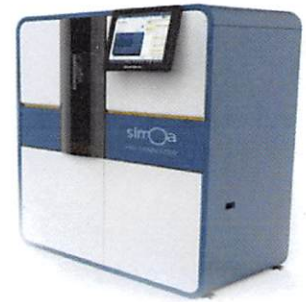


Product Number:	101552
Lot Number:	501472
Expiration:	5-Jul-2019
Platform(s):	HD-1 and SR-X



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

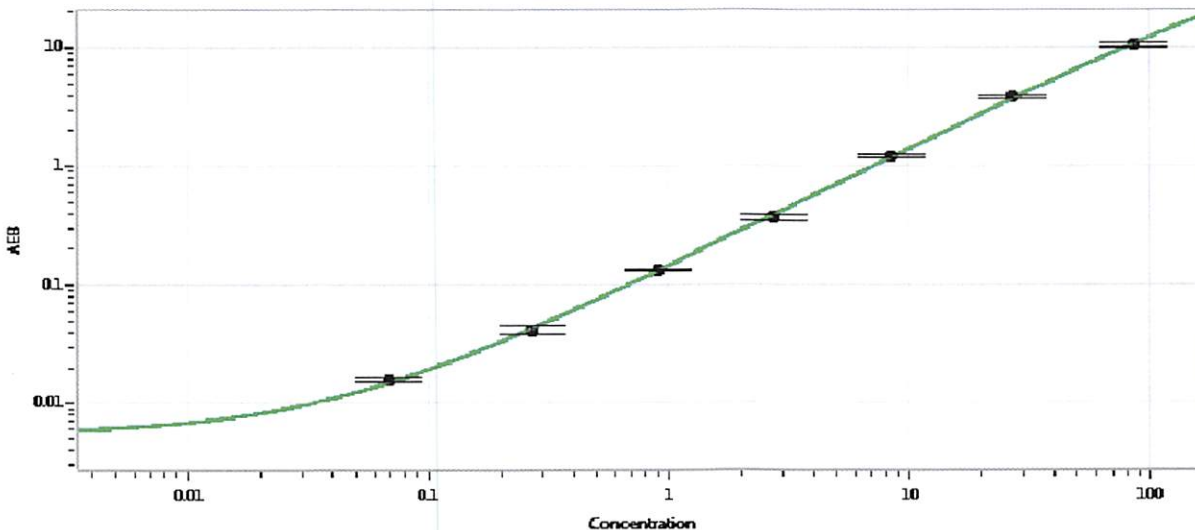
Component	Part Number	Lot Number
Bead Reagent	101554	829203
Detector Reagent	101553	829012
SBG Reagent	101555	828201
Sample Diluent	101556	827802
Calibrators	101559	819901
Controls	101560-101561	72650
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Human Total Tau Advantage SR-X 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	72650	1.16	0.532-1.60	
Control 2	72650	65.5	27.4-82.1	
Panel 1	MF-013017-02	0.887	0.251-1.23	
Panel 2	MF-050317-13	1.42	0.326-2.04	
Panel 3	MF-050417-01	82.2	38.2-115	
Panel 4	MF-050417-02	216	99.8-299	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.270
D	0.903
E	2.76
F	8.48
G	26.8
H	85.4

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date