

Product Number:	101876
Lot Number:	501238
Expiration:	2-Aug-2018
Platform(s):	HD-1



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101879	81241
Detector Reagent	101877	81242
SBG Reagent	101881	81014
Sample Diluent	101882	81080
Calibrators	101884	62150
Controls	101885-101886	62151
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials

Tau


	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	62151	1.88	0.961-2.88
Control 2	62151	108	59.9-180
Panel 2	MF-010818-02	187	138-206
Panel 3	MF-010818-03	3.27	2.78-4.17

Abeta 42

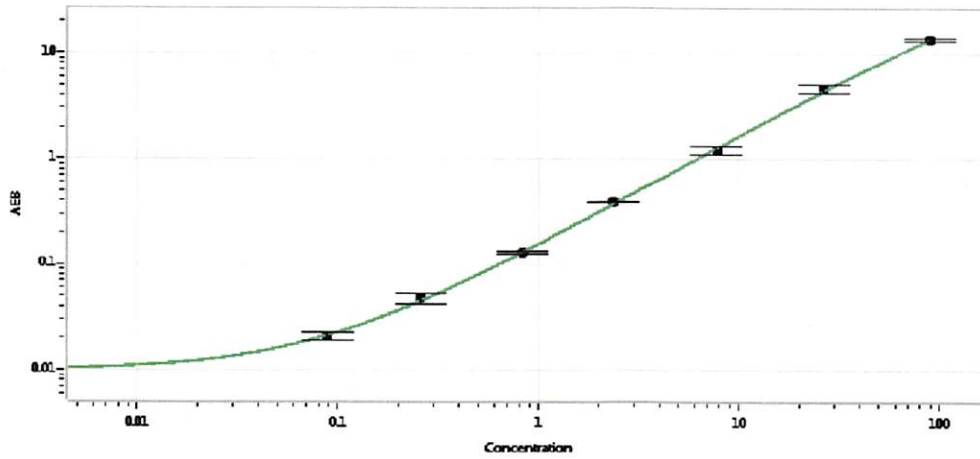
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	62151	0.588	0.391-1.16
Control 2	62151	45.9	42.9-64.3
Panel 2	MF-010818-02	14.2	11.0-19.3
Panel 3	MF-010818-03	161	111-263

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

James Drummey
Senior Director, Manufacturing Operations
On behalf of:

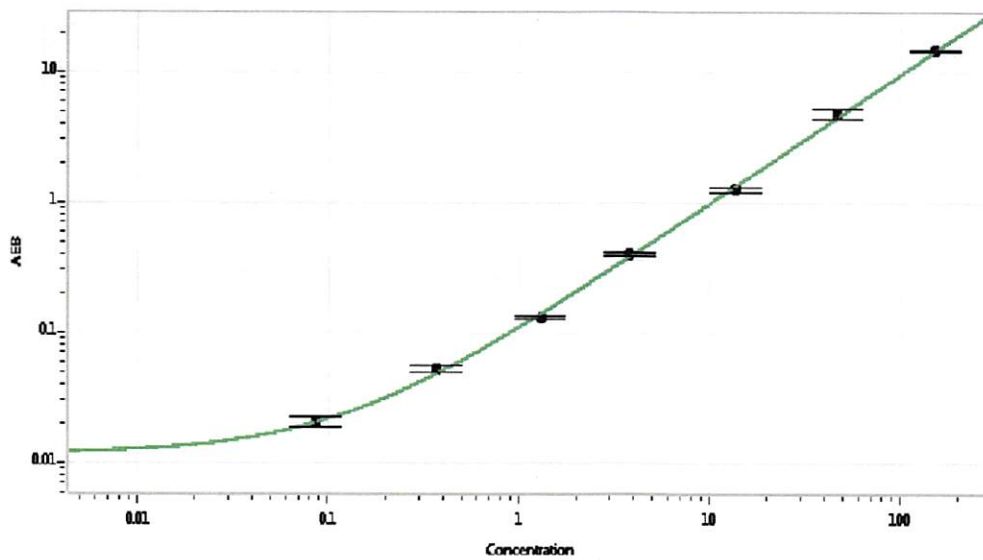
Review/Approval	Sivasankari Murugan	QC Supervisor		May 17, 2018
	Name	Title		

Tau Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.089
C	0.260
D	0.831
E	2.37
F	7.75
G	26.2
H	88.7

Abeta 42



Calibrator Levels (pg/mL)	
A	0.000
B	0.086
C	0.373
D	1.30
E	3.79
F	13.6
G	46.8
H	152