

Product Number:	101995
Lot Number:	502634
Expiration:	16-Oct-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101998	032402
Detector Reagent	101996	032806
SBG Reagent	102000	032602
Sample Diluent	102002	032805
Calibrators	103221	034402
Control 1	103222	032903
Control 2	103223	032903
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	032903	1.62	0.745-2.23
Control 2	032903	59.9	30.5-91.5
Panel 1	003802	1.55	0.935-2.42
Panel 2	003802	146	106-194
Panel 3	003802	1.32	0.552-1.68

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	032903	4.89	3.95-5.92
Control 2	032903	138	92.4-164
Panel 1	003802	4.36	2.66-6.02
Panel 2	003802	3.81	2.78-5.20
Panel 3	003802	108	88.7-140

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	032903	15.4	8.70-20.3
Control 2	032903	218	143-335
Panel 1	003802	60.7	47.2-101
Panel 2	003802	32.4	26.4-50.9
Panel 3	003802	73.4	50.4-128

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

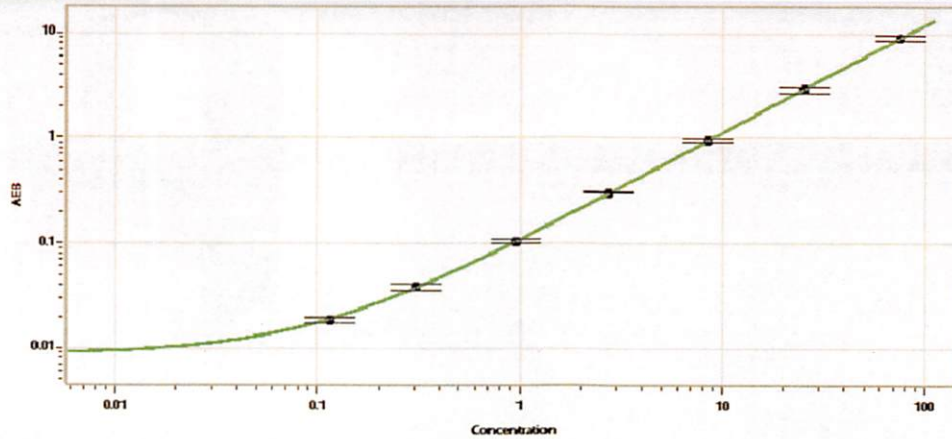
Review/Approval

Junpil Park
Name

QC Supervisor
Title

 16 Dec 2020
Signature/Date

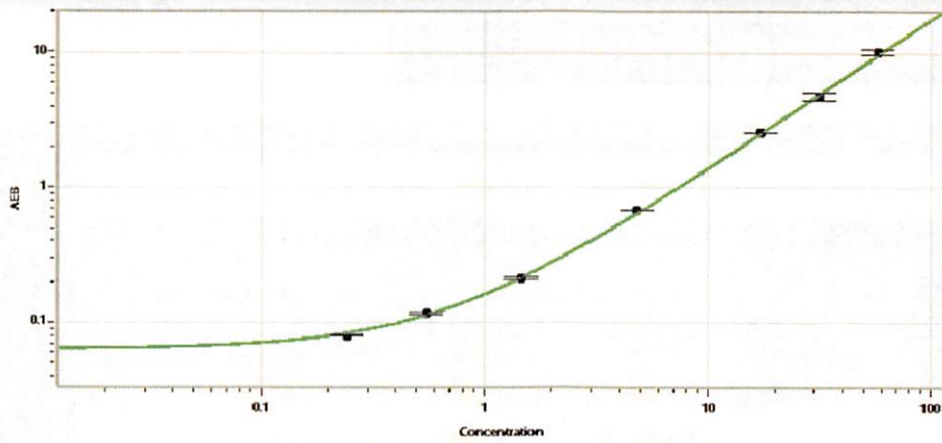
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.114
C	0.305
D	0.964
E	2.73
F	8.49
G	25.3
H	75.6

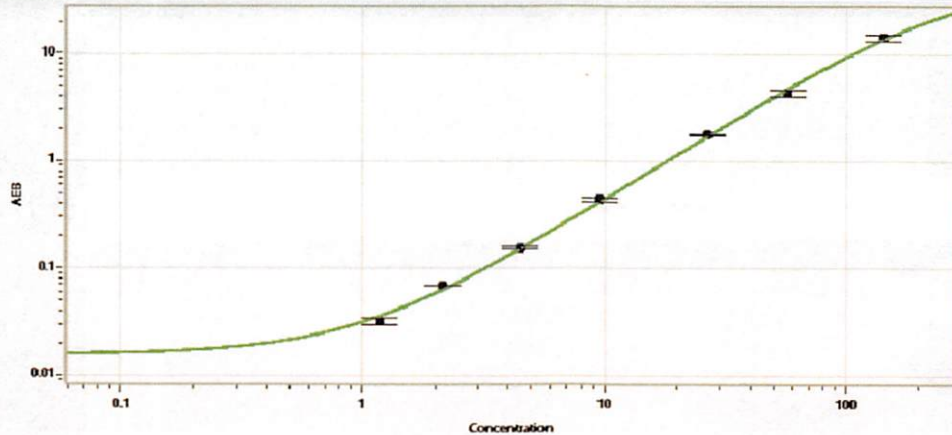
Abeta 42



Calibrator Levels (pg/mL)

A	0.000
B	0.244
C	0.545
D	1.47
E	4.84
F	17.2
G	31.5
H	57.6

Abeta 40



Calibrator Levels (pg/mL)

A	0.000
B	1.18
C	2.14
D	4.49
E	9.58
F	26.4
G	57.0
H	142