

Product Number:	101995
Lot Number:	503205
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	132009
Detector Reagent	101996	131909
SBG Reagent	102000	132010
Sample Diluent	102002	123412
Calibrators	103221	131504
Control 1	103222	131505
Control 2	103223	131505
SBG Diluent (SRX)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	131505	3.55	1.96-4.20
Control 2	131505	128	70.6-181
Panel 1	124607	2.91	2.55-3.83
Panel 2	124607	220	176-264
Panel 3	124607	4.43	3.49-5.92

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	131505	5.77	4.26-6.71
Control 2	131505	137	90.8-187
Panel 1	124607	5.77	4.54-7.59
Panel 2	124607	5.43	4.50-7.18
Panel 3	124607	89.8	79.1-126

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	131505	16.8	12.0-18.0
Control 2	131505	226	176-289
Panel 1	124607	138	113-170
Panel 2	124607	97.6	81.8-123
Panel 3	124607	156	134-203

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

Review/Approval

Junpil Park

QC Supervisor

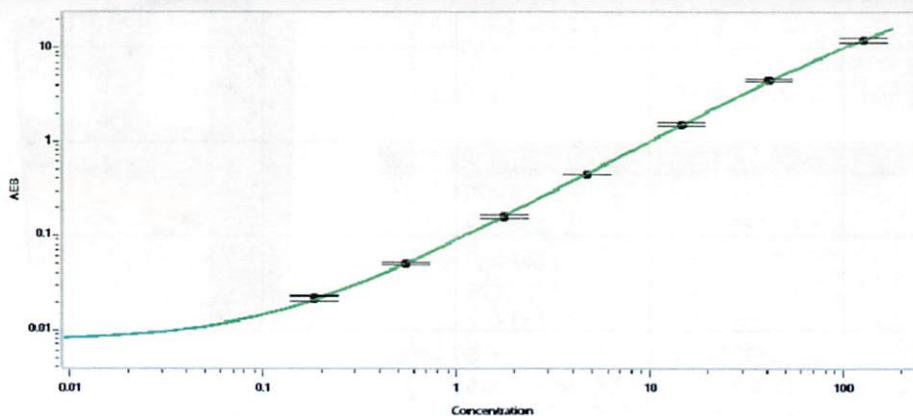
02 Dec 2021

Name

Title

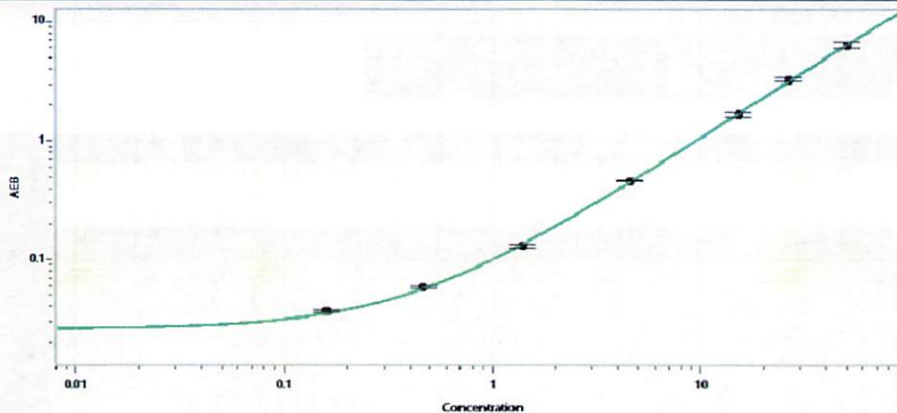
Signature/Date

**Tau Calibration Curve**



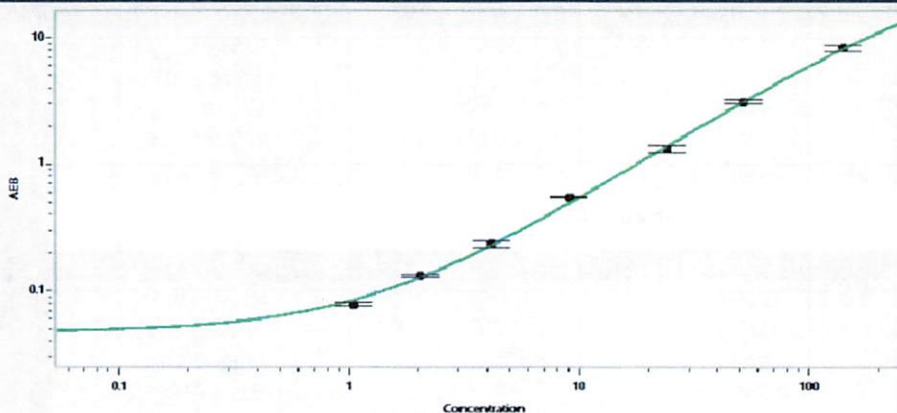
Calibrator Levels (pg/mL)	
A	0.000
B	0.182
C	0.551
D	1.74
E	4.67
F	14.5
G	40.8
H	127

**Abeta 42**



Calibrator Levels (pg/mL)	
A	0.000
B	0.159
C	0.465
D	1.39
E	4.55
F	15.2
G	26.3
H	49.9

**Abeta 40**



Calibrator Levels (pg/mL)	
A	0.000
B	1.05
C	2.03
D	4.13
E	9.07
F	24.2
G	52.0
H	140