

Product Number:	101995
Lot Number:	502161
Expiration:	29-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	935210
Detector Reagent	101996	935301
SBG Reagent	102000	934017
Sample Diluent	102002	934016
Calibrators	103221	934702 & 935702
Controls	103222/103223	934703
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	934703	2.81	2.38-3.70
Control 2	934703	109	90.4-151
Panel 1	MF-062018-04	0.644	0.453-2.36
Panel 2	MF-010818-02	170	151-226
Panel 3	MF-010818-03	2.77	1.59-6.18

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	934703	4.90	3.61-6.99
Control 2	934703	111	94.4-142
Panel 1	MF-062018-04	8.83	4.73-10.24
Panel 2	MF-010818-02	12.5	6.06-14.5
Panel 3	MF-010818-03	131	98.3-147

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	934703	14.8	12.2-18.3
Control 2	934703	223	187-281
Panel 1	MF-062018-04	106	92.4-158
Panel 2	MF-010818-02	120	105-196
Panel 3	MF-010818-03	83	69.0-125

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

Review/Approval

Patrizia Stadler Assoc. Director, QC

Name

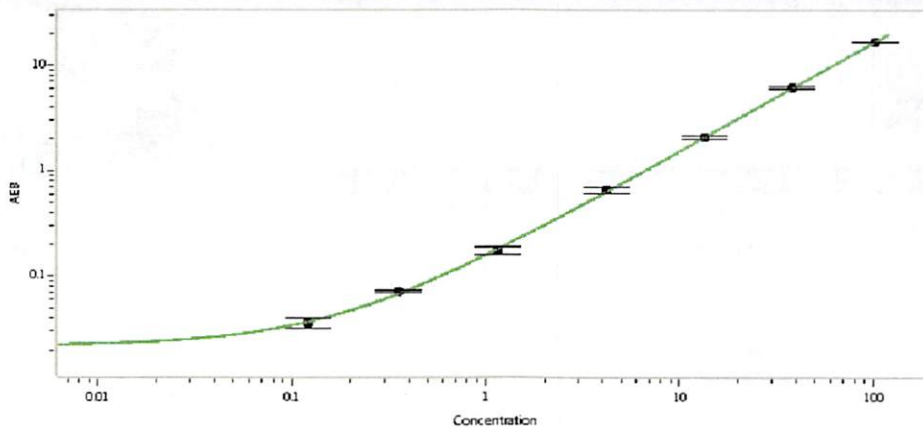
Title

*p. stadler*

30 Dec 2019

Signature/Date

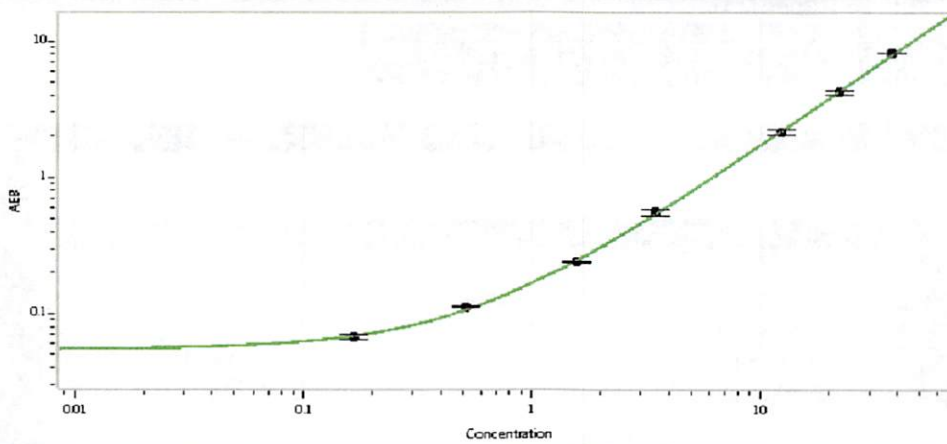
**Tau Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.122
C	0.353
D	1.16
E	4.18
F	13.40
G	38.2
H	102.0

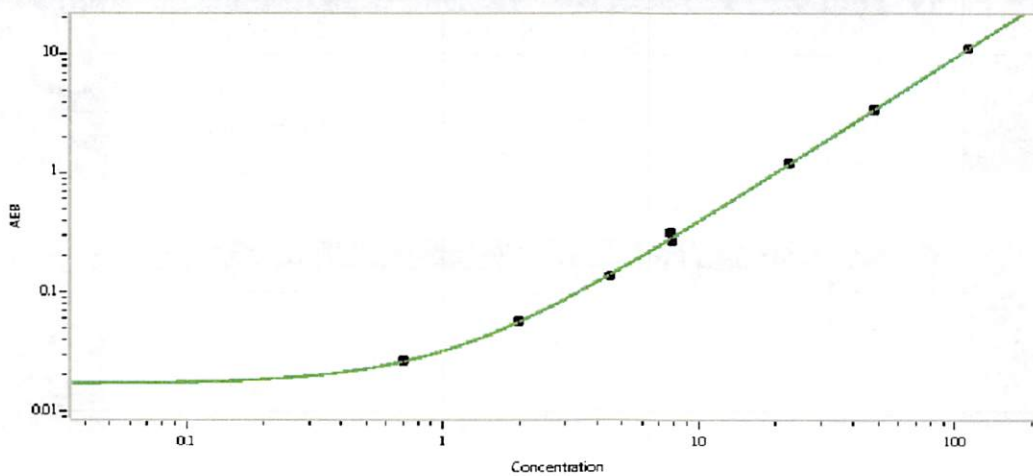
**Abeta 42**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.168
C	0.519
D	1.58
E	3.48
F	12.4
G	22.3
H	37.6

**Abeta 40**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.699
C	1.97
D	4.50
E	7.87
F	22.5
G	48.0
H	112