

Product Number:	101995
Lot Number:	502216
Expiration:	21-Nov-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	001513
Detector Reagent	101996	935301
SBG Reagent	102000	934017
Sample Diluent	102002	934016
Calibrators	103221	934702 / 935702
Controls	103222/103223	931002
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	931002	1.43	0.768-1.80
Control 2	931002	60.2	43.5-77.2
Panel 1	MF-062018-04	1.99	0.453-2.36
Panel 2	MF-010818-02	178	151-226
Panel 3	MF-010818-03	2.60	1.59-6.18

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931002	5.68	4.70-7.05
Control 2	931002	138	107-161
Panel 1	MF-062018-04	8.94	4.73-10.2
Panel 2	MF-010818-02	12.4	6.06-14.5
Panel 3	MF-010818-03	127	98.3-147

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931002	16.7	13.2-19.8
Control 2	931002	249	211-317
Panel 1	MF-062018-04	103	92.4-158
Panel 2	MF-010818-02	125	105-196
Panel 3	MF-010818-03	74.7	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

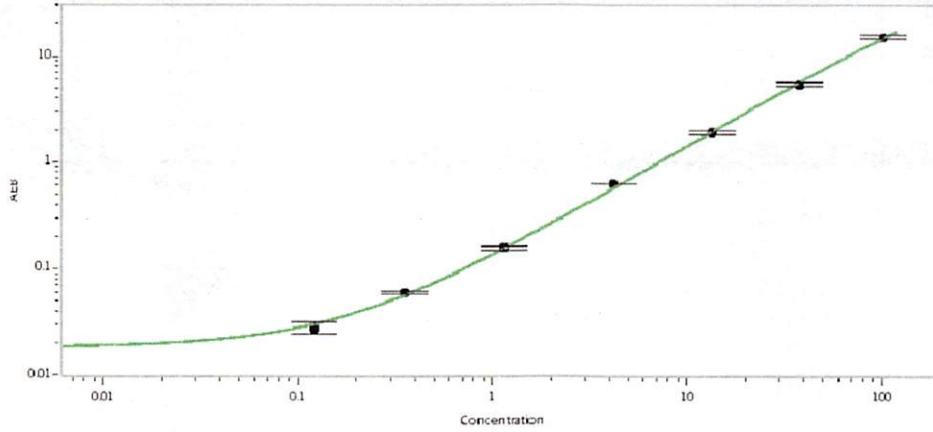
Name

Title

P. Stadler 30 Jan 2020

Signature/Date

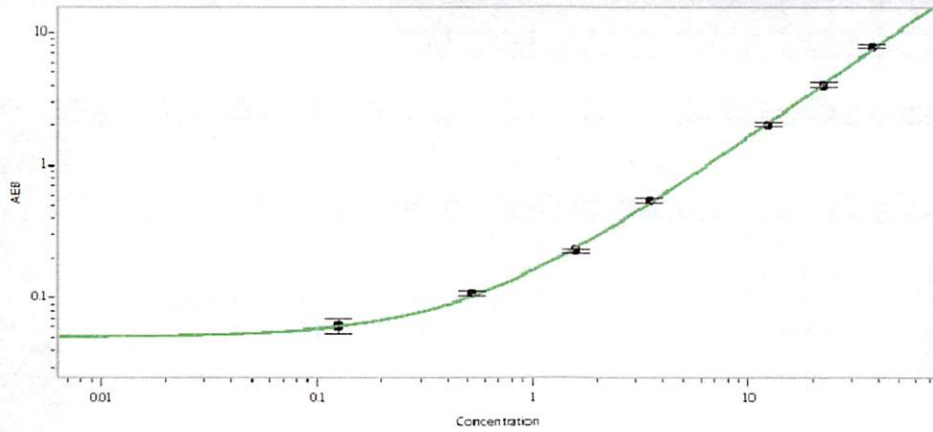
Tau Calibration Curve



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

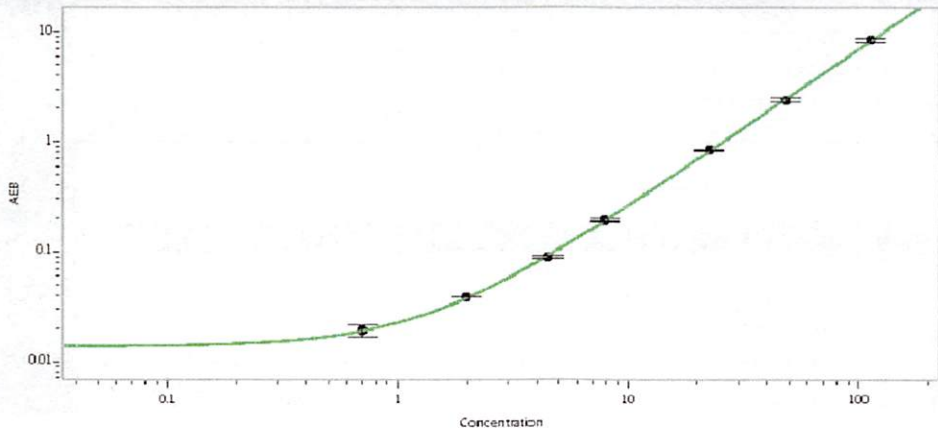
Abeta 42



Calibrator Levels (pg/mL)

A	0.000
B	0.125
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