

Product Number:	101552
Lot Number:	501245
Expiration:	30-Jan-2019
Platform(s):	HD-1 and SR-X



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

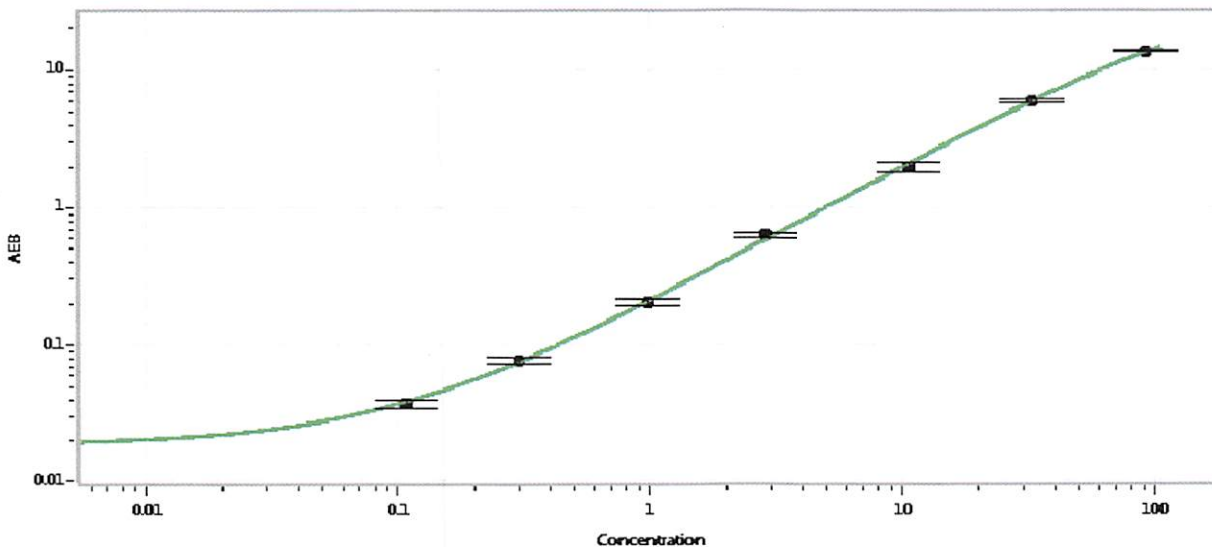
Component	Part Number	Lot Number
Bead Reagent	101554	81272
Detector Reagent	101553	81271
SBG Reagent	101555	81213
Sample Diluent	101556	81210
Calibrators	101559	80400
Controls	101560-101561	80401
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<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 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	1.10	0.787-1.40	
Control 2	80401	65.0	51.7-77.6	
Panel 1	MF-013017-02	0.722	0.251-1.23	
Panel 2	MF-050317-13	0.912	0.326-2.04	
Panel 3	MF-050417-01	83.1	38.2-115	
Panel 4	MF-050417-02	217	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]*  
Signature/Date