

| Neuro 2-Plex B Kit |             |
|--------------------|-------------|
| Product Number:    | 103520      |
| Lot Number:        | 503416      |
| Expiration:        | 26-Jan-2023 |
| Platform(s):       | HD-X & SR-X |



| Component           | Part Number | Lot Number       |
|---------------------|-------------|------------------|
| Bead Reagent        | 103510      | 204906           |
| Detector Reagent    | 103512      | 204907           |
| SBG Reagent         | 103514      | 205001           |
| Sample Diluent      | 103516      | 205002           |
| Calibrators         | 103517      | 203302           |
| Control 1           | 103518      | 203303           |
| Control 2           | 103519      | 203303           |
| SBG Diluent (SR-X)† | 103158      | N/A <sup>1</sup> |
| RGP Reagent         | 103159      | N/A <sup>1</sup> |

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

<sup>1</sup> 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) | Mean Range (pg/mL) |
|-----------|--------|----------------|--------------------|
| Control 1 | 203303 | 5.86           | 4.53-7.10          |
| Control 2 | 203303 | 536            | 435-653            |
| Panel 1   | 207003 | 14.2           | 10.7-16.1          |
| Panel 2   | 207003 | 6.44           | 4.77-7.68          |
| Panel 3   | 207003 | 126            | 101-151            |
| Panel 4   | 207003 | 321            | 273-416            |

**GFAP®**

|           | Lot    | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------|----------------|--------------------|
| Control 1 | 203303 | 91.4           | 72.2-111           |
| Control 2 | 203303 | 9231           | 7332-10997         |
| Panel 1   | 207003 | 57.1           | 49.4-83.9          |
| Panel 2   | 207003 | 63.9           | 52.9-104           |
| Panel 3   | 207003 | 757            | 684-1104           |
| Panel 4   | 207003 | 4846           | 4259-7215          |

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

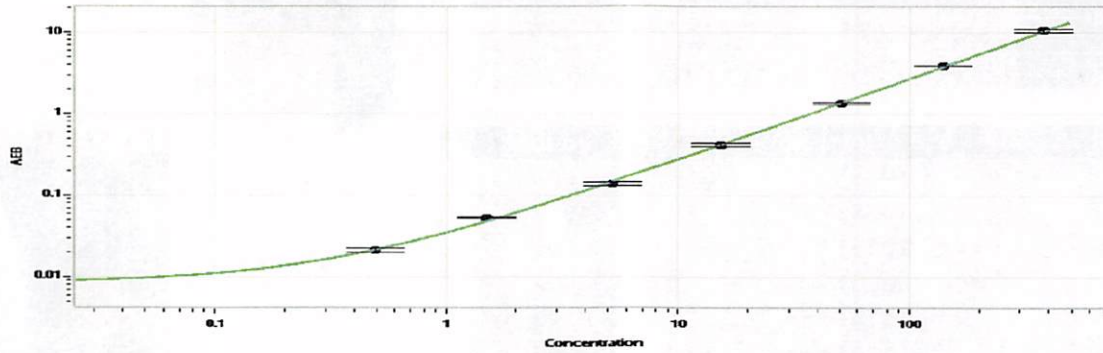
Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 29 Mar 2022  
Signature/Date

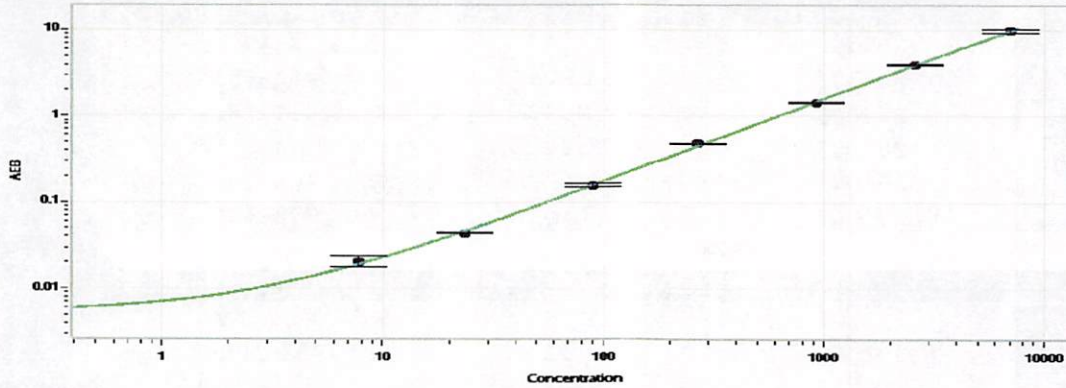
**NF-light® Calibration Curve**



**Calibrator Levels  
(pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.492 |
| C | 1.50  |
| D | 5.20  |
| E | 15.3  |
| F | 50.7  |
| G | 140   |
| H | 377   |

**GFAP® Calibration Curve**



**Calibrator Levels  
(pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 7.75  |
| C | 23.6  |
| D | 89.5  |
| E | 268   |
| F | 925   |
| G | 2601  |
| H | 7092  |