

Product Number:	101995
Lot Number:	501736
Expiration:	19-Jun-2019
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101998	817905
Detector Reagent	101996	817906
SBG Reagent (HD-1)	102000	817010
Sample Diluent	102002	817011
Calibrators	103221	903702/904203
Controls	103222/103223	903703
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

**Release Materials**

**Tau**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	903703	1.03	0.893-1.38
Control 2	903703	46.6	36.0-54.0
Panel 1	MF-062018-04	1.54	1.16-2.04
Panel 2	MF-010818-02	182	156-234
Panel 3	MF-010818-03	4.10	1.30-5.98

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	903703	4.60	3.62-5.43
Control 2	903703	88.5	67.8-101
Panel 1	MF-062018-04	5.86	5.28-7.92
Panel 2	MF-010818-02	9.08	7.07-10.7
Panel 3	MF-010818-03	124	96.1-150

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	903703	14.4	12.2-20.9
Control 2	903703	221	181-272
Panel 1	MF-062018-04	109	102-158
Panel 2	MF-010818-02	130	107-200
Panel 3	MF-010818-03	84.0	71.5-124

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

Review/Approval

Patrizia Stadler Assoc. Director, QC

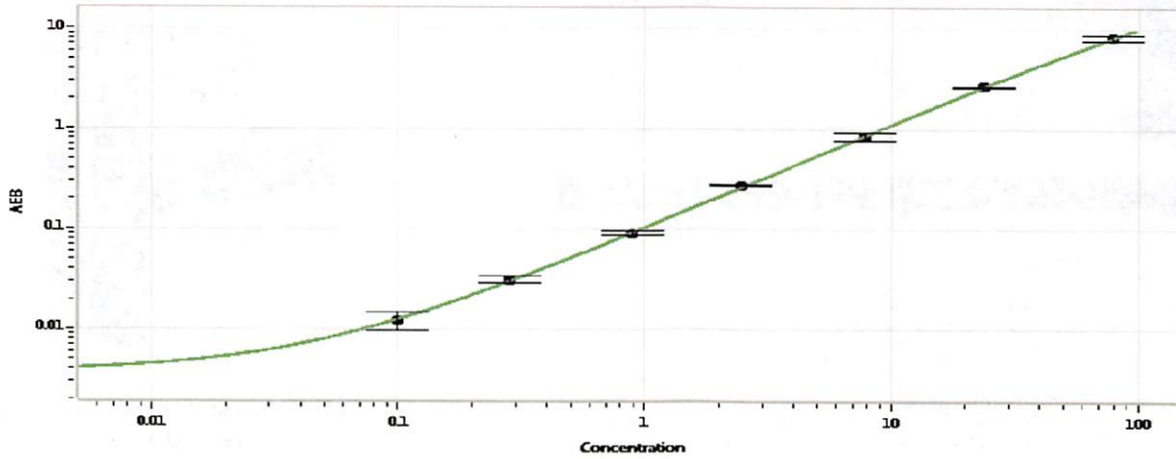
*P. Stadler* 26 Mar 2019

Name

Title

Signature/Date

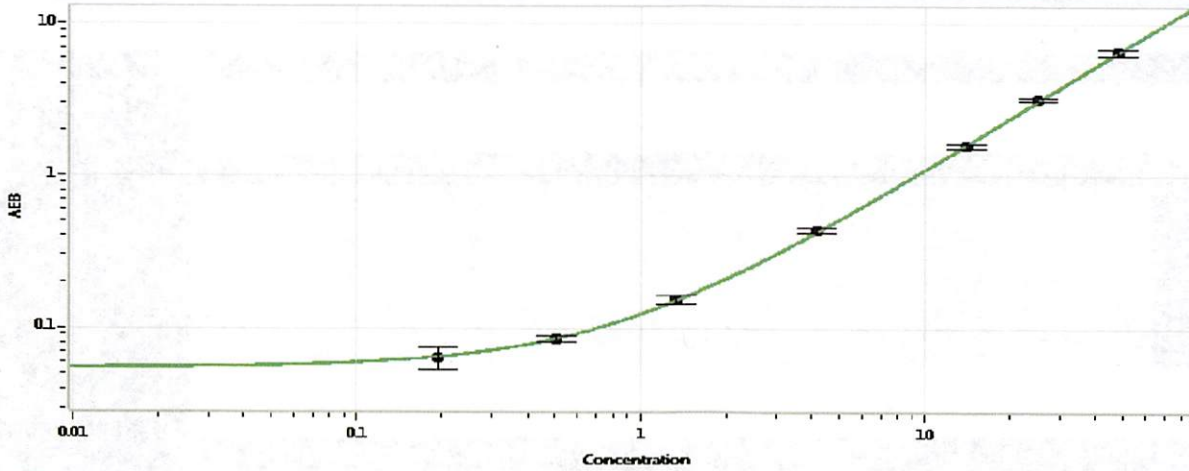
**Tau Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.100
C	0.283
D	0.886
E	2.44
F	7.82
G	23.8
H	79.0

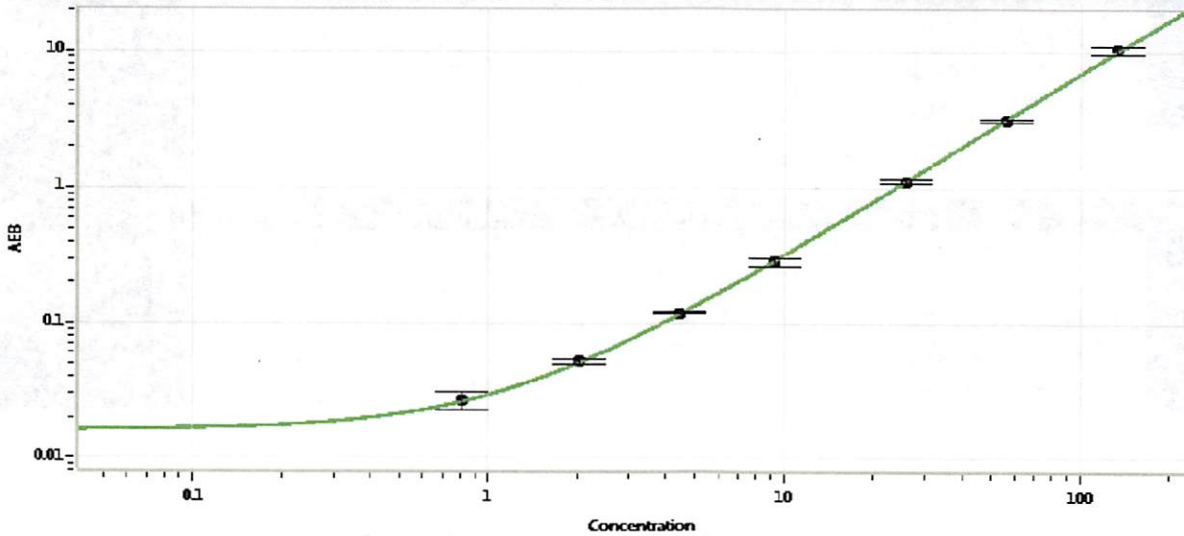
**Abeta 42**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.194
C	0.507
D	1.34
E	4.20
F	14.1
G	25.1
H	47.9

**Abeta 40**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.814
C	2.04
D	4.45
E	9.27
F	25.9
G	56.1
H	133