

Product Number:	101995
Lot Number:	502179
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	102000	928302
Sample Diluent	102002	928301
Calibrators	103221	934702 / 935702
Controls	103222/103223	934703
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	934703	3.36	2.43-3.80
Control 2	934703	120	93.2-148
Panel 1	MF-062018-04	1.87	0.453-2.36
Panel 2	MF-010818-02	192	151-226
Panel 3	MF-010818-03	4.58	1.59-6.18

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	934703	5.39	3.82-6.83
Control 2	934703	113	93.6-140
Panel 1	MF-062018-04	9.30	4.73-10.2
Panel 2	MF-010818-02	12.9	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	15.4	12.2-18.4
Control 2	934703	229	186-279
Panel 1	MF-062018-04	133	92.4-158
Panel 2	MF-010818-02	141	105-196
Panel 3	MF-010818-03	94.1	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

P. Stadler

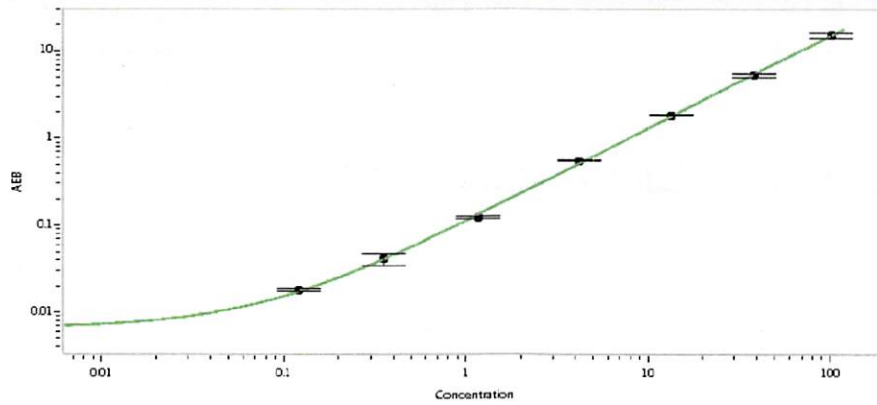
09 Jan 2020

Name

Title

Signature/Date

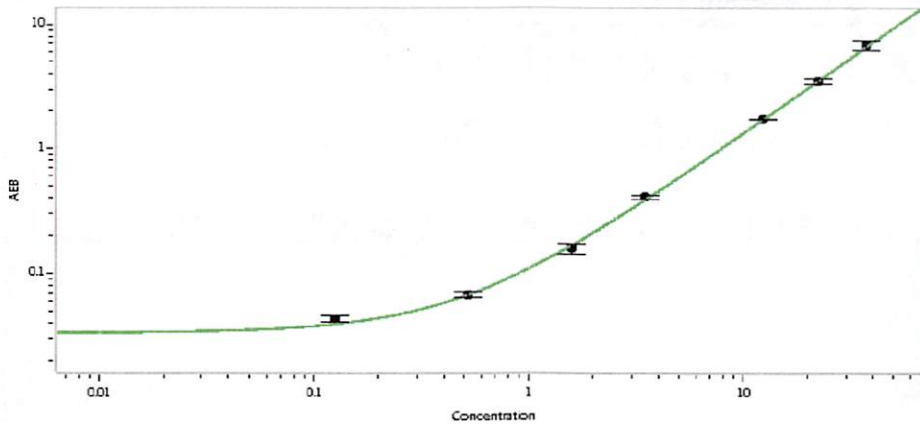
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.122
C	0.353
D	1.16
E	4.18
F	13.4
G	38.2
H	102

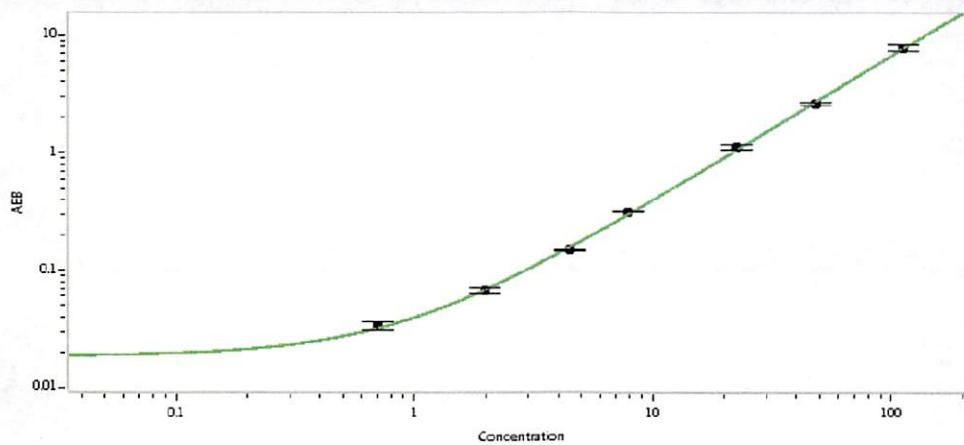
Abeta 42



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