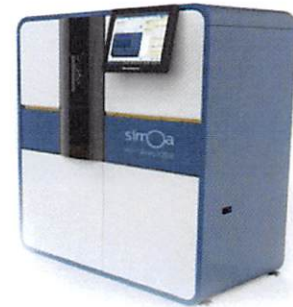


Product Number:	101995
Lot Number:	501509
Expiration:	19-Jun-2019
Platform(s):	HD-1 & SR-X



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	102000	817010
Sample Diluent	102002	817011
Calibrators	103221	820001
Controls	102005-102006	80172
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0083).

**Release Materials**

**Tau**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80172	2.25	1.64-2.46
Control 2	80172	92.4	70.8-108
Panel 1	MF-062018-04	1.84	1.29-1.96
Panel 2	MF-010818-02	168	156-234
Panel 3	MF-010818-03	4.73	1.30-5.98

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80172	5.79	4.74-7.11
Control 2	80172	118	106-159
Panel 1	MF-062018-04	6.27	5.32-7.98
Panel 2	MF-010818-02	8.58	7.07-10.7
Panel 3	MF-010818-03	116	96.1-150

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80172	15.7	13.8-20.7
Control 2	80172	225	194-291
Panel 1	MF-062018-04	114	106-159
Panel 2	MF-010818-02	126	107-200
Panel 3	MF-010818-03	90.1	71.5-124

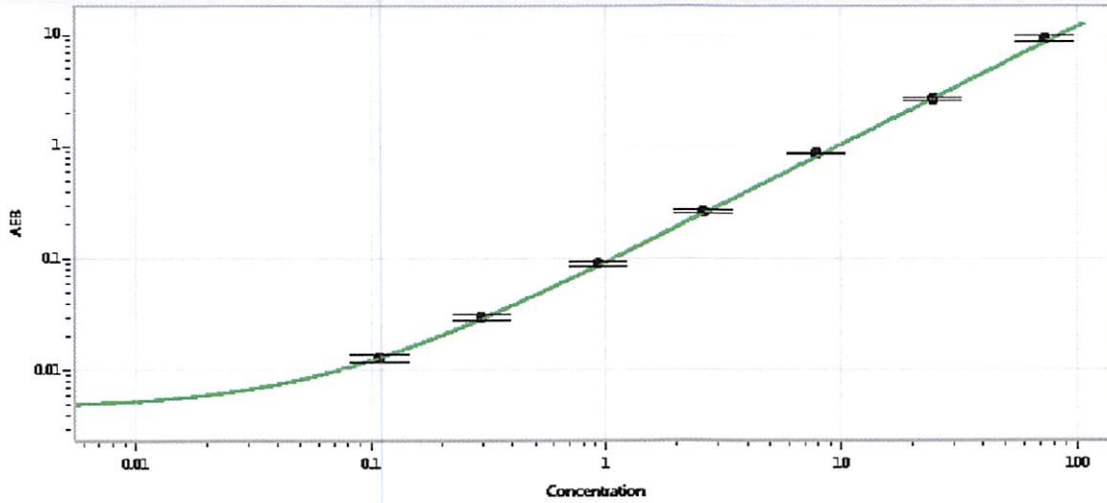
Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Sivasankari Murugar QC Supervisor  
Name Title

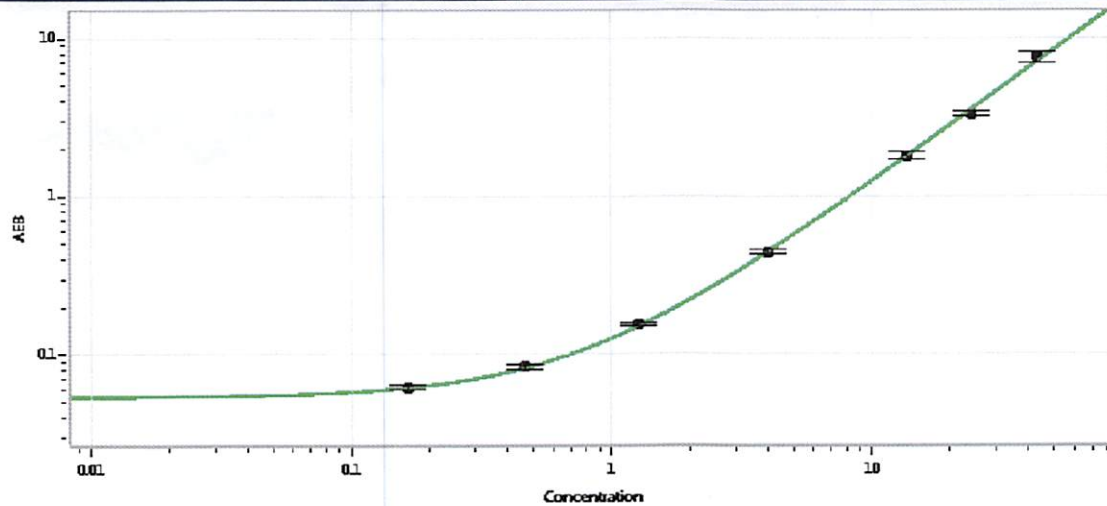
*[Signature]* 20 Nov 2018  
Signature/Date

**Tau Calibration Curve**



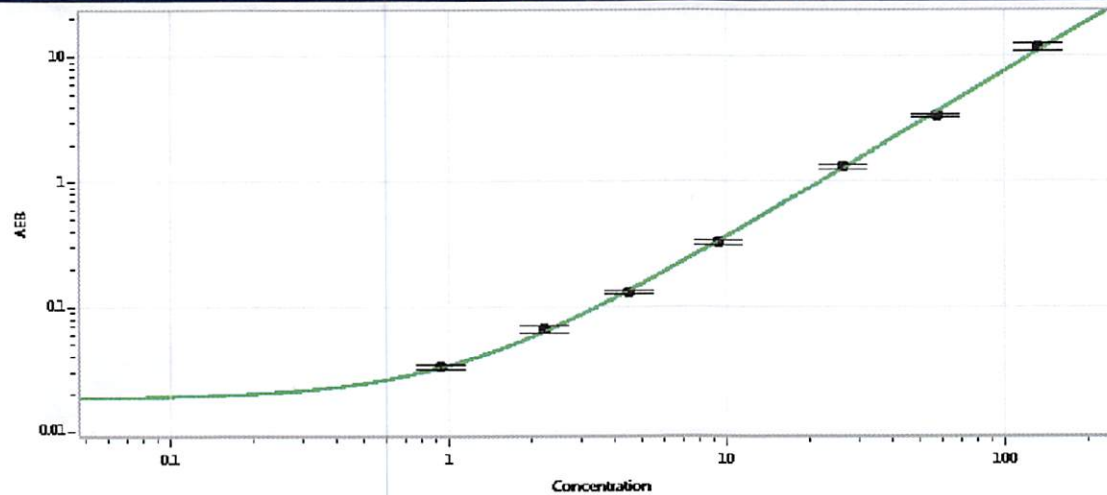
Calibrator Levels (pg/mL)	
A	0.000
B	0.109
C	0.294
D	0.919
E	2.59
F	7.88
G	24.7
H	73.5

**Abeta 42 Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.462
D	1.27
E	4.03
F	13.7
G	24.2
H	43.1

**Abeta 40 Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.938
C	2.22
D	4.48
E	9.43
F	26.3
G	57.3
H	133