

Product Number:	101995
Lot Number:	501742
Expiration:	15-Mar-2020
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	910703
Detector Reagent	101996	910704
SBG Reagent (HD-1)	102000	907403
Sample Diluent	102002	907402
Calibrators	103221	911201
Controls	103222/103223	911202
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	911202	1.92	1.47-2.21
Control 2	911202	73.6	58.3-87.5
Panel 1	MF-062018-04	1.82	1.16-2.04
Panel 2	MF-010818-02	174	156-234
Panel 3	MF-010818-03	4.43	1.30-5.98

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911202	4.67	3.77-5.66
Control 2	911202	102	78.9-118
Panel 1	MF-062018-04	6.69	5.28-7.92
Panel 2	MF-010818-02	8.60	7.07-10.7
Panel 3	MF-010818-03	114	96.1-150

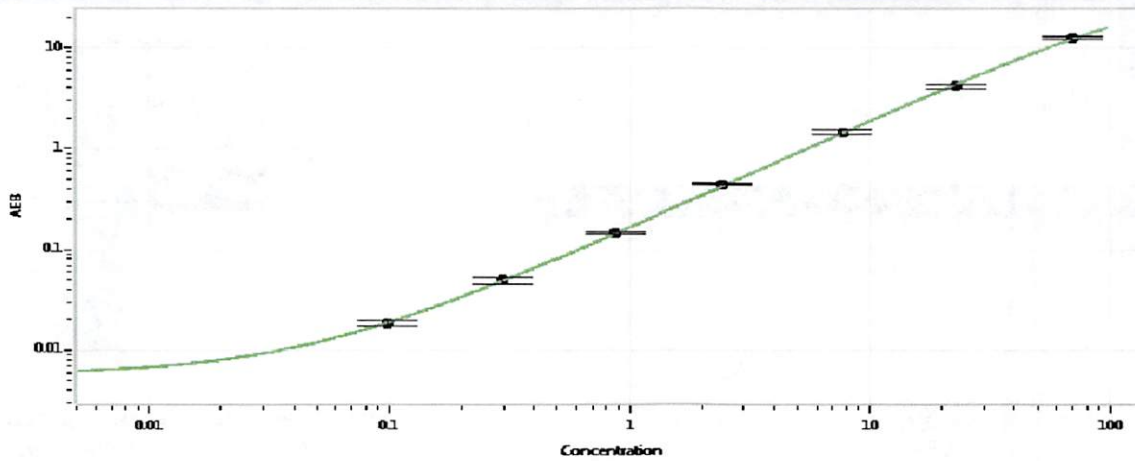
**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911202	14.7	11.6-17.4
Control 2	911202	238	182-273
Panel 1	MF-062018-04	150	102-158
Panel 2	MF-010818-02	170	107-200
Panel 3	MF-010818-03	104	71.5-124

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

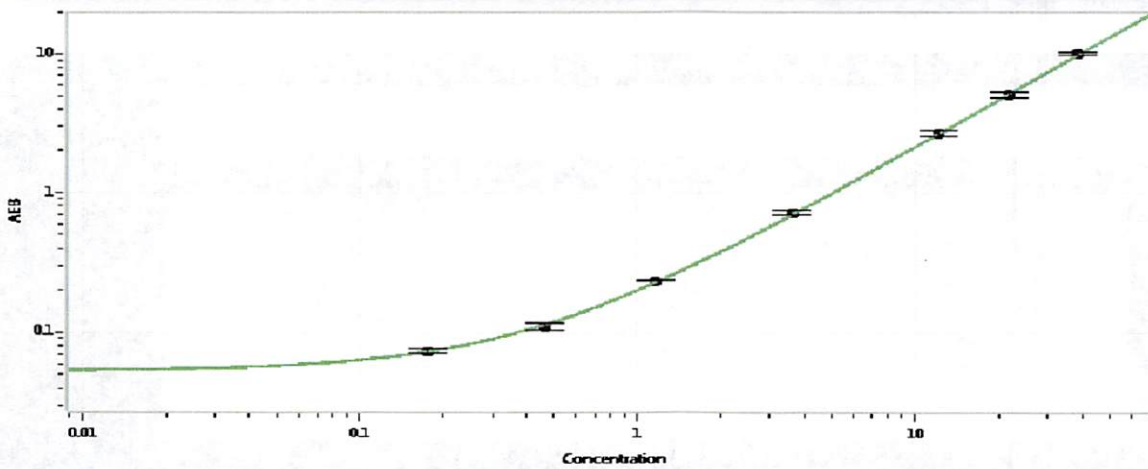
Review/Approval	Patrizia Stadler Assoc. Director, QC	<i>P. Stadler</i>	29 Apr 2019
	Name	Title	Signature/Date

**Tau Calibration Curve**



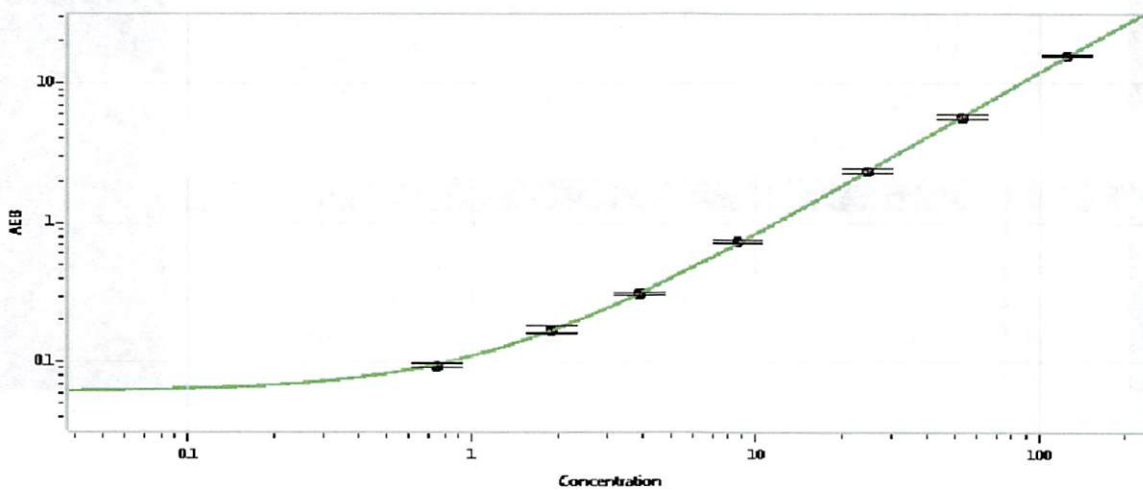
Calibrator Levels (pg/mL)	
A	0.000
B	0.099
C	0.296
D	0.879
E	2.42
F	7.70
G	23.0
H	70.3

**Abeta 42**



Calibrator Levels (pg/mL)	
A	0.000
B	0.176
C	0.463
D	1.16
E	3.60
F	12.2
G	21.8
H	38.2

**Abeta 40**



Calibrator Levels (pg/mL)	
A	0.000
B	0.752
C	1.90
D	3.88
E	8.62
F	24.6
G	52.9
H	124