

Product Number:	101995
Lot Number:	501208
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	102000	80571
Sample Diluent	102002	80572
Calibrators	102004	80850/80941
Controls	102005-102006	80851
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 SRX platform - refer to the SR-X Kit Instructions (KI-0083).

**Release Materials**

**Tau**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80851	2.19	1.48-2.22
Control 2	80851	83.6	67.5-101
Panel 1	MF-010818-01	1.04	0.286-2.01
Panel 2	MF-010818-02	191	156-233
Panel 3	MF-010818-03	3.59	1.23-6.14

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80851	4.65	3.76-5.99
Control 2	80851	101	85.5-132
Panel 1	MF-010818-01	1.55	0.539-3.91
Panel 2	MF-010818-02	9.12	6.97-10.8
Panel 3	MF-010818-03	128	96.8-145

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80851	20.7	16.3-24.5
Control 2	80851	276	234-351
Panel 1	MF-010818-01	46.8	26.1-63.3
Panel 2	MF-010818-02	150	105-203
Panel 3	MF-010818-03	99.4	70.1-125

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

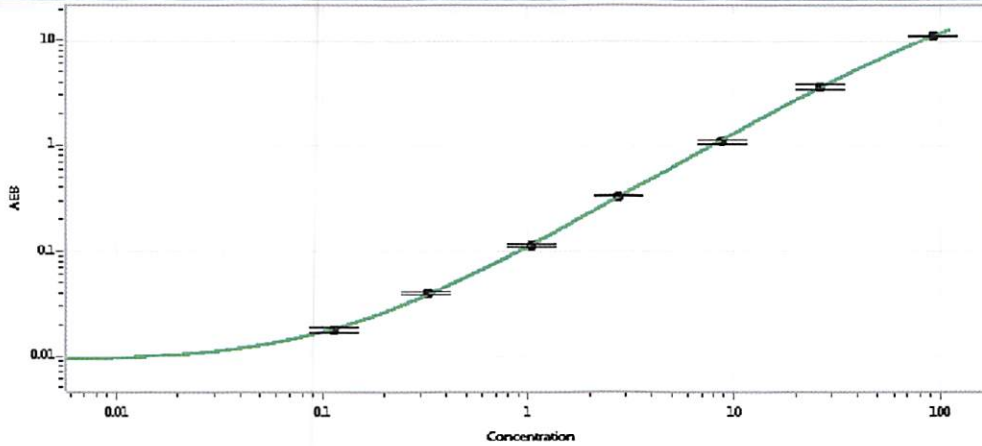
Review/Approval

Sivasankari Murugan QC Supervisor  
Name Title

*[Signature]* 18 APR 2018  
Signature/Date

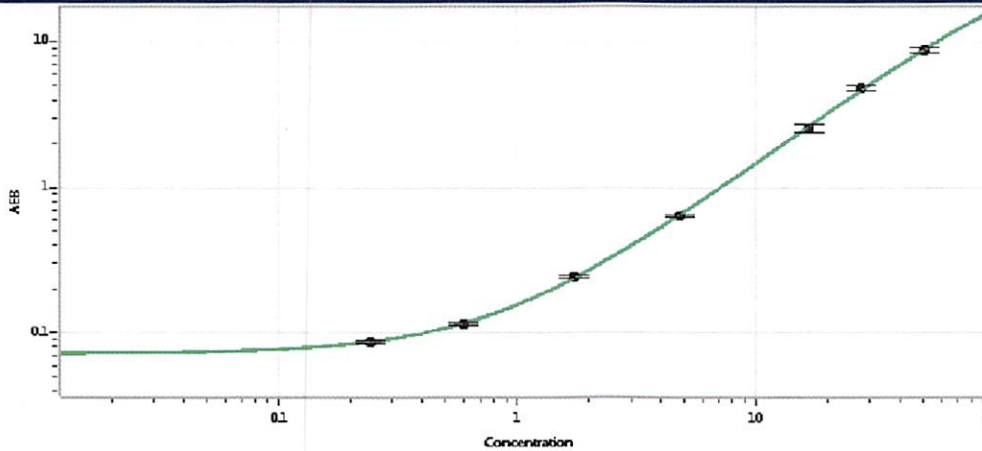


**Tau Calibration Curve**



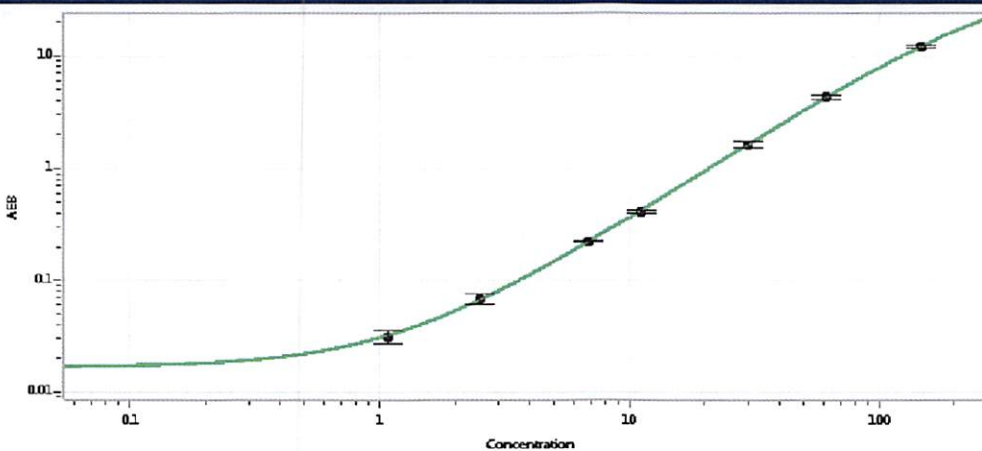
Calibrator Levels (pg/mL)	
A	0.000
B	0.114
C	0.324
D	1.04
E	2.75
F	8.78
G	26.3
H	92.4

**Abeta 42**



Calibrator Levels (pg/mL)	
A	0.000
B	0.242
C	0.594
D	1.73
E	4.84
F	16.7
G	27.6
H	51.1

**Abeta 40**



Calibrator Levels (pg/mL)	
A	0.000
B	1.08
C	2.52
D	6.81
E	11.1
F	30.1
G	62.0
H	147