

<b>Product Number:</b>	101995
<b>Lot Number:</b>	504191
<b>Expiration:</b>	1-Jun-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101998	333604
Detector Reagent	101996	333507
SBG Reagent	102000	333508
Sample Diluent	102002	333511
Calibrators	103221	333408
Control 1	103222	333409
Control 2	103223	333409
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 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

 Name	QC MANAGER Title	 Signature/Date	28ME2024
---	---------------------	---	----------

**Release Materials\***

**Tau**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	333409	1.47	1.09 - 1.64
Control 2	333409	63.6	48.7 - 73.1
Panel 1	319110A	3.43	2.74 - 4.11
Panel 2	208403	138	110 - 167
Panel 3	208403	2.36	1.61 - 2.84

**Abeta 42**

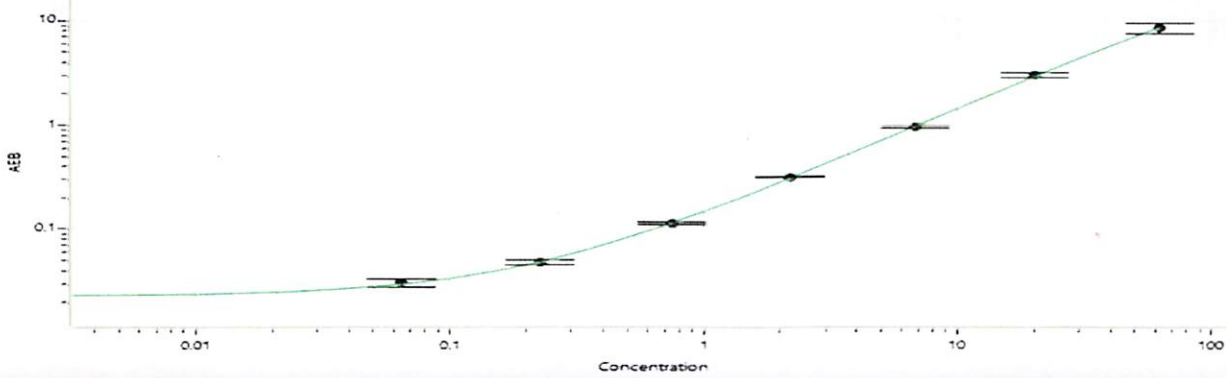
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	333409	4.14	2.42 - 4.68
Control 2	333409	97.6	68.5 - 112
Panel 1	319110A	2.92	2.30 - 3.45
Panel 2	208403	9.98	5.69 - 11.3
Panel 3	208403	132	80.4 - 155

**Abeta 40**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	333409	11.9	7.38 - 13.1
Control 2	333409	172	142 - 213
Panel 1	319110A	78.9	63.4 - 95.0
Panel 2	208403	130	105 - 185
Panel 3	208403	79.0	60.6 - 113

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

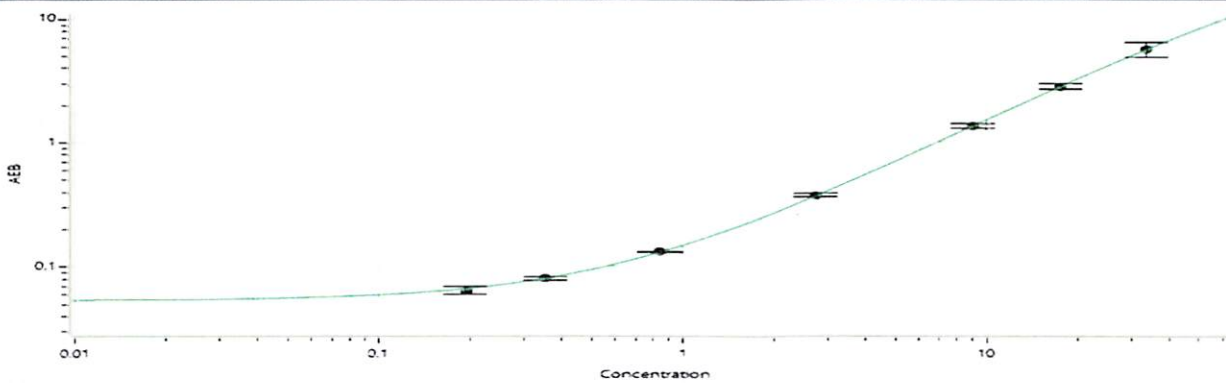
**Tau Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.064
C	0.227
D	0.757
E	2.19
F	6.79
G	20.2
H	61.9

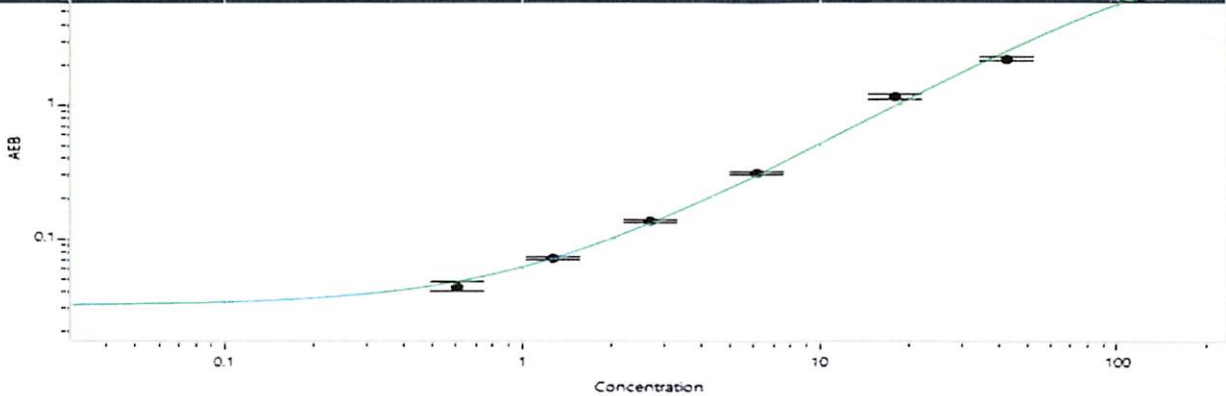
**Abeta 42 Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.193
C	0.353
D	0.841
E	2.73
F	9.01
G	17.6
H	33.7

**Abeta 40 Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.604
C	1.27
D	2.70
E	6.15
F	18.0
G	42.6
H	117