

Product Number:	101995
Lot Number:	503099
Expiration:	5-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	122305
Detector Reagent	101996	131608
SBG Reagent	102000	122310
Sample Diluent	102002	122311
Calibrators	103221	122204
Control 1	103222	122205
Control 2	103223	122205
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	122205	2.44	1.91 - 2.87
Control 2	122205	90.3	78.1 - 125
Panel 1	124607	1.32	1.13-1.70
Panel 2	124607	216	191-329
Panel 3	124607	2.16	1.81-2.71

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	122205	4.37	3.67 - 5.50
Control 2	122205	104	80.9 - 150
Panel 1	124607	6.31	4.96-10.6
Panel 2	124607	5.48	4.66-7.85
Panel 3	124607	81.0	57.8-180

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	122205	14.5	10.1 - 15.6
Control 2	122205	202	166 - 248
Panel 1	124607	125	107-169
Panel 2	124607	90.3	78.6-121
Panel 3	124607	137	114-210

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

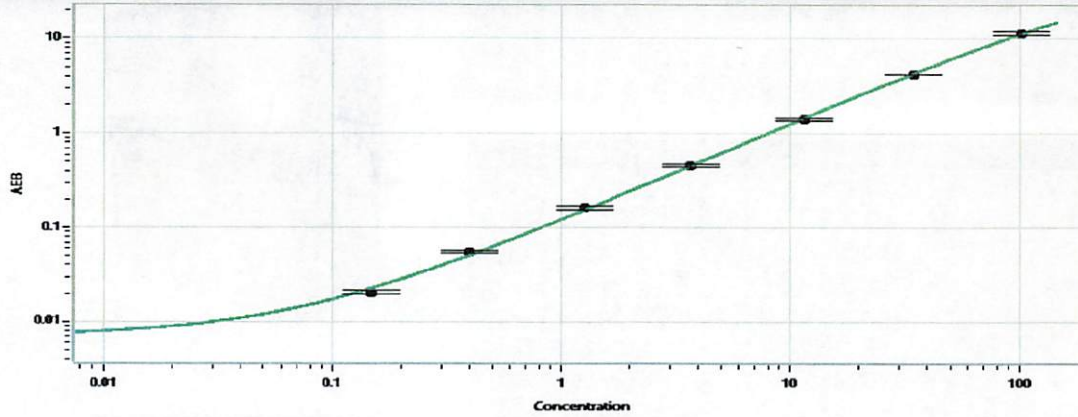
Review/Approval

Junpil Park
Name

QC Supervisor
Title

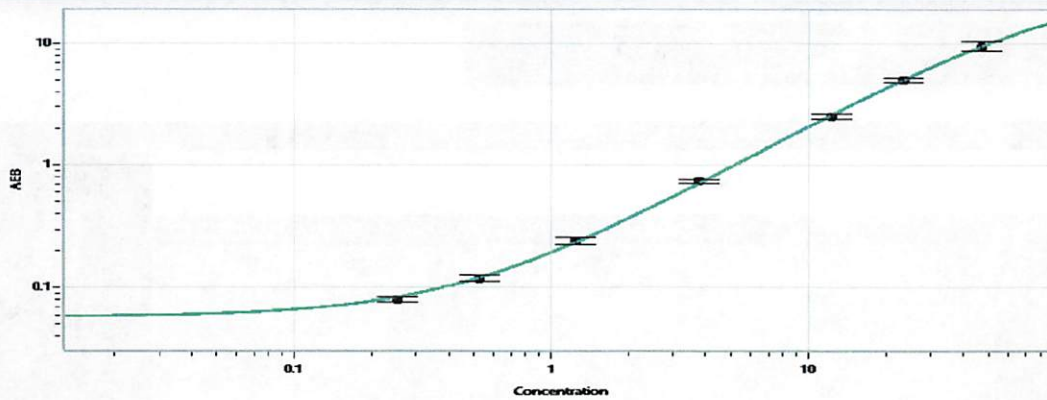
 23 NOV 2021
Signature/Date

Tau Calibration Curve



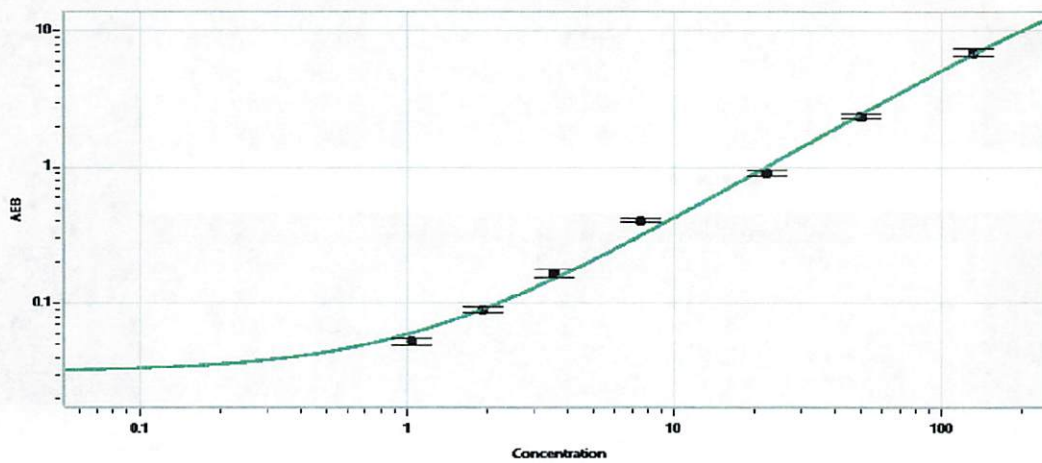
Calibrator Levels (pg/mL)	
A	0.000
B	0.148
C	0.399
D	1.27
E	3.66
F	11.6
G	34.3
H	102

Abeta 42



Calibrator Levels (pg/mL)	
A	0.000
B	0.254
C	0.526
D	1.24
E	3.76
F	12.3
G	23.5
H	46.6

Abeta 40



Calibrator Levels (pg/mL)	
A	0.000
B	1.04
C	1.93
D	3.57
E	7.54
F	22.1
G	49.6
H	132