

Product Number:	101552
Lot Number:	501576
Expiration:	18-Sep-2019
Platform(s):	HD-1 and SR-X



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

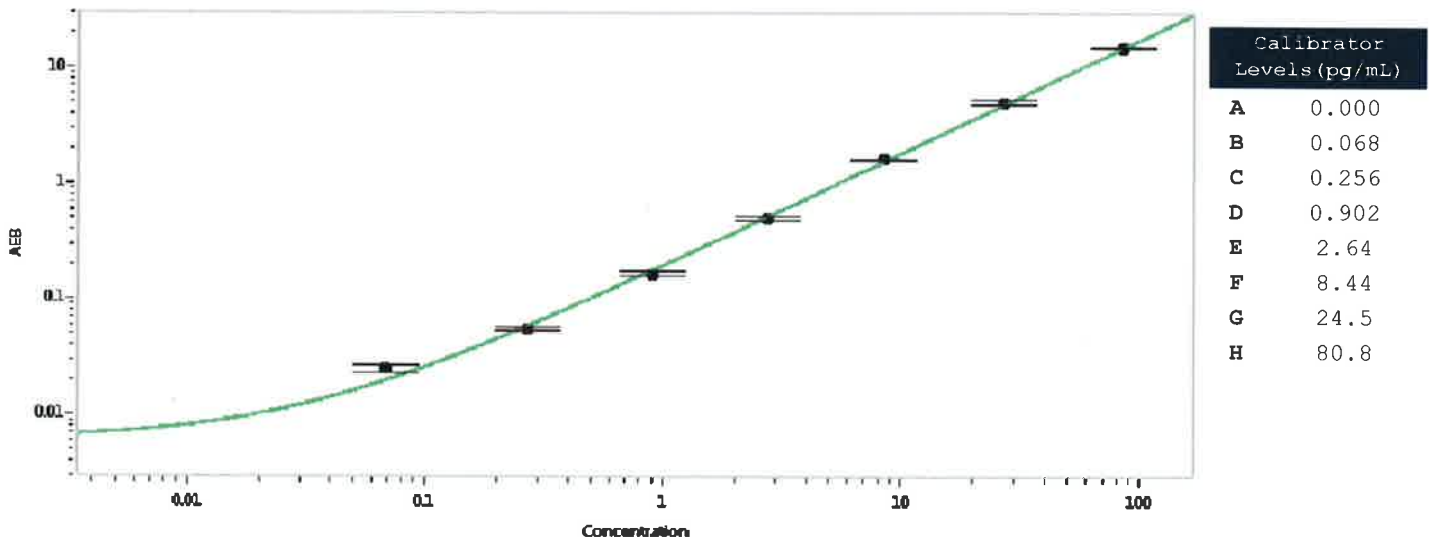
Component	Part Number	Lot Number
Bead Reagent	101554	836101
Detector Reagent	101553	835204
SBG Reagent	101555	835404
Sample Diluent	101556	835205
Calibrators	101559	900801
Controls	101560-101561	900905
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ 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	900905	1.25	1.02-1.54	
Control 2	900905	58.9	47.6-71.4	
Panel 1	MF-013017-02	0.828	0.251-1.23	
Panel 2	MF-050317-13	1.19	0.326-2.04	
Panel 3	MF-050417-01	71.2	38.2-115	
Panel 4	MF-050417-02	175	99.8-299	

Calibration Curve



Review/Approval: Patrizia Stadler Associate Director, QC
 Name: Patrizia Stadler Title: QC Signature/Date: P. Stadler 15 Jan 2019