

Product Number:	101995
Lot Number:	503240
Expiration:	22-Aug-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	132701
Detector Reagent	101996	134203
SBG Reagent	102000	132702
Sample Diluent	102002	123412
Calibrators	103221	133508
Control 1	103222	133509
Control 2	103223	133509
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	133509	3.09	2.47 - 3.71
Control 2	133509	136	109 - 163
Panel 1	124607	3.46	2.55 - 3.83
Panel 2	124607	230	176 - 264
Panel 3	124607	5.05	3.49 - 5.92

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	133509	5.15	3.10 - 8.19
Control 2	133509	131	111 - 167
Panel 1	124607	6.35	4.54 - 7.59
Panel 2	124607	6.08	4.50 - 7.18
Panel 3	124607	106	79.1 - 126

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	133509	15.7	12.5 - 18.7
Control 2	133509	228	189 - 283
Panel 1	124607	133	113 - 170
Panel 2	124607	99.8	81.8 - 123
Panel 3	124607	173	134 - 203

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

Review/Approval

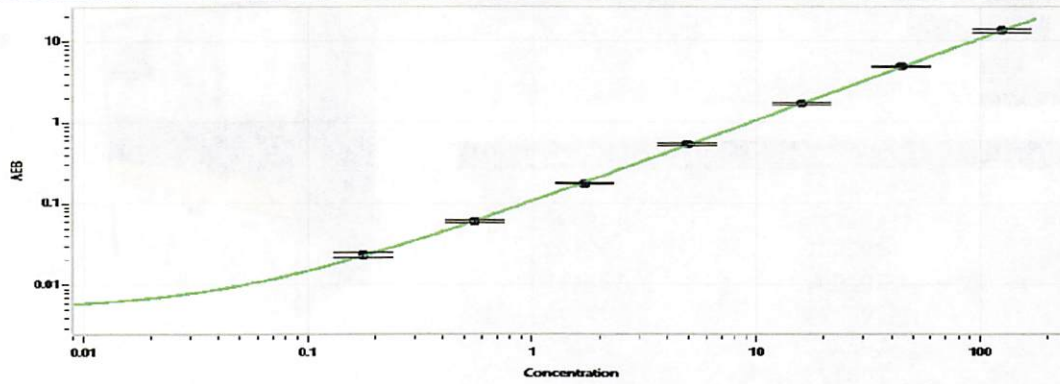
Wanku To.
Name

QC

Title

[Signature] 07/01/2022
Signature/Date

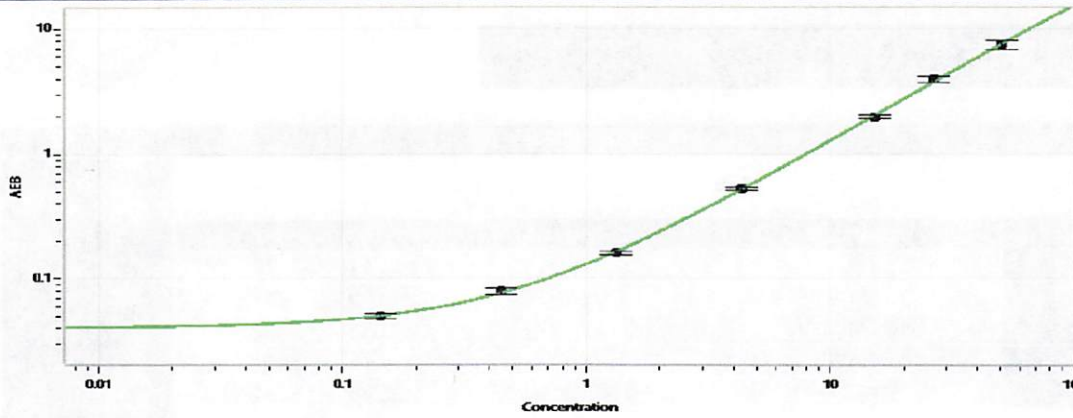
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.178
C	0.557
D	1.72
E	4.90
F	15.9
G	44.2
H	125

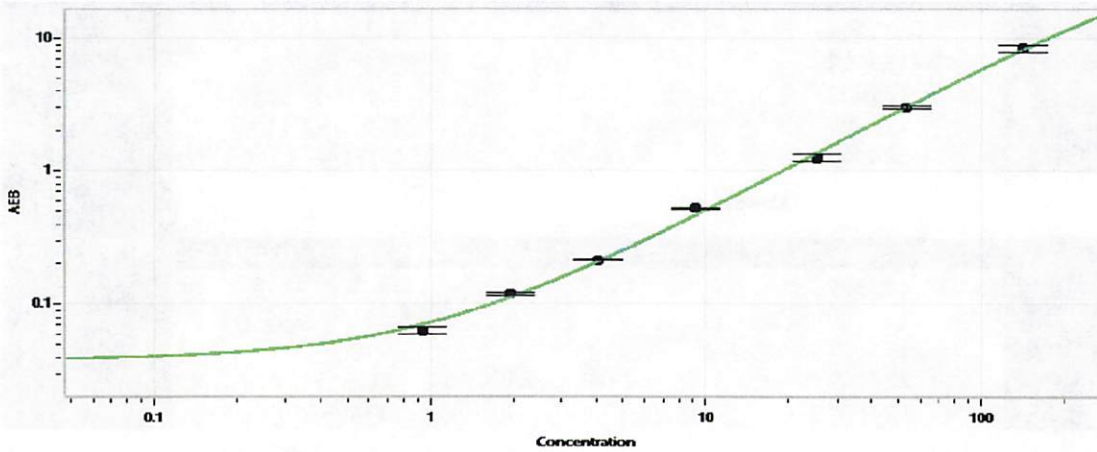
Abeta 42



Calibrator Levels (pg/mL)

A	0.000
B	0.144
C	0.450
D	1.32
E	4.30
F	15.2
G	26.4
H	50.6

Abeta 40



Calibrator Levels (pg/mL)

A	0.000
B	0.94
C	1.96
D	4.07
E	9.15
F	25.3
G	53.1
H	142