

Product Number:	101995
Lot Number:	501169
Expiration:	26-Aug-2018
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	80660
Detector Reagent	101996	80750
SBG Reagent (HD-1)	102000	80571
Sample Diluent	102002	80572
Calibrators	102004	73560
Controls	102005-102006	80172
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	80172	2.19	1.62-2.43
Control 2	80172	94.2	70.3-106
Panel 1	MF-010818-01	1.15	0.286-2.01
Panel 2	MF-010818-02	189	156-233
Panel 3	MF-010818-03	3.76	1.23-6.14

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80172	5.64	4.70-7.05
Control 2	80172	134	105-158
Panel 1	MF-010818-01	1.12	0.539-3.91
Panel 2	MF-010818-02	8.55	6.97-10.8
Panel 3	MF-010818-03	120	96.8-145

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80172	17.2	13.7-20.6
Control 2	80172	251	193-289
Panel 1	MF-010818-01	46.1	26.1-63.3
Panel 2	MF-010818-02	145	105-203
Panel 3	MF-010818-03	95.6	70.1-125

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

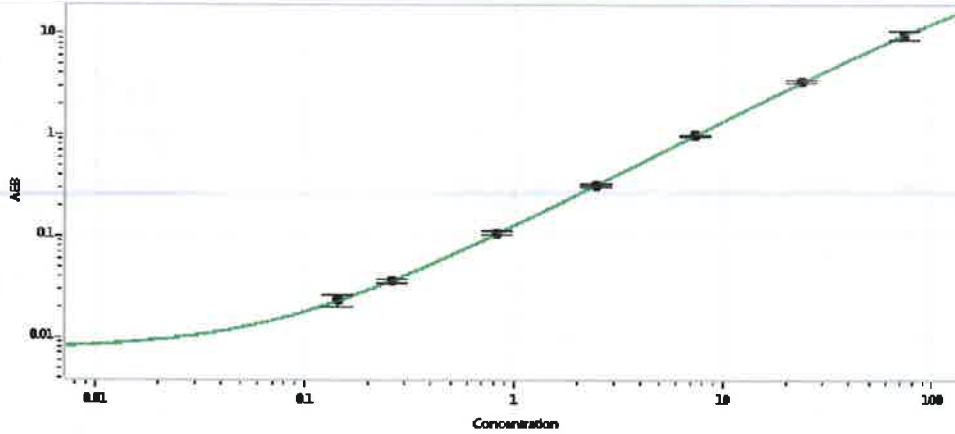
Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
21 MAR 2018

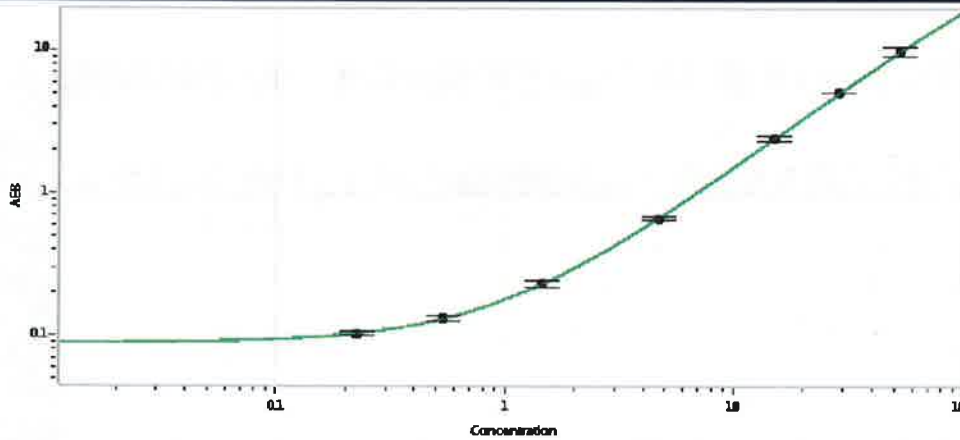
**Tau Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.143
C	0.262
D	0.828
E	2.47
F	7.37
G	23.8
H	73.1

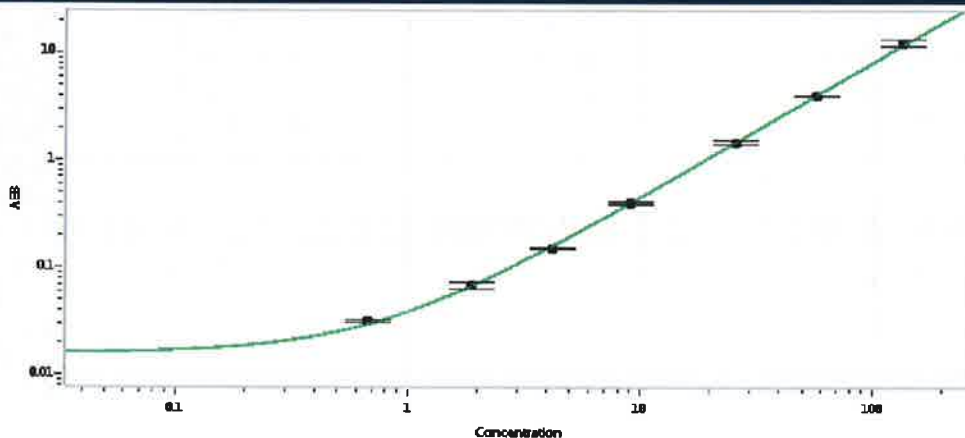
**Abeta 42**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.225
C	0.537
D	1.44
E	4.67
F	15.1
G	28.8
H	53.1

**Abeta 40**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.673
C	1.90
D	4.22
E	9.06
F	25.9
G	57.6
H	136