

<b>Product Number:</b>	101995
<b>Lot Number:</b>	504136
<b>Expiration:</b>	23-Aug-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101998	405407
Detector Reagent	101996	405508
SBG Reagent	102000	408803
Sample Diluent	102002	405406
Calibrators	103221	412804
Control 1	103222	405802
Control 2	103223	405802
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

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

† Specific to use on the SRX platform - refer to Neuro-3 plex A Kit Instructions SR-X (KI-0083).



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

Review/Approval

Cody Carlson  
Name

QC Manager  
Title

Cody Carlson 22 MAY 2024  
Signature/Date

**Release Materials\***

**Tau**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Control 1</b>	405802	2.20	1.87 - 2.81
<b>Control 2</b>	405802	109	92.6 - 139
<b>Panel 1</b>	319110	3.04	2.70 - 4.05
<b>Panel 2</b>	208403	134	110 - 167
<b>Panel 3</b>	208403	1.86	1.61 - 2.84

**Abeta 42**

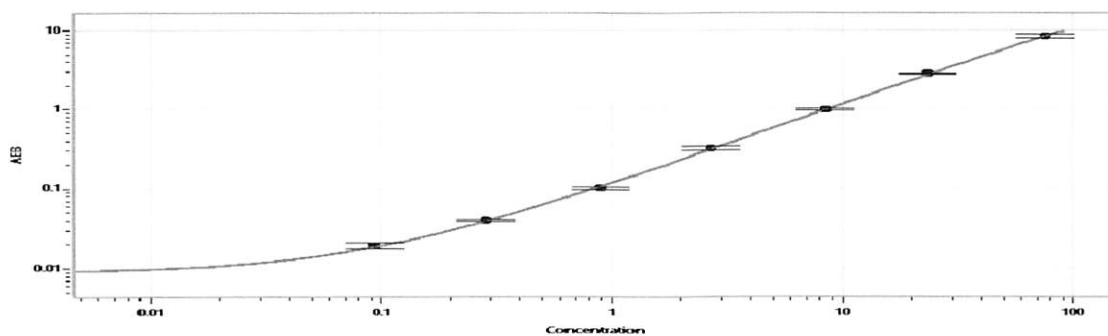
	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Control 1</b>	405802	2.87	2.63 - 3.95
<b>Control 2</b>	405802	85.4	70.8 - 106
<b>Panel 1</b>	319110	2.63	2.02 - 3.44
<b>Panel 2</b>	208403	8.74	5.69 - 11.3
<b>Panel 3</b>	208403	116	80.4 - 155

**Abeta 40**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Control 1</b>	405802	8.74	6.70 - 10.1
<b>Control 2</b>	405802	169	140 - 210
<b>Panel 1</b>	319110	85.9	58.4 - 91.4
<b>Panel 2</b>	208403	144	105 - 185
<b>Panel 3</b>	208403	90.7	60.6 - 113

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

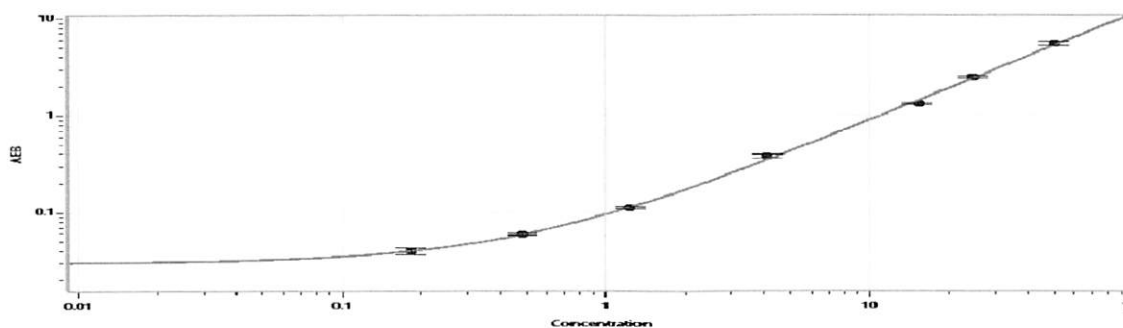
**Tau Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.093
C	0.285
D	0.902
E	2.69
F	8.50
G	23.5
H	76.0

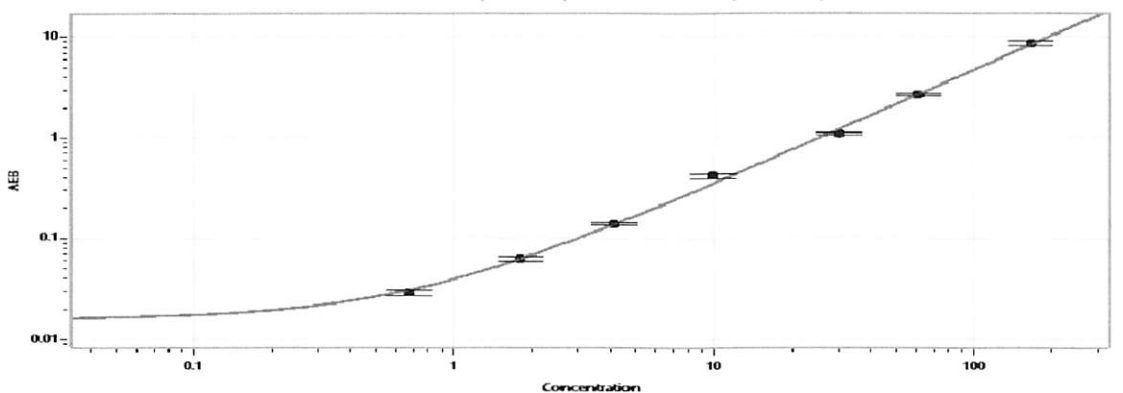
**Abeta 42 Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.182
C	0.481
D	1.23
E	4.10
F	15.3
G	24.8
H	50.2

**Abeta 40 Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.671
C	1.80
D	4.11
E	9.85
F	30.1
G	61.0
H	167