

|                 |                  |
|-----------------|------------------|
| Product Number: | 101995           |
| Lot Number:     | 502473           |
| Expiration:     | 24-Mar-2021      |
| Platform(s):    | HD-1/HD-X & SR-X |

| Component          | Part Number | Lot Number       |
|--------------------|-------------|------------------|
| Bead Reagent       | 101998      | 019708           |
| Detector Reagent   | 101996      | 026912           |
| SBG Reagent        | 102000      | 019809           |
| Sample Diluent     | 102002      | 019801           |
| Calibrators        | 103221      | 027402           |
| Control 1          | 103222      | 931002           |
| Control 2          | 103223      | 931002           |
| SBG Diluent (SRX)† | 103158      | N/A <sup>1</sup> |
| RGP Reagent        | 103159      | N/A <sup>1</sup> |

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

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 3-Plex A SR-X Kit Instructions (KI-0083).



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

**Release Materials**

**Tau**

|           | Lot    | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------|----------------|--------------------|
| Control 1 | 931002 | 1.77           | 0.695-2.11         |
| Control 2 | 931002 | 68.5           | 45.4-79.6          |
| Panel 1   | 003802 | 1.63           | 0.942-2.48         |
| Panel 2   | 003802 | 137            | 104-190            |
| Panel 3   | 003802 | 1.29           | 0.520-1.64         |

**Abeta 42**

|           | Lot    | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------|----------------|--------------------|
| Control 1 | 931002 | 5.34           | 4.59-6.89          |
| Control 2 | 931002 | 132            | 107-160            |
| Panel 1   | 003802 | 4.78           | 2.54-6.08          |
| Panel 2   | 003802 | 4.44           | 2.67-5.20          |
| Panel 3   | 003802 | 111            | 88.4-138           |

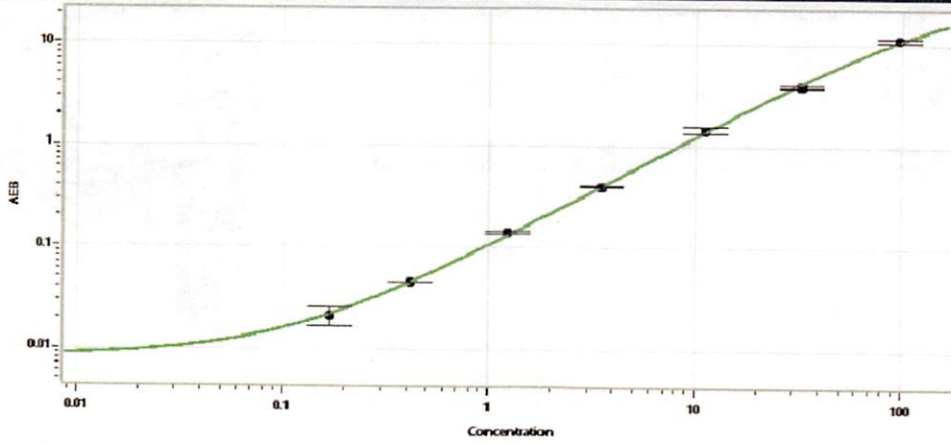
**Abeta 40**

|           | Lot    | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------|----------------|--------------------|
| Control 1 | 931002 | 17.2           | 13.3-20.0          |
| Control 2 | 931002 | 238            | 206-308            |
| Panel 1   | 003802 | 67.1           | 44.5-103           |
| Panel 2   | 003802 | 34.6           | 25.1-51.1          |
| Panel 3   | 003802 | 79.6           | 46.9-127           |

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

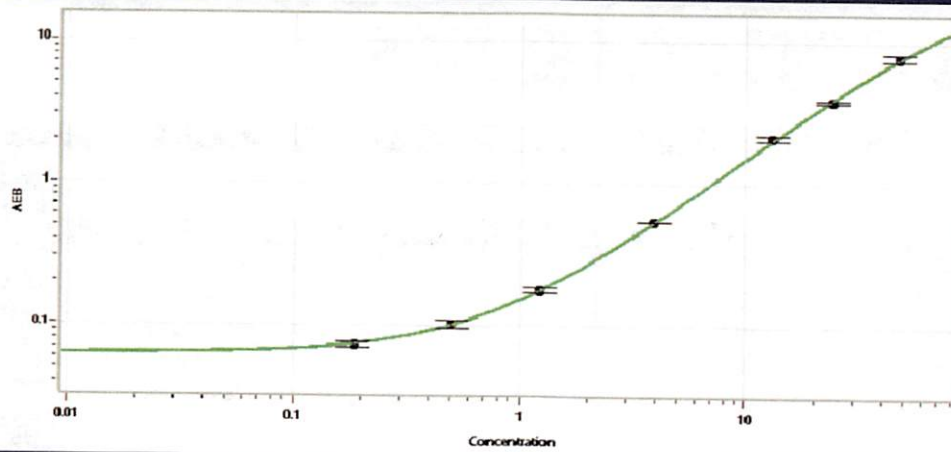
|                 |              |            |                |              |
|-----------------|--------------|------------|----------------|--------------|
| Review/Approval | Shivani Goel | QC Manager | Shivani Goel   | 15 Oct 2020. |
|                 | Name         | Title      | Signature/Date |              |

**Tau Calibration Curve**



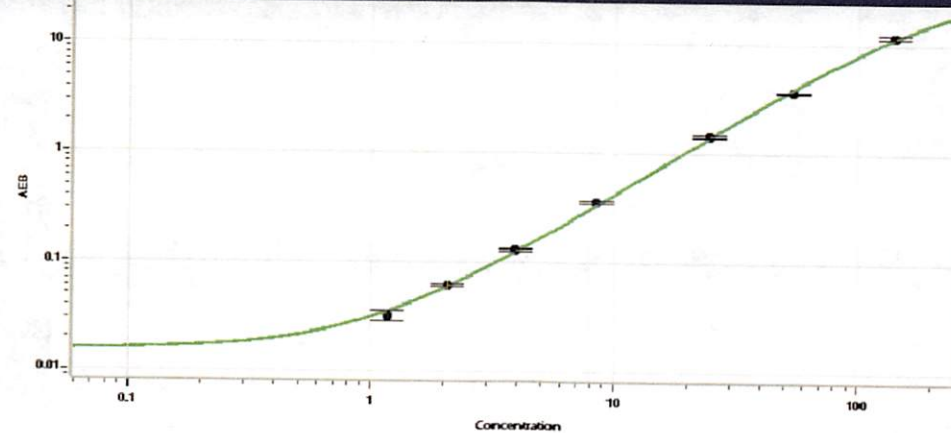
| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.170 |
| C                         | 0.413 |
| D                         | 1.24  |
| E                         | 3.52  |
| F                         | 11.2  |
| G                         | 32.7  |
| H                         | 98.2  |

**Abeta 42**



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.186 |
| C                         | 0.501 |
| D                         | 1.23  |
| E                         | 3.91  |
| F                         | 13.0  |
| G                         | 24.0  |
| H                         | 47.3  |

**Abeta 40**



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 1.17  |
| C                         | 2.08  |
| D                         | 3.93  |
| E                         | 8.36  |
| F                         | 24.5  |
| G                         | 54.0  |
| H                         | 141   |