

Product Number:	101995
Lot Number:	502013
Expiration:	30-Jul-2020
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101998	921704
Detector Reagent	101996	921708
SBG Reagent (HD-1)	102000	921203
Sample Diluent	102002	921107
Calibrators	103221	923502 & 924001
Controls	103222/103223	923503
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	923503	1.91	1.09-2.17
Control 2	923503	74.8	41.4-82.9
Panel 1	MF-062018-04	1.77	1.18-2.03
Panel 2	MF-010818-02	179	153-230
Panel 3	MF-010818-03	4.63	1.36-6.23

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	923503	5.39	2.85-6.77
Control 2	923503	116	66.5-139
Panel 1	MF-062018-04	7.69	5.32-7.98
Panel 2	MF-010818-02	11.2	6.52-12.2
Panel 3	MF-010818-03	134	98.0-147

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	923503	10.7	7.38-12.7
Control 2	923503	178	127-200
Panel 1	MF-062018-04	140	104-159
Panel 2	MF-010818-02	154	123-191
Panel 3	MF-010818-03	108	71.4-125

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

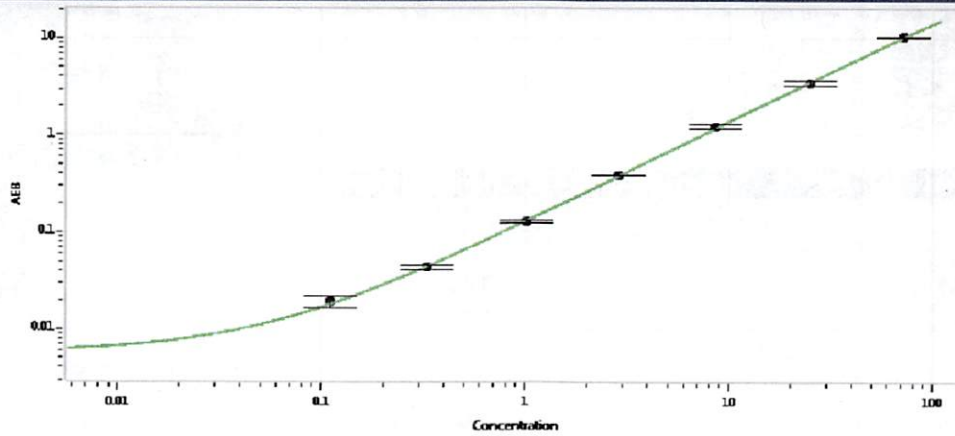
23 Sep 2019

Name

Title

Signature/Date

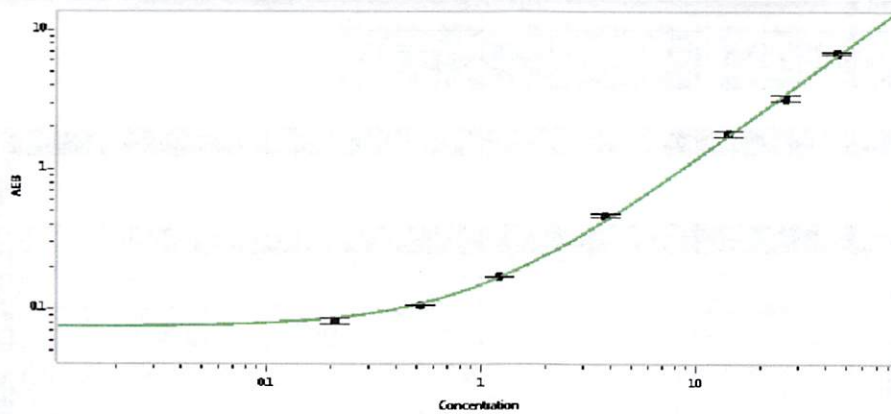
Tau Calibration Curve



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

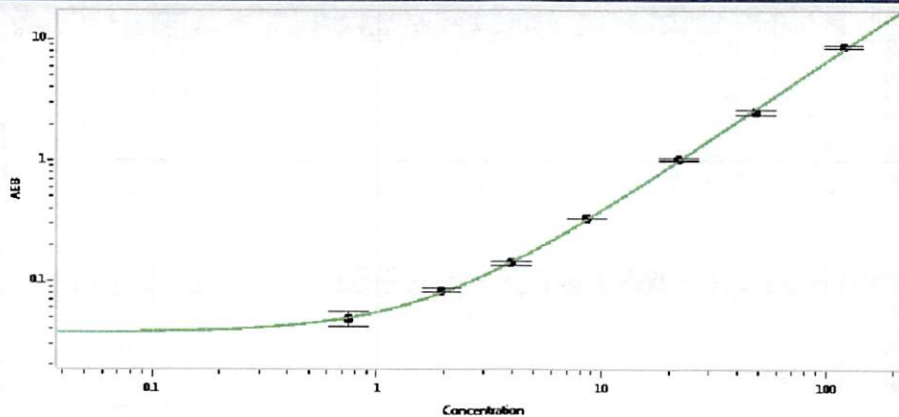
Abeta 42



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