

Product Number:	101995
Lot Number:	503203
Expiration:	22-Aug-2022
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101998	129125
Detector Reagent	101996	128904
SBG Reagent	102000	128905
Sample Diluent	102002	123412
Calibrators	103221	129007
Control 1	103222	129008
Control 2	103223	129008
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	129008	3.05	2.36 - 3.54
Control 2	129008	122	97.2 - 146
Panel 1	124607	3.24	1.28 - 4.59
Panel 2	124607	225	179 - 268
Panel 3	124607	4.67	2.47 - 6.40


**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	129008	6.30	4.94 - 7.40
Control 2	129008	160	132 - 201
Panel 1	124607	5.80	4.58 - 8.08
Panel 2	124607	5.35	4.77 - 7.19
Panel 3	124607	104	84.8 - 127

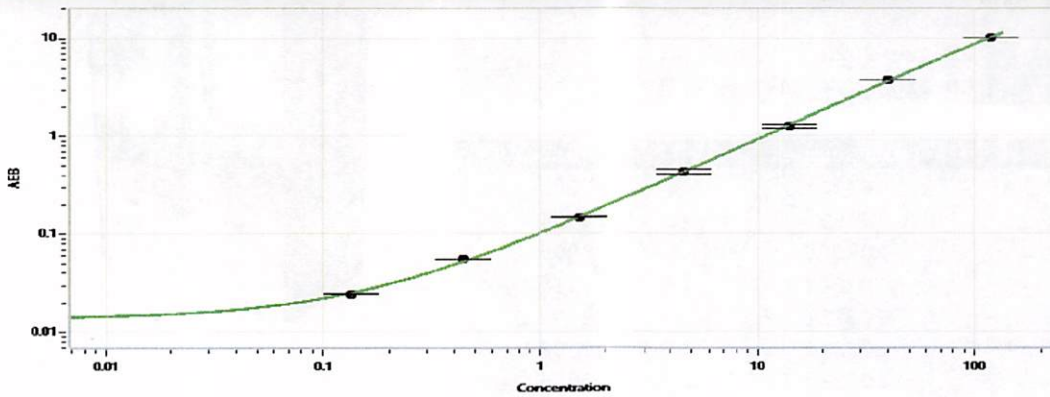
**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	129008	15.4	11.7 - 17.6
Control 2	129008	233	199 - 298
Panel 1	124607	132	115 - 172
Panel 2	124607	98.7	83.1 - 125
Panel 3	124607	166	138 - 207

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

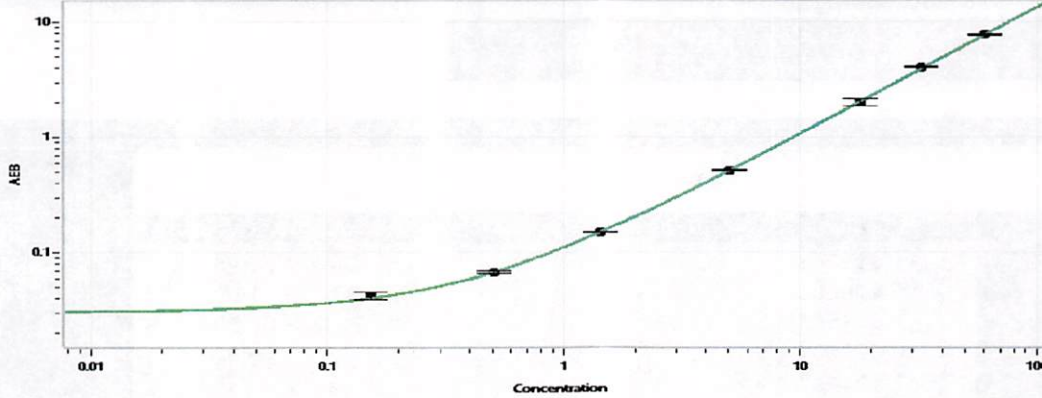
Review/Approval	Junpil Park	QC Supervisor		10 Nov 2021
	Name	Title	Signature/Date	

**Tau Calibration Curve**



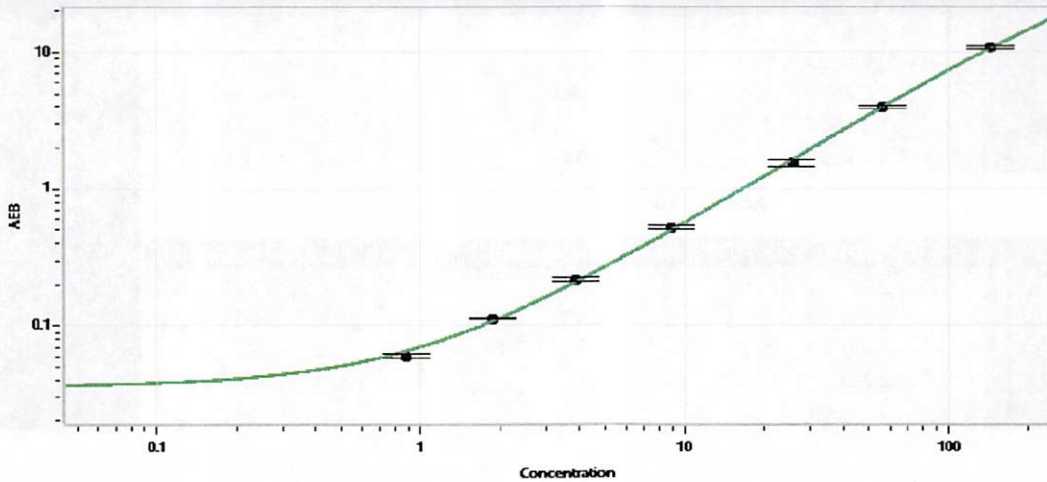
Calibrator Levels (pg/mL)	
A	0.000
B	0.136
C	0.445
D	1.51
E	4.56
F	14.0
G	39.6
H	119

**Abeta 42**



Calibrator Levels (pg/mL)	
A	0.000
B	0.152
C	0.506
D	1.43
E	5.07
F	18.0
G	32.5
H	60.5

**Abeta 40**



Calibrator Levels (pg/mL)	
A	0.000
B	0.885
C	1.87
D	3.87
E	8.90
F	25.5
G	56.0
H	144