

|                 |               |
|-----------------|---------------|
| Product Number: | 104670        |
| Lot Number:     | 504575        |
| Expiration:     | 10-Sep-2026   |
| Platform(s):    | HD-X and SR-X |

| Component                 | Part Number | Lot Number       |
|---------------------------|-------------|------------------|
| Bead Reagent              | 104652      | 509013           |
| Detector Reagent          | 104654      | 509301           |
| SBG Reagent               | 104656      | 509302           |
| Plasma Sample Diluent     | 104722      | 509019           |
| CSF Sample Diluent        | 104723      | 509018           |
| Calibrator Diluent        | 104657      | 506918           |
| Control Diluent           | 104721      | 506919           |
| Calibrator Concentrate    | 104666      | 402948           |
| Control 1 Concentrate     | 104668      | 402949           |
| Control 2 Concentrate     | 104669      | 402950           |
| RGP 2.0 Reagent           | 104321      | N/A <sup>1</sup> |
| Activation Buffer Reagent | 104324      | 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

*Dhruvi Patel*

Name

*Doc Control II, QA*

Title

*SDP 22 May 2025*

Signature/Date

**Release Materials\***

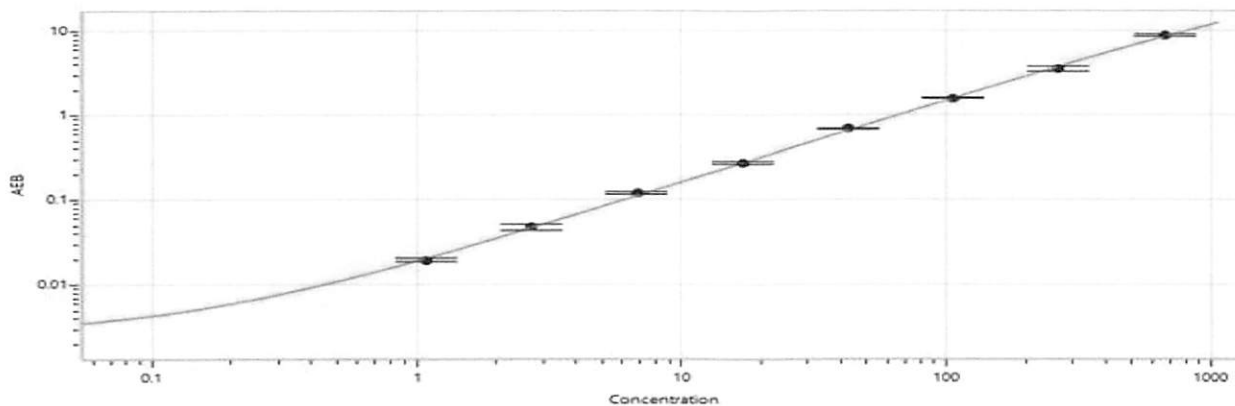
**GFAP®**

|           | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|--------|----------------|---------------------|
| Control 1 | 402949 | 81.6           | 62.5 - 116          |
| Control 2 | 402950 | 1167           | 809 - 1503          |

\* 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 | 1.09  |
| C | 2.73  |
| D | 6.81  |
| E | 17.1  |
| F | 42.6  |
| G | 106   |
| H | 266   |
| I | 666   |

**Release Materials\***

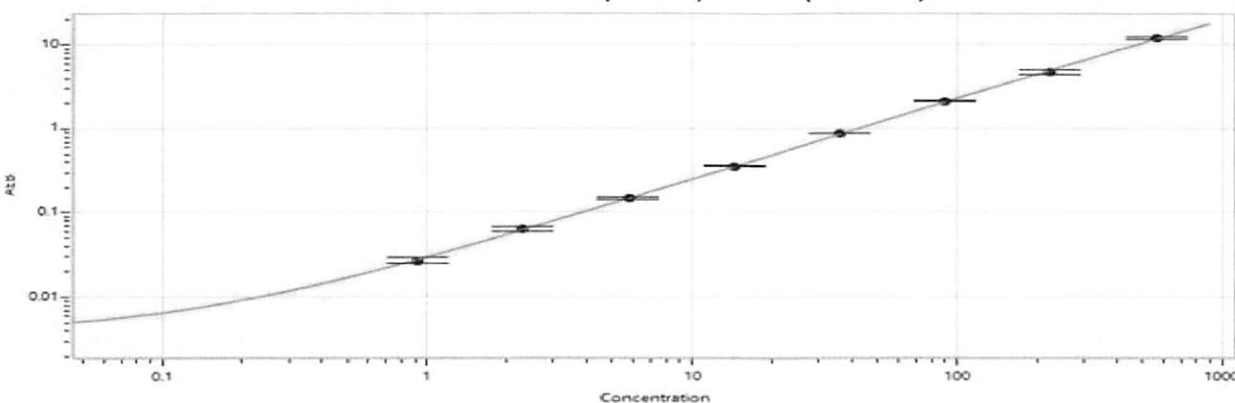
**NF-Light®**

|           | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
|-----------|--------|----------------|---------------------|
| Control 1 | 402949 | 75.1           | 58.3 - 108          |
| Control 2 | 402950 | 1020           | 753 - 1398          |

\* 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 | 0.921 |
| C | 2.30  |
| D | 5.75  |
| E | 14.4  |
| F | 35.9  |
| G | 89.9  |
| H | 224   |
| I | 562   |