

|                 |                  |
|-----------------|------------------|
| Product Number: | 103345           |
| Lot Number:     | 503684           |
| Expiration:     | 9-Nov-2023       |
| Platform(s):    | HD-1/HD-X & SR-X |

| Component           | Part Number | Lot Number       |
|---------------------|-------------|------------------|
| Bead Reagent        | 103338      | 231311           |
| Detector Reagent    | 103339      | 235310           |
| SBG Reagent         | 103340      | 231505           |
| Sample Diluent      | 103341      | 231506           |
| Calibrators         | 103342      | 225803           |
| Control 1           | 103343      | 217802           |
| Control 2           | 103344      | 217802           |
| 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 B SR-X Kit Instructions (KI-0123).



Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

#### Release Materials

##### Tau

|           | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|--------|----------------|---------------------|
| Control 1 | 217802 | 2.40           | 1.62-2.77           |
| Control 2 | 217802 | 81.8           | 62.4-107            |
| Panel 3   | 206807 | 27.5           | 20.2-30.3           |
| Panel 4   | 206807 | 108            | 79.2-119            |

##### NF-light®

|           | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|--------|----------------|---------------------|
| Control 1 | 217802 | 4.91           | 3.69-5.69           |
| Control 2 | 217802 | 440            | 311-572             |
| Panel 3   | 206807 | 169            | 130-195             |
| Panel 4   | 206807 | 788            | 617-926             |

##### GFAP®

|           | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|--------|----------------|---------------------|
| Control 1 | 217802 | 135            | 90.6-166            |
| Control 2 | 217802 | 10054          | 7164-12142          |
| Panel 3   | 206807 | 995            | 655-1090            |
| Panel 4   | 206807 | 4963           | 3366-5512           |

##### UCH-L1®

|           | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|--------|----------------|---------------------|
| Control 1 | 217802 | 115            | 84.4-135            |
| Control 2 | 217802 | 10139          | 7961-11942          |
| Panel 3   | 206807 | 1214           | 863-1443            |
| Panel 4   | 206807 | 8188           | 6163-9265           |

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

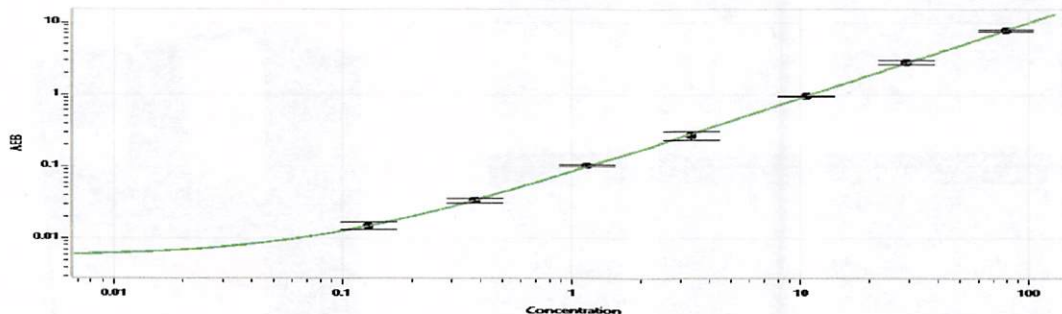
Review/Approval

*Sonia Khilku*  
Name

*QC Manager*  
Title

*Sonia Khilku / 24 FEB 2023*  
Signature/Date

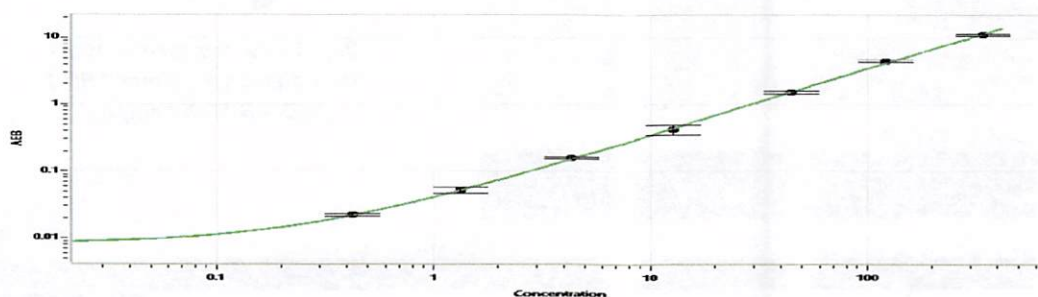
**Tau Calibration Curve**



**Calibrator Levels  
(pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.131 |
| C | 0.375 |
| D | 1.17  |
| E | 3.35  |
| F | 10.6  |
| G | 29.1  |
| H | 78.5  |

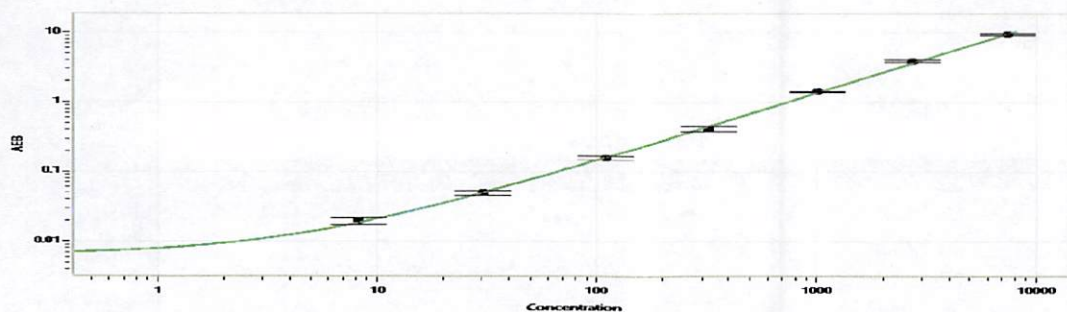
**NF-light® Calibration Curve**



**Calibrator Levels  
(pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.427 |
| C | 1.33  |
| D | 4.36  |
| E | 12.7  |
| F | 44.9  |
| G | 122   |
| H | 340   |

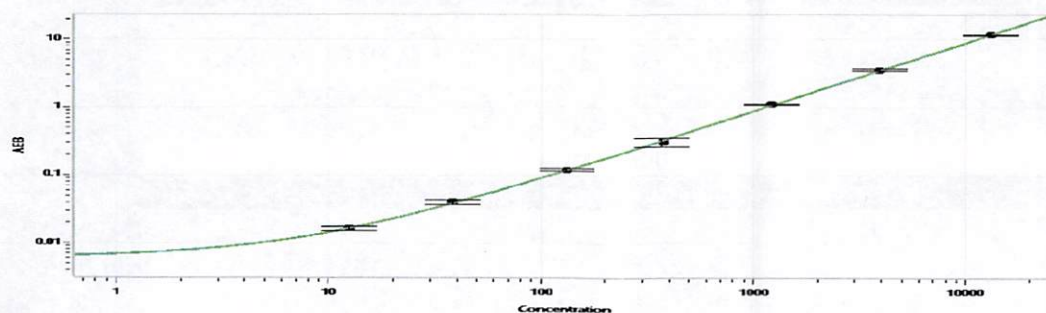
**GFAP® Calibration Curve**



**Calibrator Levels  
(pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 8.22  |
| C | 30.3  |
| D | 109   |
| E | 320   |
| F | 1012  |
| G | 2741  |
| H | 7457  |

**UCHL1® Calibration Curve**



**Calibrator Levels  
(pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 12.5  |
| C | 38.4  |
| D | 132   |
| E | 373   |
| F | 1223  |
| G | 3926  |
| H | 13117 |