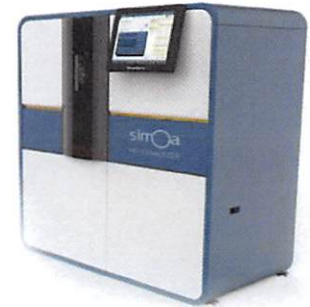


Product Number:	101552
Lot Number:	501341
Expiration:	26-Mar-2019
Platform(s):	HD-1 and SR-X



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

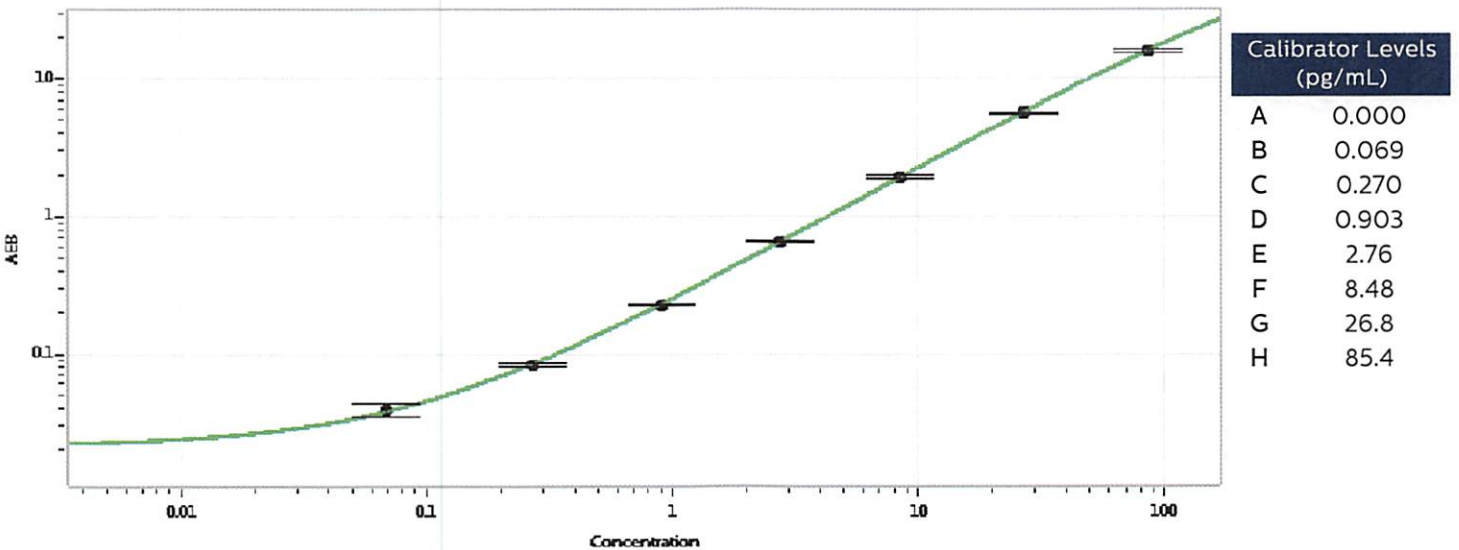
Component	Part Number	Lot Number
Bead Reagent	101554	81784
Detector Reagent	101553	818601
SBG Reagent	101555	81780
Sample Diluent	101556	81770
Calibrators	101559	819901
Controls	101560-101561	72650
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<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	72560	1.04	0.532-1.60	
Control 2	72560	51.7	27.4-82.1	
Panel 1	MF-013017-02	0.527	0.251-1.23	
Panel 2	MF-050317-13	0.894	0.326-2.04	
Panel 3	MF-050417-01	55.4	38.2-115	
Panel 4	MF-050417-02	172	99.8-299	

Calibration Curve



Review/Approval: Sivasankari Murugan (Name), QC Supervisor (Title), [Signature] 11 AUG 2018 (Signature/Date)