

Product Number:	101995
Lot Number:	501427
Expiration:	21-Mar-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	832310
Detector Reagent	101996	832311
SBG Reagent	102000	826410
Sample Diluent	102002	826501
Calibrators	103221	820001
Controls	102005-102006	80172
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ 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	2.47	1.63-2.49
Control 2	80172	96.5	70.8-108
Panel 1	MF-062018-04	1.74	1.29-2.00
Panel 2	MF-010818-02	173	156-234
Panel 3	MF-010818-03	4.37	1.30-5.98

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80172	6.10	4.74-7.11
Control 2	80172	125	106-159
Panel 1	MF-062018-04	6.34	5.28-7.93
Panel 2	MF-010818-02	8.74	7.07-10.7
Panel 3	MF-010818-03	108	96.1-150

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80172	17.7	13.8-20.7
Control 2	80172	247	194-291
Panel 1	MF-062018-04	133	106-159
Panel 2	MF-010818-02	151	107-200
Panel 3	MF-010818-03	102	71.5-124

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

Review/Approval

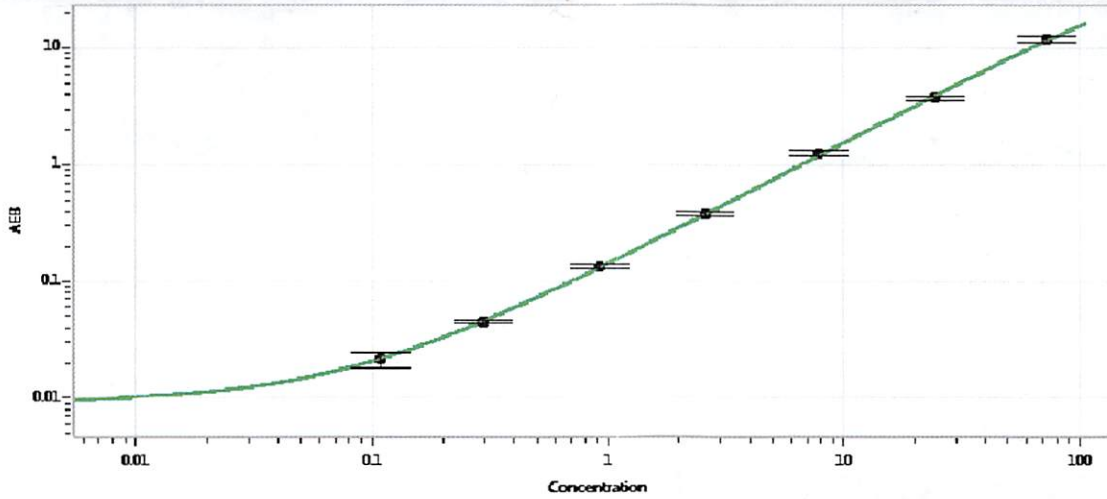
Sivasankari Murugan QC Supervisor

Name

Title

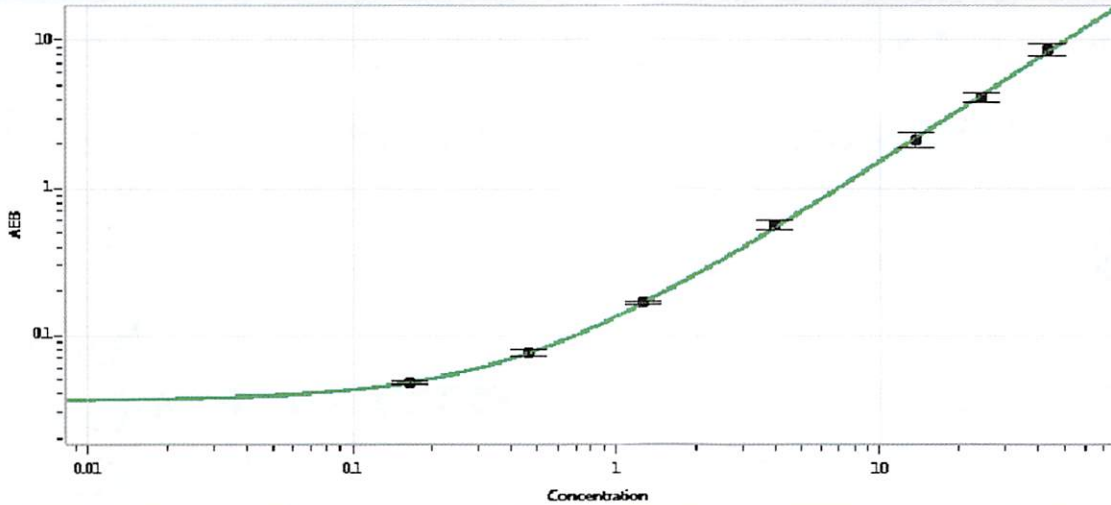
[Signature] 03DEC2018
Signature/Date

Tau Calibration Curve



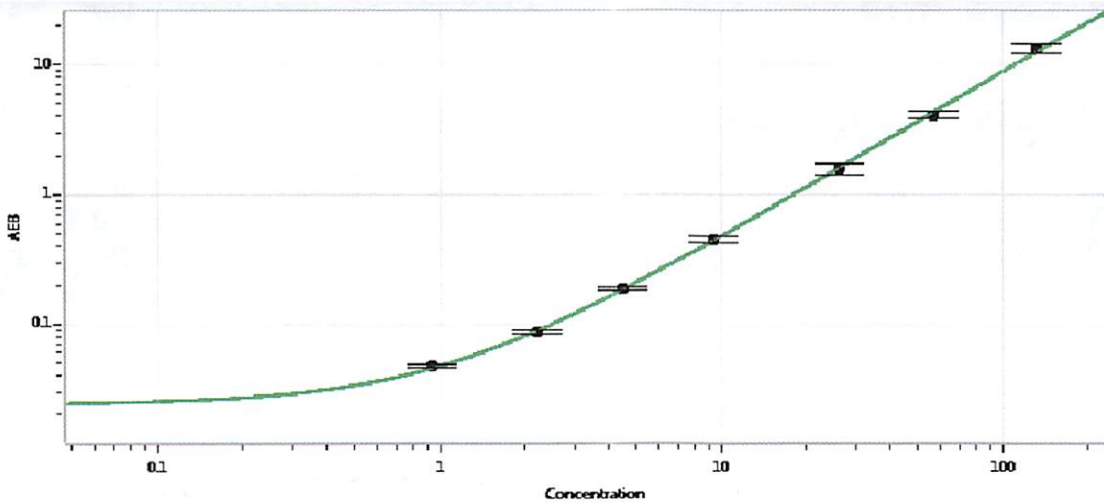
Calibrator Levels (pg/mL)	
A	0.000
B	0.109
C	0.294
D	0.919
E	2.59
F	7.88
G	24.7
H	73.5

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