

|                        |               |
|------------------------|---------------|
| <b>Product Number:</b> | 104670        |
| <b>Lot Number:</b>     | 504607        |
| <b>Expiration:</b>     | 19-Nov-2026   |
| <b>Platform(s):</b>    | HD-X and SR-X |

| Component                 | Part Number | Lot Number       | Control Lot Number |
|---------------------------|-------------|------------------|--------------------|
| Bead Reagent              | 104652      | 513905           | N/A                |
| Detector Reagent          | 104654      | 514908           | N/A                |
| SBG Reagent               | 104656      | 514909           | N/A                |
| Plasma Sample Diluent     | 104722      | 514706           | N/A                |
| CSF Sample Diluent        | 104723      | 514309           | N/A                |
| Calibrator Diluent        | 104657      | 514310           | N/A                |
| Control Diluent           | 104721      | 514311           | N/A                |
| Calibrator Concentrate    | 104666      | 317341           | N/A                |
| Control 1 Concentrate     | 104668      | 317342           | N/A                |
| Control 2 Concentrate     | 104669      | 317343           | 317343A            |
| RGP 2.0 Reagent           | 104321      | N/A <sup>1</sup> | N/A <sup>1</sup>   |
| Activation Buffer Reagent | 104324      | N/A <sup>1</sup> | N/A <sup>1</sup>   |

<sup>1</sup> Reagents are not Kit Lot Specific

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



Review/Approval

Dharmendra Patel      DOC Control II, QA      [Signature] / 24 Jun 2025

Name                      Title                      Signature/Date

**Release Materials\***

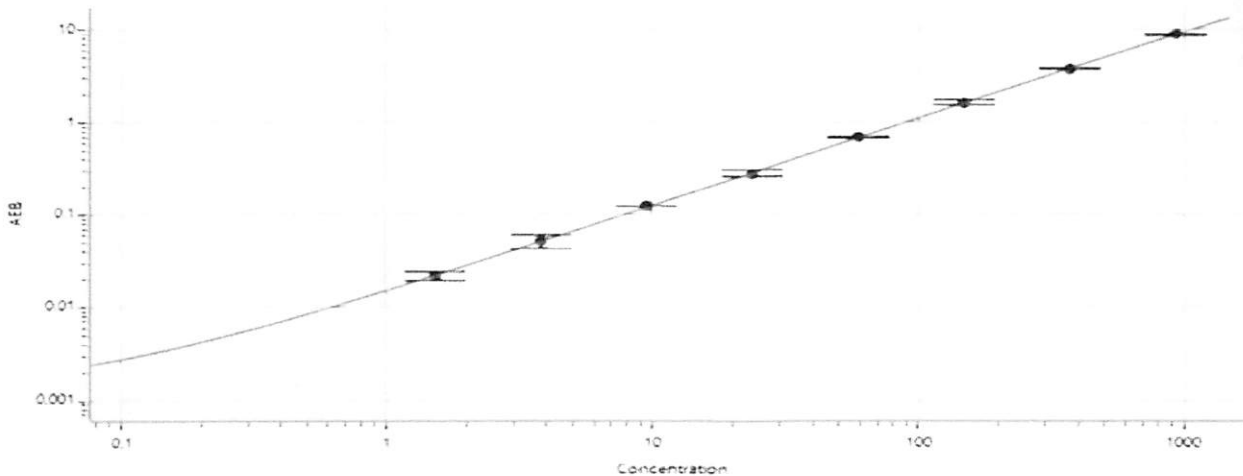
**GFAP®**

|           | Lot     | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|---------|----------------|---------------------|
| Control 1 | 317342  | 116            | 84.9 - 158          |
| Control 2 | 317343A | 1639           | 1080 - 2005         |

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

|                                                   |      |
|---------------------------------------------------|------|
| Calibrator Concentrate Reconstitution Volume (uL) | 1144 |
| Control 1 Concentrate Reconstitution Volume (uL)  | 1144 |
| Control 2 Concentrate Reconstitution Volume (uL)  | 1144 |

**Calibration Curve**



**Calibrator Levels (pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 1.52  |
| C | 3.80  |
| D | 9.48  |
| E | 23.8  |
| F | 59.5  |
| G | 149   |
| H | 372   |
| I | 929   |

**Release Materials\***

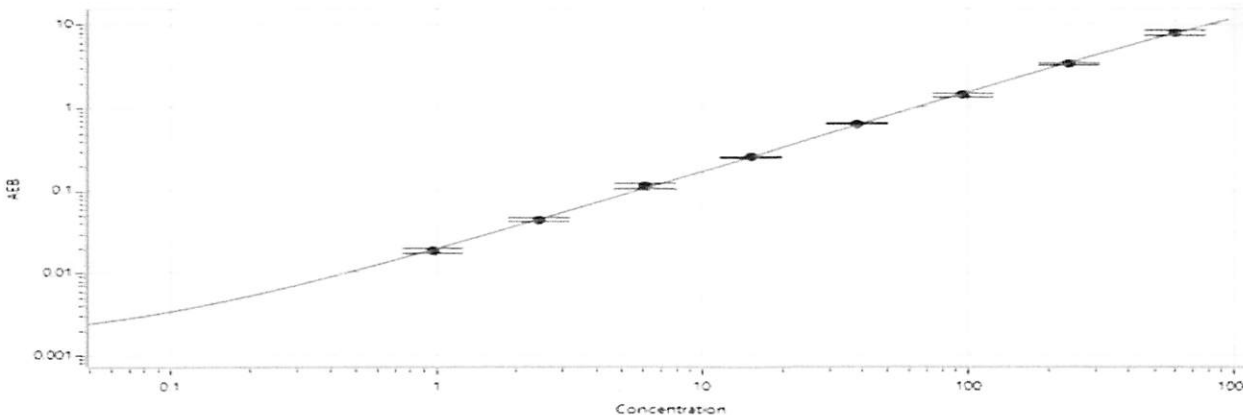
**NF-Light®**

|           | Lot     | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|---------|----------------|---------------------|
| Control 1 | 317342  | 89.8           | 62.6 - 116          |
| Control 2 | 317343A | 1135           | 808 - 1500          |

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

|                                                   |      |
|---------------------------------------------------|------|
| Calibrator Concentrate Reconstitution Volume (uL) | 1144 |
| Control 1 Concentrate Reconstitution Volume (uL)  | 1144 |
| Control 2 Concentrate Reconstitution Volume (uL)  | 1144 |

**Calibration Curve**



**Calibrator Levels (pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.975 |
| C | 2.44  |
| D | 6.09  |
| E | 15.3  |
| F | 38.2  |
| G | 95.2  |
| H | 238   |
| I | 596   |