

<b>Product Number:</b>	101995
<b>Lot Number:</b>	504224
<b>Expiration:</b>	27-Dec-2024
<b>Platform(s):</b>	SR-X and HD-1/HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	101998	317809	317809A
Detector Reagent	101996	318102	318102A
SBG Reagent	102000	318101	318101A
Sample Diluent	102002	317904	317904A
Calibrators	103221	317710	317710A
Control 1	103222	315103	315103A
Control 2	103223	315103	315103A
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>	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).

Review/Approval

Cody Carlson

QC Manager

Name

Title

 18 APR 2024  
Signature/Date

**Release Materials\***

**Tau**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315103	1.18	1.02 - 1.54
Control 2	315103	52.5	39.6 - 60.9
Panel 1	319110	3.03	2.74 - 4.11
Panel 2	208403	143	109 - 169
Panel 3	208403	1.94	1.57 - 2.82

**Abeta 42**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315103	2.36	1.67 - 3.23
Control 2	315103	57.4	39.5 - 76.7
Panel 1	319110	2.46	2.28 - 3.43
Panel 2	208403	8.38	5.86 - 10.6
Panel 3	208403	113	79.2 - 150

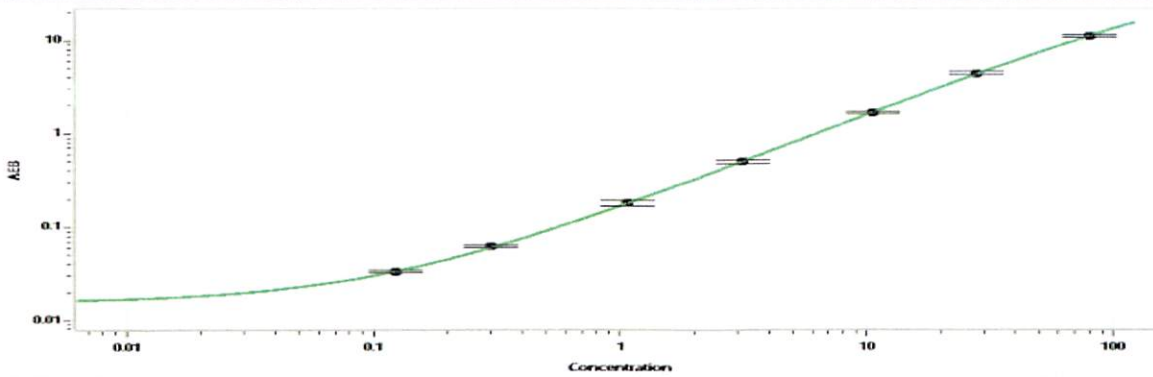
**Abeta 40**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315103	7.84	5.88 - 9.43
Control 2	315103	127	90.2 - 175
Panel 1	319110	73.7	63.4 - 95.2
Panel 2	208403	138	109 - 187
Panel 3	208403	76.1	62.0 - 115

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



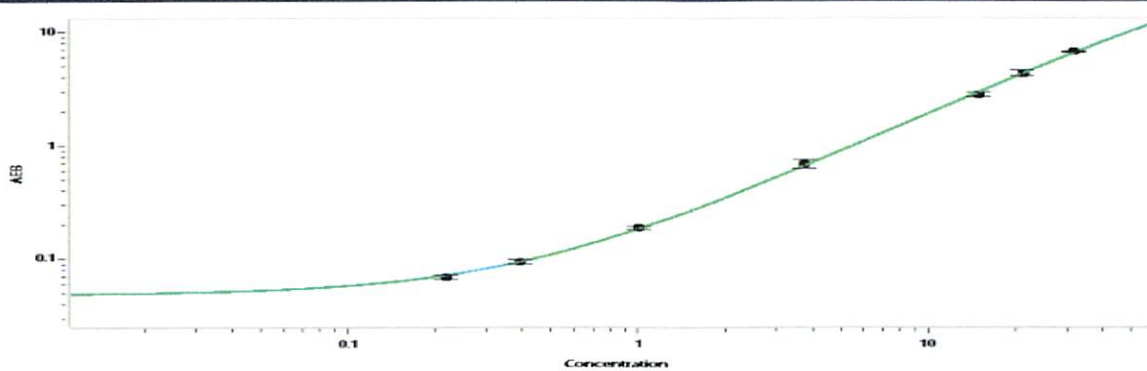
**Tau Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.123
C	0.300
D	1.07
E	3.16
F	10.5
G	27.8
H	80.0

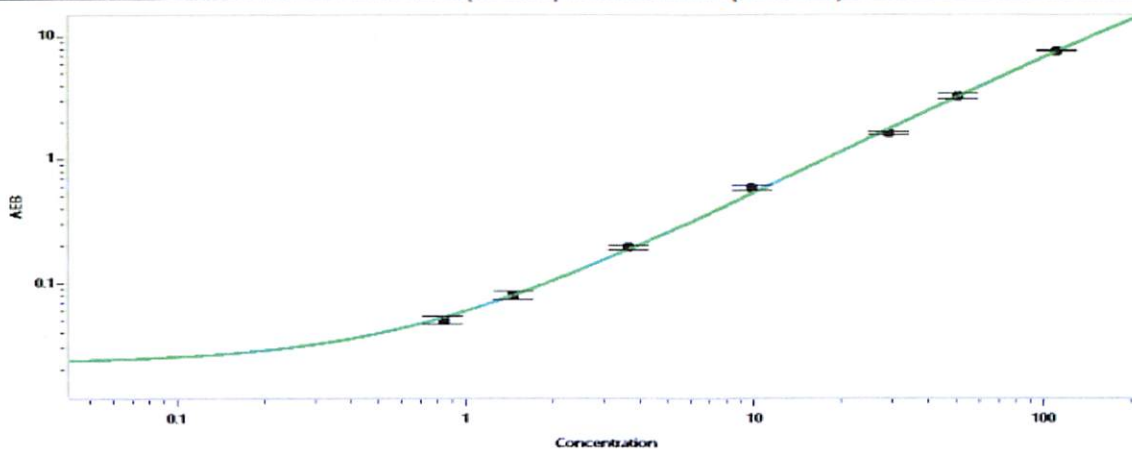
**Abeta 42 Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.219
C	0.390
D	1.01
E	3.75
F	14.9
G	21.2
H	31.9

**Abeta 40 Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.835
C	1.45
D	3.67
E	9.75
F	29.2
G	50.6
H	111