

|                    |                    |
|--------------------|--------------------|
| Product Number:    | 103671             |
| Lot Number:        | 502559             |
| Paired Reagent Kit | 502542             |
| Expiration:        | 21-Sep-2021        |
| Platform(s) :      | HD-1 / HD-X & SR-X |

| Component                       | Part Number | Lot Number |
|---------------------------------|-------------|------------|
| N4PE Plasma Sample Diluent      | 103659      | 026708     |
| N4PE CSF Sample Diluent         | 103727      | 026802     |
| Calibrator Diluent              | 103661      | 026507     |
| Abeta 40 Calibrator Concentrate | 101734      | 927002     |
| Abeta 42 Calibrator Concentrate | 103217      | 026801     |
| GFAP Calibrator Concentrate     | 103622      | 027001     |
| NF-light Calibrator Concentrate | 102619      | 026716     |

| Calibrator Concentrate<br>Value Assignment (pg/mL) |        |
|----------------------------------------------------|--------|
| Abeta 40                                           | 3592   |
| Abeta 42                                           | 2184   |
| GFAP                                               | 149367 |
| NF-Light                                           | 9343   |

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 16 Oct 2020  
Signature/Date