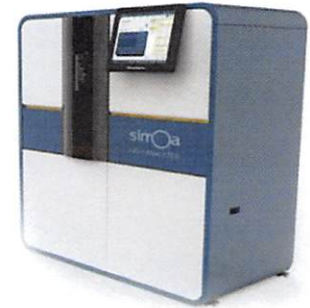


Product Number:	101552
Lot Number:	501132
Expiration:	4-Nov-2018
Platform(s):	HD-1 and SR-X



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

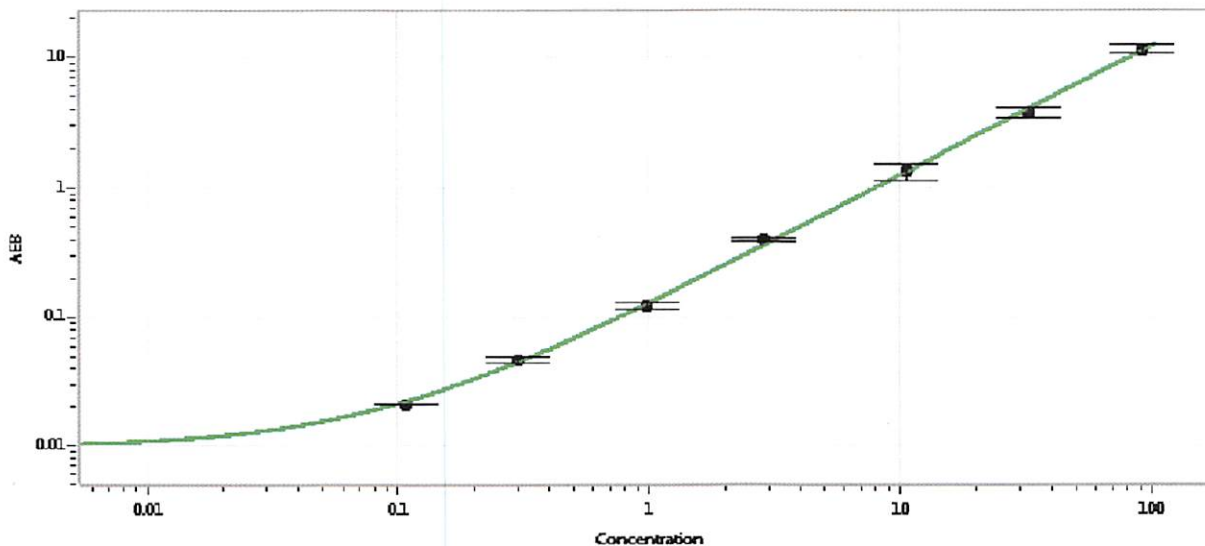
Component	Part Number	Lot Number
Bead Reagent	101554	80380
Detector Reagent	101553	80350
SBG Reagent (HD-1)†	101555	80351
Sample Diluent	101556	80451
Calibrators	101559	80400
Controls	101560-101561	80401
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

† Specific for use on the SRX platform - refer to the Simoa™ Human Total Tau Advantage SRX Kit Instructions (KI-0034).

Release Materials				Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	80401	1.07	0.860-1.37	
Control 2	80401	57.6	52.1-78.1	
Panel 1	MF-013017-02	0.710	0.251-1.23	
Panel 2	MF-050317-13	1.14	0.326-2.04	
Panel 3	MF-050417-01	64.5	38.2-115	
Panel 4	MF-050417-02	213	99.8-299	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.107
C	0.301
D	0.982
E	2.84
F	10.6
G	32.3
H	92.3

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
26 FEB 2018
Signature/Date