

Product Number:	101995
Lot Number:	503104
Expiration:	17-Aug-2022
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101998	122905
Detector Reagent	101996	126609
SBG Reagent	102000	122919
Sample Diluent	102002	124608
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).



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

Release Materials

Tau

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	122205	2.58	1.90-2.85
Control 2	122205	103	83.2-125
Panel 1	124607	1.71	1.10-1.93
Panel 2	124607	231	207-320
Panel 3	124607	3.07	1.50-3.47

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	122205	4.87	3.71-5.56
Control 2	122205	118	83.5-154
Panel 1	124607	6.85	5.49-10.3
Panel 2	124607	5.65	4.85-7.75
Panel 3	124607	104	89.3-160

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	122205	13.5	10.0-15.0
Control 2	122205	196	167-250
Panel 1	124607	130	111-167
Panel 2	124607	97.3	81.4-122
Panel 3	124607	165	135-203

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

Review/Approval

Junpil Park

QC Supervisor

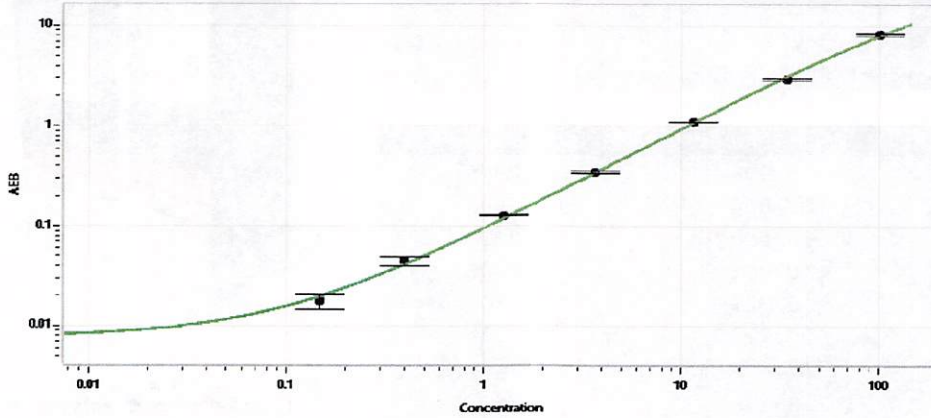
Name

Title

Signature/Date

 15 Oct 2021

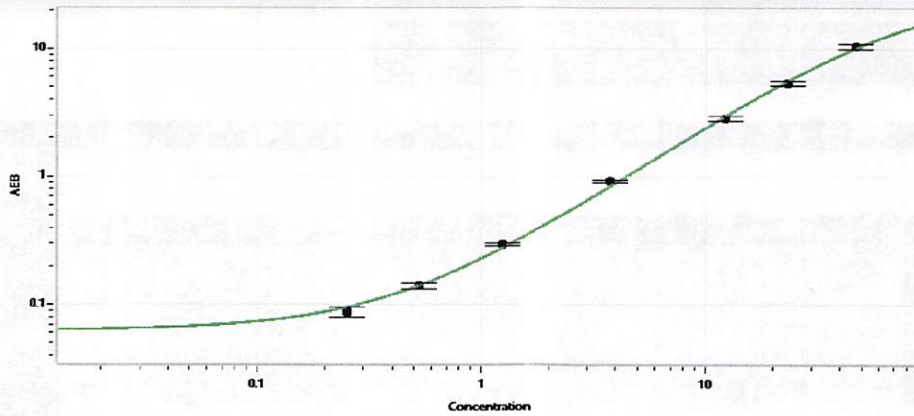
Tau Calibration Curve



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

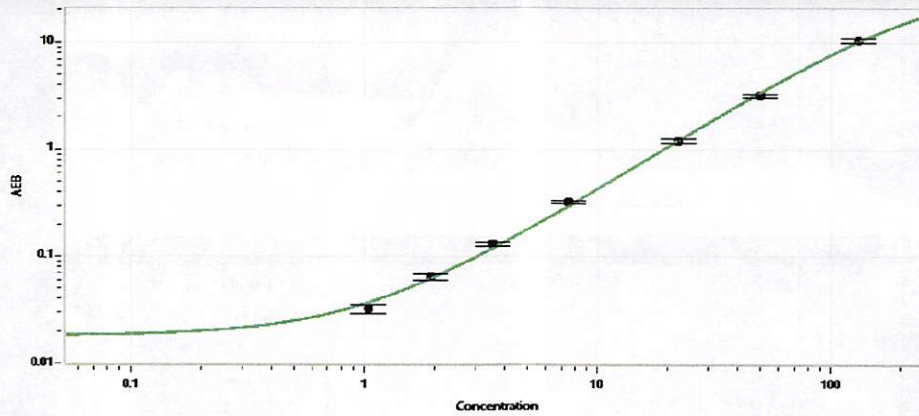
Abeta 42



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