

|                    |                    |
|--------------------|--------------------|
| Product Number:    | 103671             |
| Lot Number:        | 503972             |
| Paired Reagent Kit | 503893             |
| Expiration:        | 20-Apr-2024        |
| Platform(s):       | HD-1 / HD-X & SR-X |



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

| Component                       | Part Number | Lot Number | Control Number |
|---------------------------------|-------------|------------|----------------|
| N4PE Plasma Sample Diluent      | 103659      | 318704     | 318704A        |
| N4PE CSF Sample Diluent         | 103727      | 323502     | N/A            |
| Calibrator Diluent              | 103661      | 319917     | 319917A        |
| Abeta 40 Calibrator Concentrate | 101734      | 310203     | 310203A        |
| Abeta 42 Calibrator Concentrate | 103217      | 211005     | N/A            |
| GFAP Calibrator Concentrate     | 103622      | 219910     | N/A            |
| NF-light Calibrator Concentrate | 102619      | 329214     | N/A            |

| Calibrator Concentrate<br>Value Assignment (pg/mL) |       |
|----------------------------------------------------|-------|
| Abeta 40                                           | 2478  |
| Abeta 42                                           | 2125  |
| GFAP                                               | 82774 |
| NF-Light                                           | 15619 |

Review/Approval

*Sonia Khilgar*

Name

QC

Title

*Sonia Khilgar 26 OCT 2023*

Signature/Date