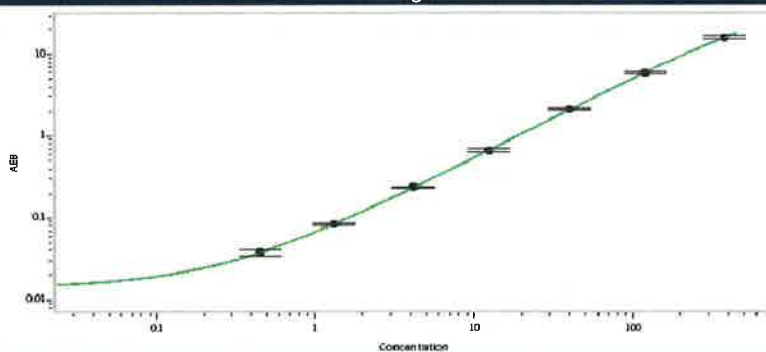
Tau Calibration Curve



**Calibrator Levels
(pg/mL)**

| | |
|---|-------|
| A | 0.000 |
| B | 0.070 |
| C | 0.204 |
| D | 0.679 |
| E | 1.99 |
| F | 6.34 |
| G | 19.7 |
| H | 61.3 |

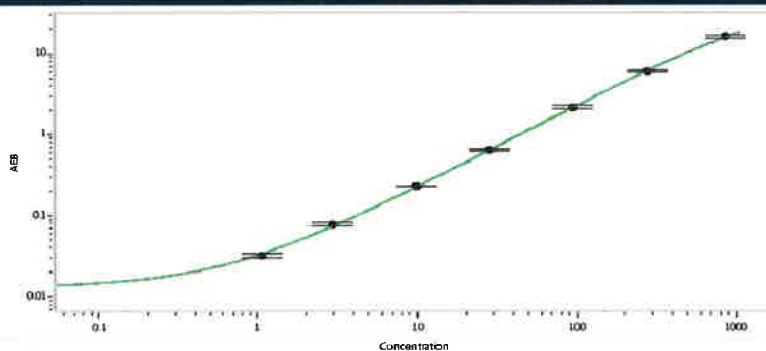
NF-light® Calibration Curve



**Calibrator Levels
(pg/mL)**

| | |
|---|-------|
| A | 0.000 |
| B | 0.455 |
| C | 1.31 |
| D | 4.12 |
| E | 12.2 |
| F | 40.0 |
| G | 120 |
| H | 376 |

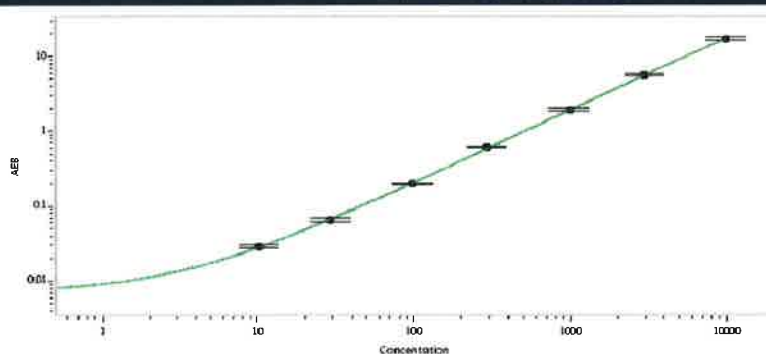
GFAP® Calibration Curve



**Calibrator Levels
(pg/mL)**

| | |
|---|-------|
| A | 0.000 |
| B | 1.06 |
| C | 2.93 |
| D | 9.77 |
| E | 28.2 |
| F | 93.6 |
| G | 277 |
| H | 855 |

UCHL1® Calibration Curve



**Calibrator Levels
(pg/mL)**

| | |
|---|-------|
| A | 0.000 |
| B | 10.1 |
| C | 28.9 |
| D | 97.8 |
| E | 288 |
| F | 981 |
| G | 2981 |
| H | 9782 |