

Product Number:	101995
Lot Number:	502804
Expiration:	11-Mar-2022
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101998	107005
Detector Reagent	101996	107006
SBG Reagent	102000	107009
Sample Diluent	102002	107007
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).

Release Materials

Tau

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114607	2.59	2.07-3.10
Control 2	114607	126	109-163
Panel 1	003802	1.67	0.896-2.74
Panel 2	003802	141	106-186
Panel 3	003802	1.36	0.731-1.75

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114607	4.27	3.65-5.48
Control 2	114607	130	108-193
Panel 1	003802	3.42	2.66-5.91
Panel 2	003802	3.28	2.74-5.10
Panel 3	003802	88.6	77.7-141

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114607	12.4	9.65-14.5
Control 2	114607	190	175-263
Panel 1	003802	59.5	43.6-101
Panel 2	003802	29.3	22.6-50.9
Panel 3	003802	69.4	46.9-123

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

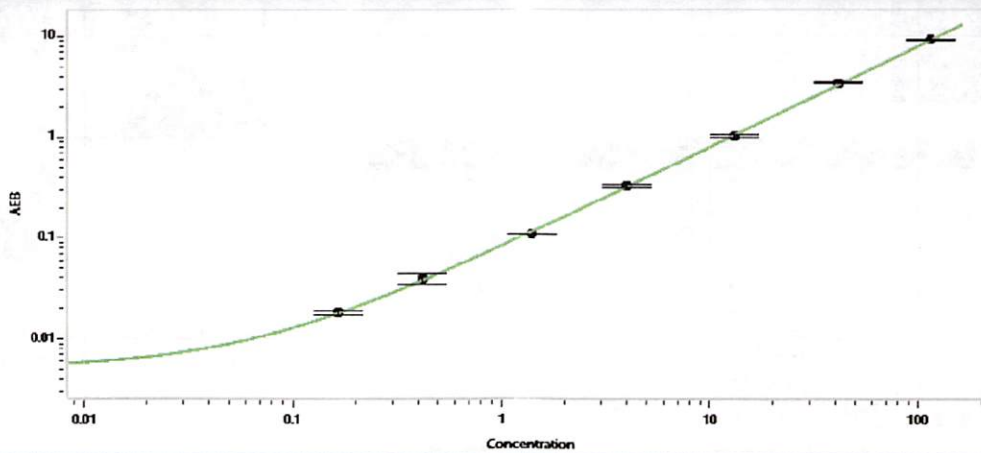
Review/Approval

Jennifer Park
Name

QC Supervisor
Title

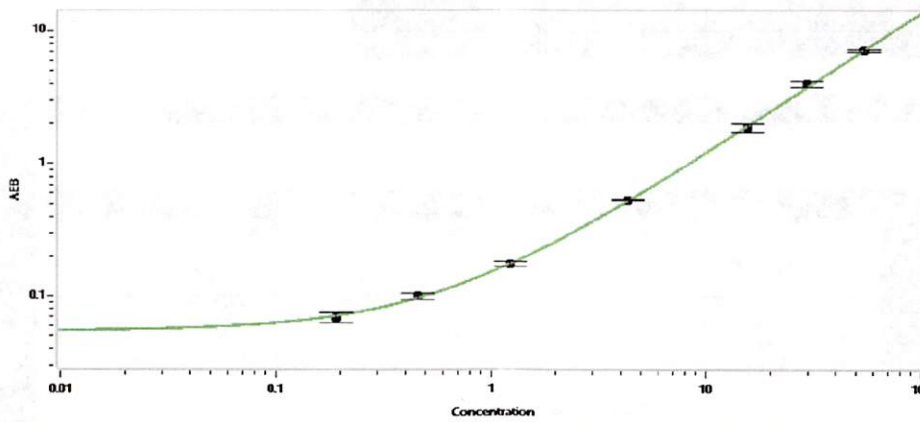
[Signature] 04 Jun 2021
Signature/Date

Tau Calibration Curve



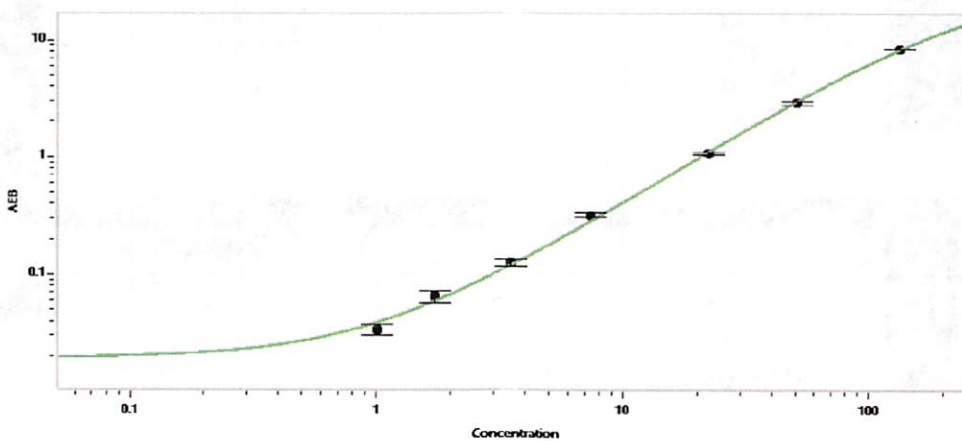
Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.420
D	1.40
E	4.00
F	13.2
G	42.0
H	115

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