

Product Number:	101995
Lot Number:	502838
Expiration:	21-Mar-2022
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101998	108704
Detector Reagent	101996	108705
SBG Reagent	102000	108817
Sample Diluent	102002	108818
Calibrators	103221	A-E: 111902 F-H: 112717
Control 1	103222	112718
Control 2	103223	112718
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	112718	2.35	1.84-2.77
Control 2	112718	111	80.7-138
Panel 1	003802	2.27	0.896-2.74
Panel 2	003802	140	106-186
Panel 3	003802	1.57	0.731-1.75

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	112718	4.36	2.67-5.08
Control 2	112718	99.4	76.7-131
Panel 1	003802	4.19	2.67-5.91
Panel 2	003802	3.85	2.74-5.10
Panel 3	003802	90.8	77.7-141

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	112718	13.1	10.1-15.2
Control 2	112718	205	171-266
Panel 1	003802	63.4	43.6-101
Panel 2	003802	30.2	22.6-50.9
Panel 3	003802	75.6	46.9-123

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

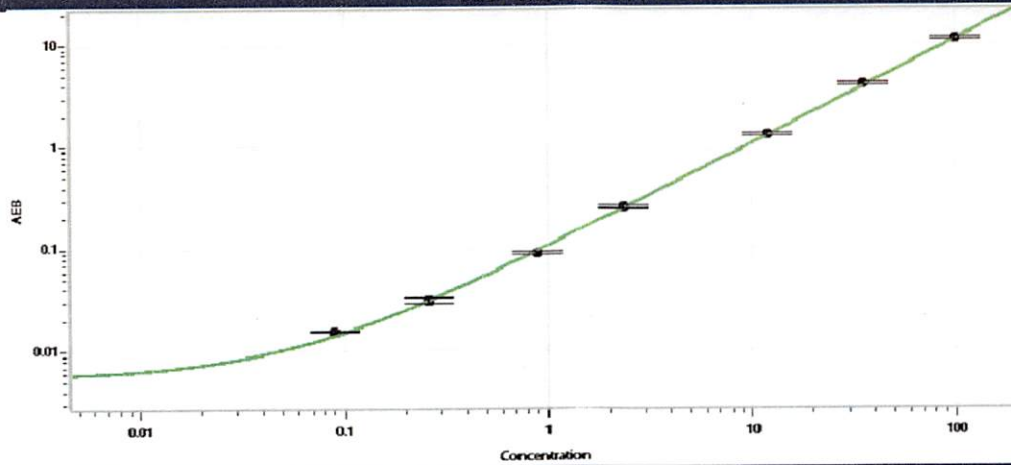
Review/Approval

Junpil Park
Name

QC Supervisor
Title

 20 May 2021
Signature/Date

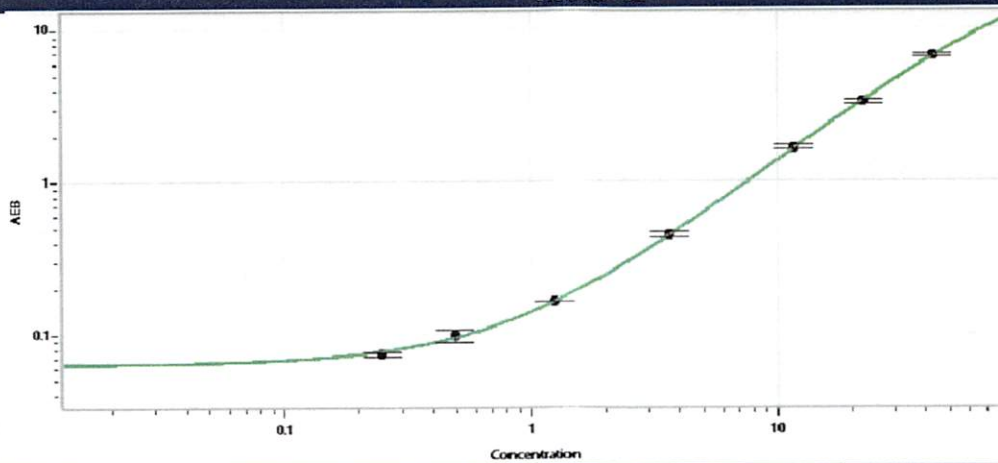
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.090
C	0.261
D	0.890
E	2.36
F	11.9
G	35.4
H	101

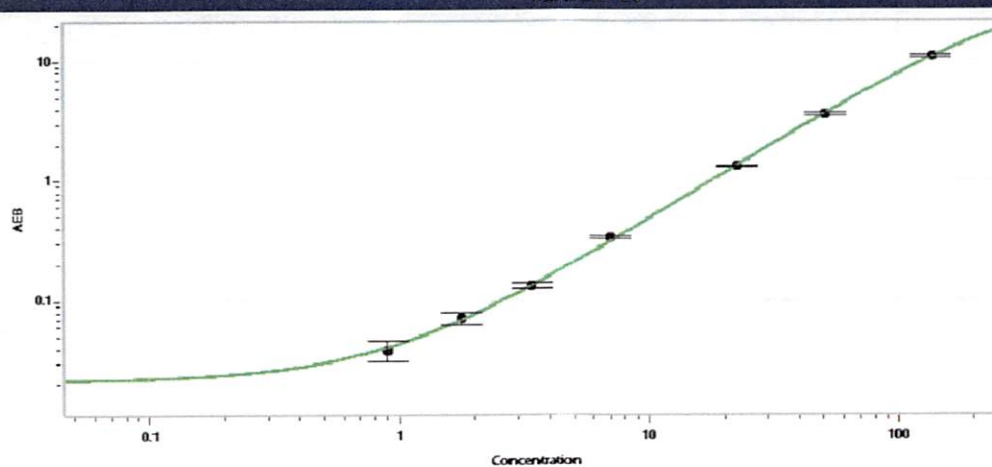
Abeta 42



Calibrator Levels (pg/mL)

A	0.000
B	0.249
C	0.491
D	1.25
E	3.64
F	11.7
G	22.2
H	42.8

Abeta 40



Calibrator Levels (pg/mL)

A	0.000
B	0.900
C	1.78
D	3.42
E	7.00
F	22.7
G	51.1
H	136