

Product Number:	101995
Lot Number:	503095
Expiration:	26-Jan-2022
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101998	113502
Detector Reagent	101996	121713
SBG Reagent	102000	108817
Sample Diluent	102002	108818
Calibrators	103221	114606
Control 1	103222	114607
Control 2	103223	114607
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ 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	114607	2.64	2.08 - 3.11
Control 2	114607	119	106 - 159
Panel 1	003802	1.79	0.896 - 2.74
Panel 2	003802	138	106 - 186
Panel 3	003802	1.45	0.731 - 1.75

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114607	4.78	3.73 - 5.76
Control 2	114607	140	114 - 192
Panel 1	003802	3.75	2.66 - 5.91
Panel 2	003802	3.73	2.74 - 5.10
Panel 3	003802	102	77.7 - 141

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114607	13.2	8.68 - 17.2
Control 2	114607	198	176 - 264
Panel 1	003802	61.9	43.6 - 101
Panel 2	003802	30.1	22.6 - 50.9
Panel 3	003802	77.1	46.9 - 123

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

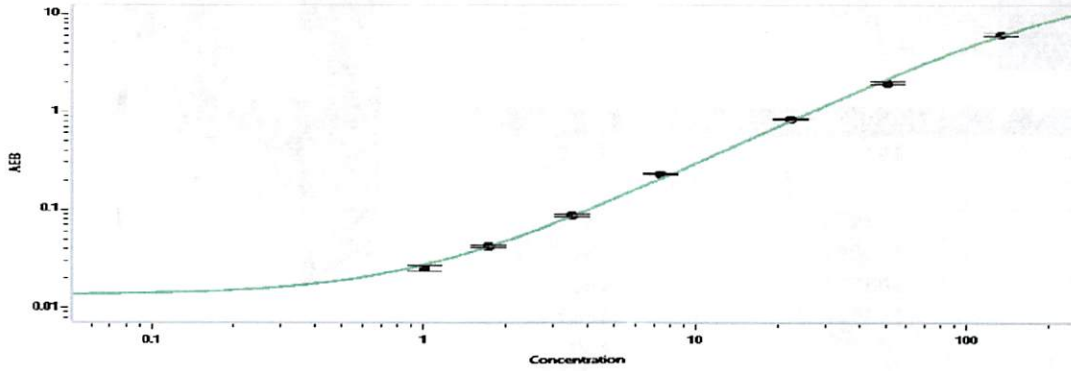
Review/Approval

Junpil Park
Name

QC Supervisor
Title

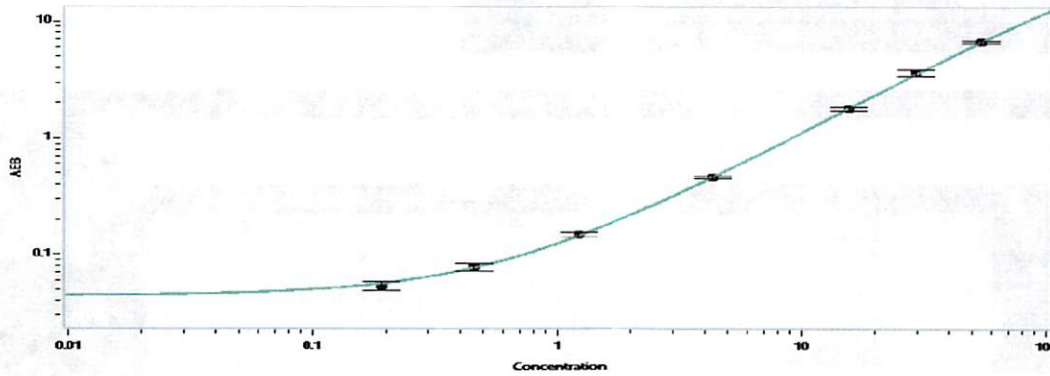
 12 Aug 2021
Signature/Date

Tau Calibration Curve



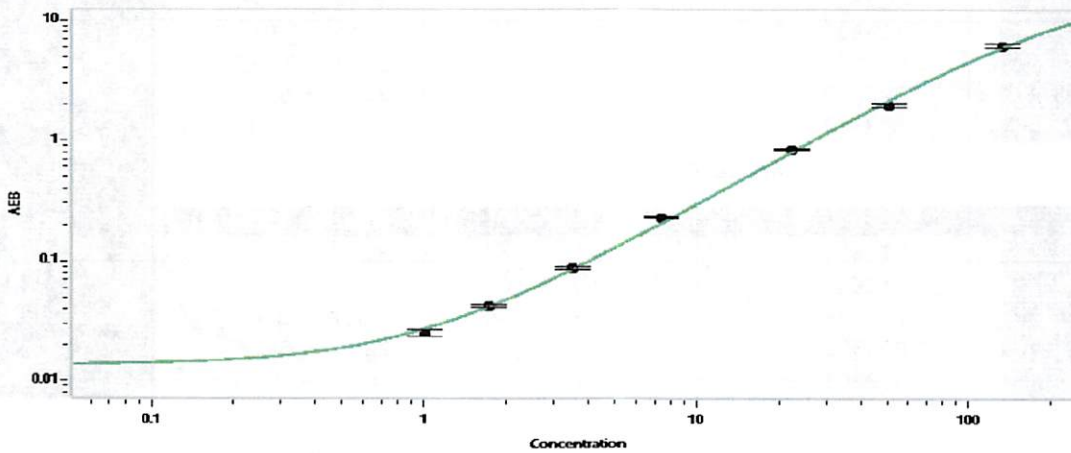
Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.420
D	1.40
E	4.00
F	13.2
G	42.0
H	115

Abeta 42



Calibrator Levels (pg/mL)	
A	0.000
B	0.191
C	0.459
D	1.22
E	4.30
F	15.7
G	29.5
H	54.4

Abeta 40



Calibrator Levels (pg/mL)	
A	0.000
B	1.01
C	1.73
D	3.52
E	7.38
F	22.4
G	50.9
H	132