

Product Number:	101995
Lot Number:	502072
Expiration:	10-Oct-2020
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	928808
Detector Reagent	101996	928801
SBG Reagent (HD-1)	102000	928302
Sample Diluent	102002	928301
Calibrators	103221	923502 & 924001
Controls	103222/103223	923503
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	923503	1.85	1.11-2.20
Control 2	923503	70.1	41.6-85.6
Panel 1	MF-062018-04	1.62	1.15-2.05
Panel 2	MF-010818-02	185	151-230
Panel 3	MF-010818-03	4.44	1.55-6.20

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	923503	5.48	2.86-7.05
Control 2	923503	119	66.1-146
Panel 1	MF-062018-04	9.02	4.63-9.51
Panel 2	MF-010818-02	13.2	6.14-14.6
Panel 3	MF-010818-03	132	98.9-148

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	923503	11.1	6.72-14.2
Control 2	923503	185	119-217
Panel 1	MF-062018-04	137	104-156
Panel 2	MF-010818-02	149	124-190
Panel 3	MF-010818-03	103	67.6-127

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

Review/Approval

Patrizia Stadler Assoc. Director, QC

*P. Stadler*

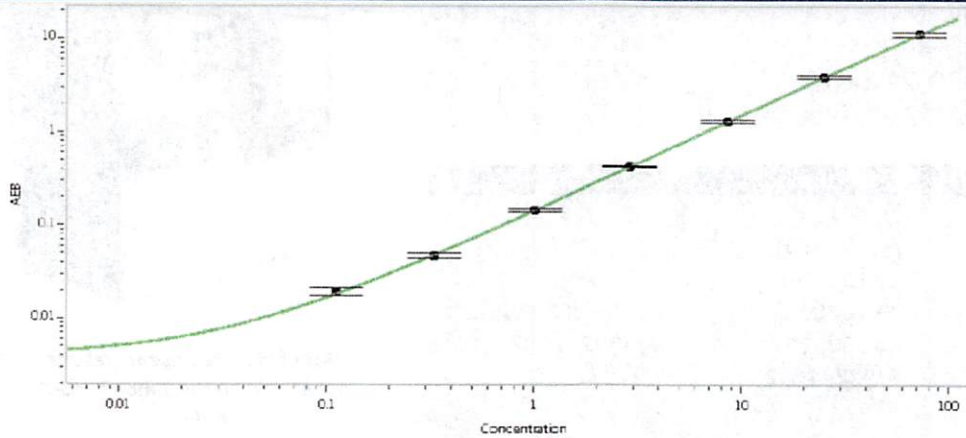
25 NOV 2019

Name

Title

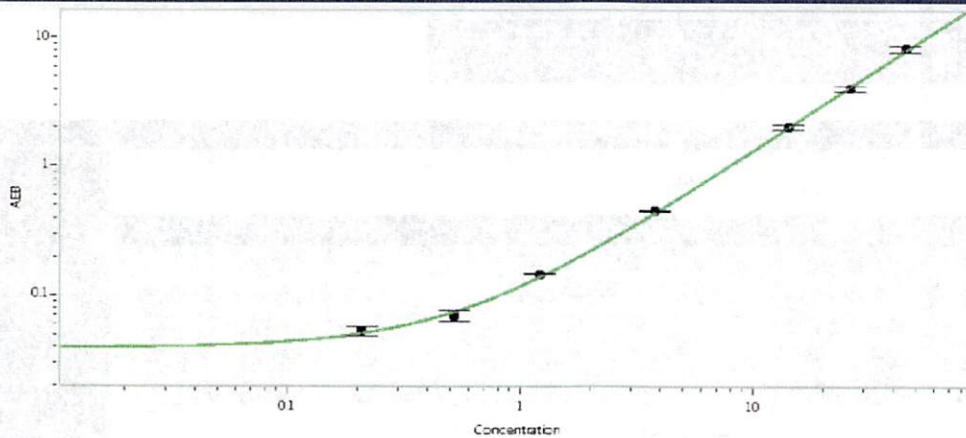
Signature/Date

**Tau Calibration Curve**



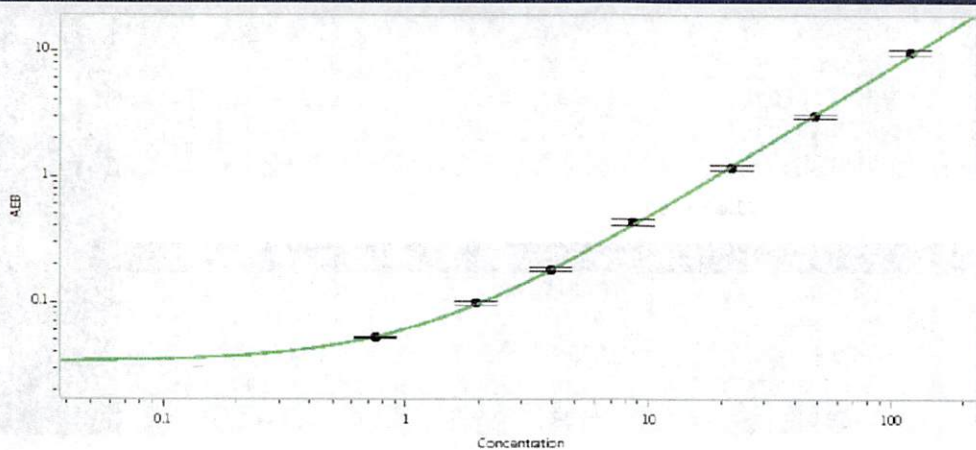
Calibrator Levels (pg/mL)	
A	0.000
B	0.111
C	0.331
D	1.02
E	2.86
F	8.51
G	24.9
H	71.9

**Abeta 42**



Calibrator Levels (pg/mL)	
A	0.000
B	0.211
C	0.524
D	1.22
E	3.83
F	14.2
G	26.3
H	45.3

**Abeta 40**



Calibrator Levels (pg/mL)	
A	0.000
B	0.748
C	1.95
D	3.97
E	8.61
F	22.1
G	48.7
H	119