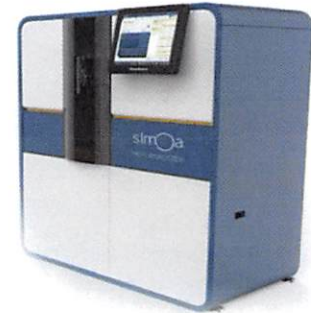


Product Number:	101995
Lot Number:	501229
Expiration:	26-Aug-2018
Platform(s):	HD-1 & SR-X



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	81142
Calibrators	102004	80850/80941
Controls	102005-102006	80851
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 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	1.94	1.49-2.23
Control 2	80851	88.6	68.0-102
Panel 1	MF-010818-01	1.05	0.317-1.96
Panel 2	MF-010818-02	199	156-234
Panel 3	MF-010818-03	3.29	1.30-5.98

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80851	4.27	3.60-5.97
Control 2	80851	104	85.9-131
Panel 1	MF-010818-01	1.36	0.379-3.87
Panel 2	MF-010818-02	8.98	7.07-10.7
Panel 3	MF-010818-03	140	96.1-150

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80851	20.0	16.3-24.4
Control 2	80851	292	234-351
Panel 1	MF-010818-01	46.9	27.2-62.6
Panel 2	MF-010818-02	149	107-200
Panel 3	MF-010818-03	103	71.5-124

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

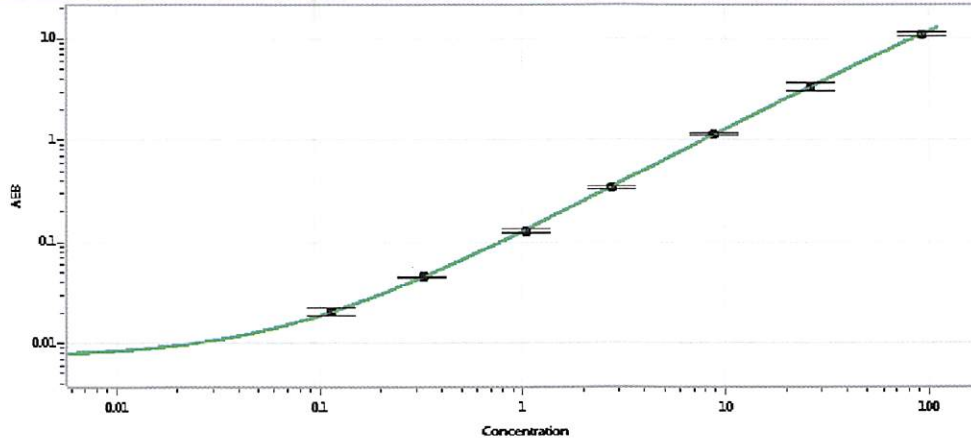
Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date 30 APR 2018

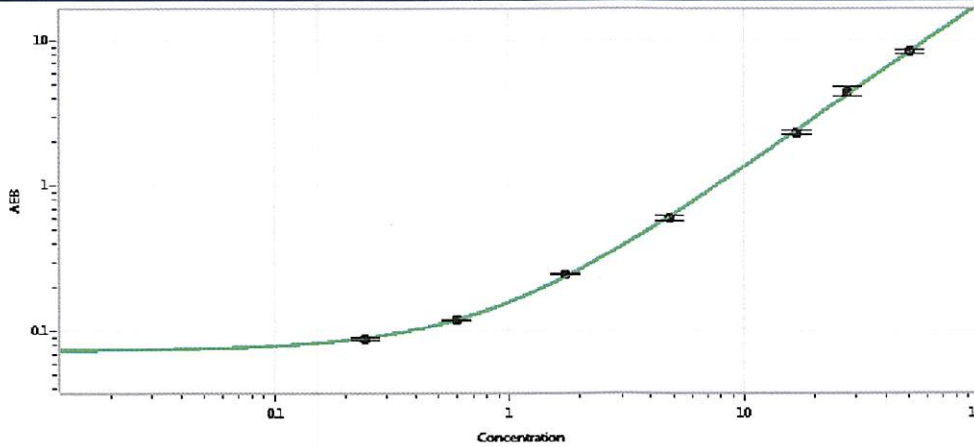
Tau Calibration Curve



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

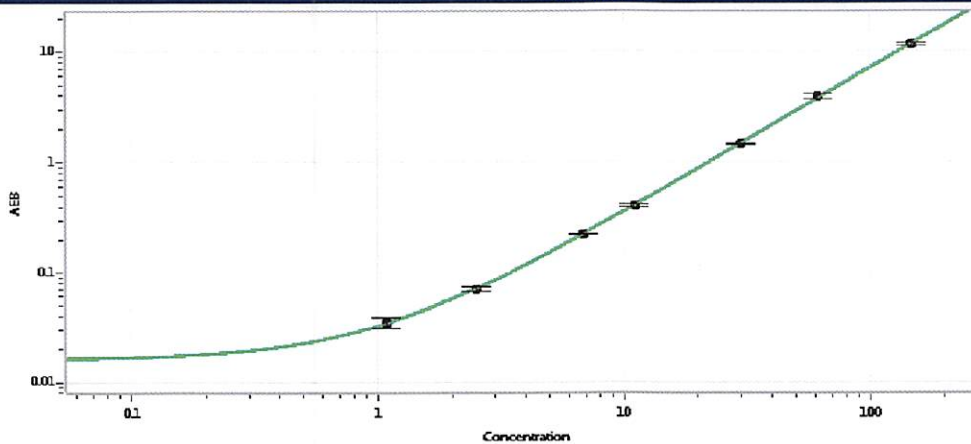
Abeta 42



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