

| | |
|-----------------|------------|
| Product Number: | 103670 |
| Lot Number: | 503819 |
| Expiration: | 3-Mar-2024 |
| Platform(s): | HD-X |



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

| Component | Part Number | Lot Number |
|-----------------------|-------------|------------------|
| Bead Reagent | 103653 | 306209 |
| Detector Reagent | 103655 | 306205 |
| SBG Reagent | 103657 | 306206 |
| Plasma Sample Diluent | 103659 | 306207 |
| CSF Sample Diluent | 103727 | 306208 |
| Calibrators | 103667 | 306505 |
| Control 1 | 103668 | 306506 |
| Control 2 | 103669 | 306506 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| Release Materials | | | |
|-------------------|--------|----------------|---------------------|
| Abeta 40 | | | |
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 306506 | 13.9 | 10.9-16.4 |
| Control 2 | 306506 | 108 | 83.4-125 |
| Panel 1 | 226304 | 22.3 | 20.8-31.2 |
| Panel 2 | 225507 | 53.6 | 44.5 - 66.8 |
| Panel 3 | 226908 | 221 | 183 - 312 |
| Abeta 42 | | | |
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 306506 | 4.71 | 3.58-5.38 |
| Control 2 | 306506 | 26.5 | 20.9 - 31.4 |
| Panel 1 | 226304 | 2.20 | 1.65 - 2.97 |
| Panel 2 | 225507 | 20.2 | 16.5 - 24.7 |
| Panel 3 | 226908 | 48.6 | 41.2 - 61.9 |
| GFAP® | | | |
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 306506 | 149 | 117-176 |
| Control 2 | 306506 | 3225 | 2454-3681 |
| Panel 1 | 226304 | 42.1 | 28.1 - 58.0 |
| Panel 2 | 225507 | 300 | 248 - 372 |
| Panel 3 | 226908 | 2648 | 2078 - 3165 |
| NF-light® | | | |
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 306506 | 19.1 | 16.2 - 24 |
| Control 2 | 306506 | 432 | 311 - 529 |
| Panel 1 | 226304 | 3.57 | 1.38 - 4.84 |
| Panel 2 | 225507 | 108 | 85.7 - 128 |
| Panel 3 | 226908 | 499 | 367 - 698 |

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

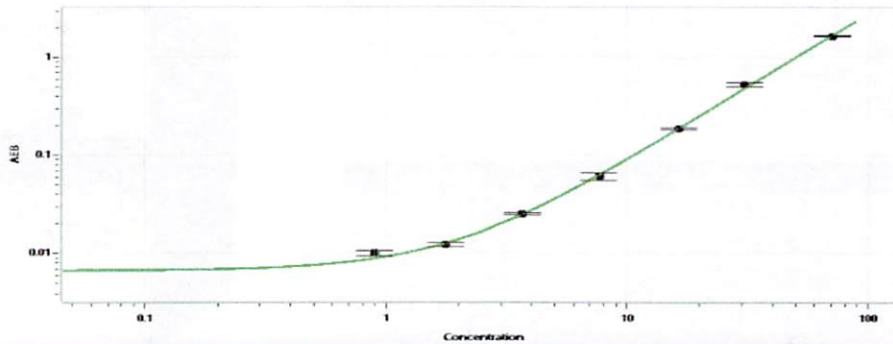
Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

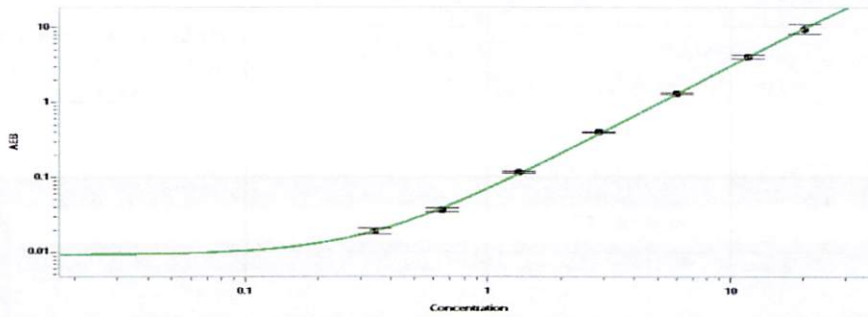
[Signature] 25 Oct 2023
Signature/Date

Abeta 40 Calibration Curve



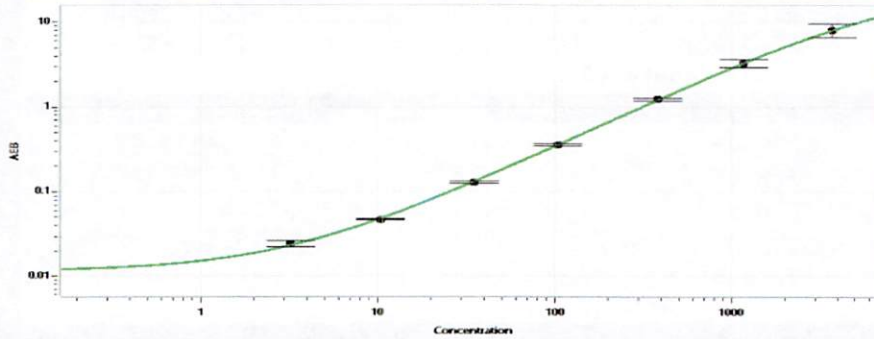
| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.897 |
| C | 1.76 |
| D | 3.67 |
| E | 7.64 |
| F | 16.3 |
| G | 30.8 |
| H | 71.3 |

Abeta 42 Calibration Curve



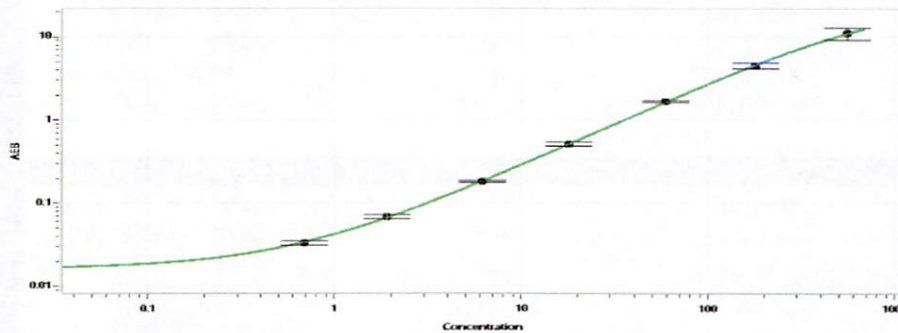
| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.343 |
| C | 0.649 |
| D | 1.34 |
| E | 2.88 |
| F | 5.96 |
| G | 11.8 |
| H | 20.2 |

GFAP® Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 3.21 |
| C | 10.4 |
| D | 34.9 |
| E | 104 |
| F | 379 |
| G | 1163 |
| H | 3709 |

NF-light® Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.691 |
| C | 1.91 |
| D | 6.17 |
| E | 17.8 |
| F | 59.2 |
| G | 180 |
| H | 552 |