

|                 |             |
|-----------------|-------------|
| Product Number: | 102153      |
| Lot Number:     | 501888      |
| Expiration:     | 23-Feb-2020 |
| Platform(s):    | HD-1 & SR-X |



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

| Component           | Part Number   | Lot Number       |
|---------------------|---------------|------------------|
| Bead Reagent        | 102141        | 919210           |
| Detector Reagent    | 102143        | 919215           |
| SBG Reagent         | 102145        | 905402           |
| Sample Diluent      | 102147        | 917503           |
| Calibrators         | 102150        | 912702           |
| Controls            | 102151-102152 | 912703           |
| SBG Diluent (SR-X)† | 103158        | N/A <sup>1</sup> |
| RGP Reagent         | 103159        | N/A <sup>1</sup> |

<sup>1</sup> 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).

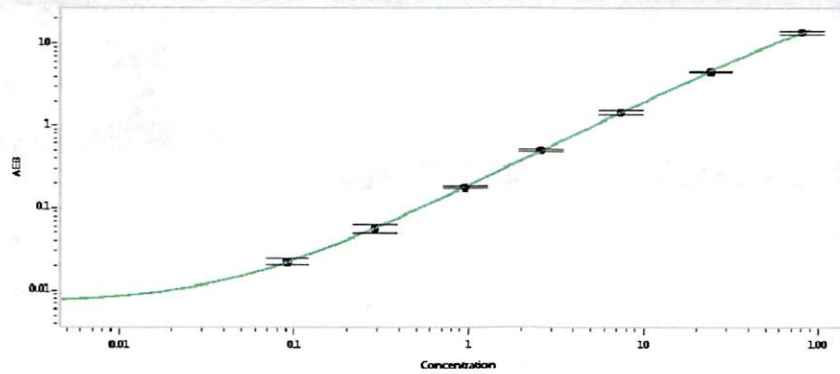
**Release Materials**

| Tau       |              |                |                    |
|-----------|--------------|----------------|--------------------|
|           | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 912703       | 1.80           | 1.37-2.47          |
| Control 2 | 912703       | 81.7           | 67.3-104           |
| Panel 1   | MF-101617-01 | 1.78           | 0.605-1.92         |
| Panel 2   | MF-101617-02 | 35.0           | 22.7-44.5          |
| Panel 3   | MF-101617-03 | 160            | 96.8-194           |
| NF-light® |              |                |                    |
|           | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 912703       | 5.33           | 4.54-6.81          |
| Control 2 | 912703       | 487            | 411-616            |
| Panel 1   | MF-101617-01 | 4.62           | 3.08-6.38          |
| Panel 2   | MF-101617-02 | 40.4           | 16.7-50.4          |
| Panel 3   | MF-101617-03 | 424            | 234-541            |
| GFAP®     |              |                |                    |
|           | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 912703       | 21.0           | 17.5-26.2          |
| Control 2 | 912703       | 1041           | 855-1282           |
| Panel 1   | MF-101617-01 | 54.3           | 40.9-62.1          |
| Panel 2   | MF-101617-02 | 190            | 111-227            |
| Panel 3   | MF-101617-03 | 2594           | 2085-3303          |
| UCHL1®    |              |                |                    |
|           | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 912703       | 123            | 93.2-161           |
| Control 2 | 912703       | 12477          | 9622-14433         |
| Panel 1   | MF-101617-01 | 51.7           | 42.9-87.8          |
| Panel 2   | MF-101617-02 | 280            | 249-503            |
| Panel 3   | MF-101617-03 | 4476           | 4232-6348          |

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

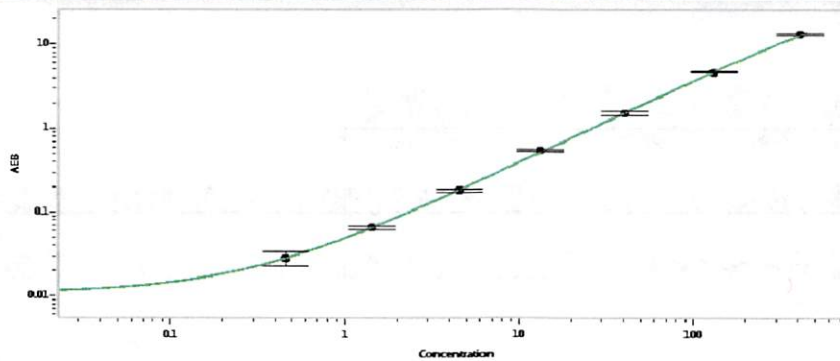
|                 |                  |                     |                   |             |
|-----------------|------------------|---------------------|-------------------|-------------|
| Review/Approval | Patrizia Stadler | Assoc. Director, QC | <i>P. Stadler</i> | 19 Jul 2019 |
|                 | <b>Name</b>      | <b>Title</b>        | <b>Signature</b>  | <b>Date</b> |

**Tau Calibration Curve**



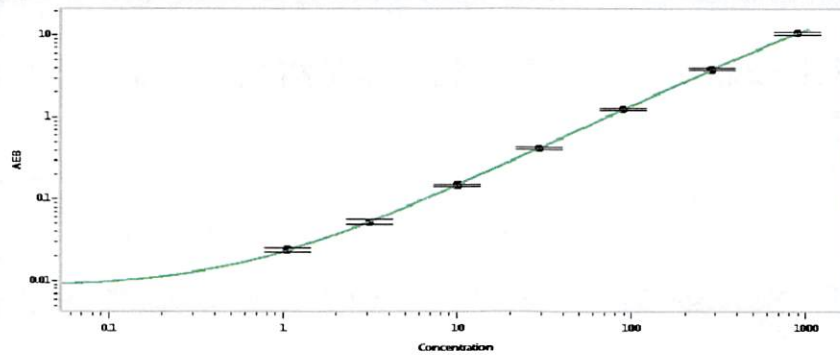
| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.092 |
| C                         | 0.290 |
| D                         | 0.952 |
| E                         | 2.56  |
| F                         | 7.43  |
| G                         | 24.2  |
| H                         | 81.4  |

**NF-light® Calibration Curve**



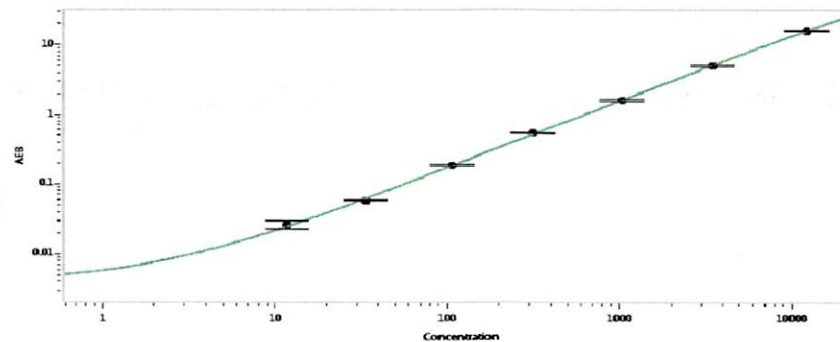
| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.457 |
| C                         | 1.44  |
| D                         | 4.54  |
| E                         | 13.1  |
| F                         | 40.8  |
| G                         | 131   |
| H                         | 412   |

**GFAP® Calibration Curve**



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 1.05  |
| C                         | 3.12  |
| D                         | 9.94  |
| E                         | 29.0  |
| F                         | 89.6  |
| G                         | 286   |
| H                         | 891   |

**UCHL1® Calibration Curve**



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 11.8  |
| C                         | 33.5  |
| D                         | 107   |
| E                         | 314   |
| F                         | 1025  |
| G                         | 3496  |
| H                         | 12235 |