

Product Number:	101552
Lot Number:	501834
Expiration:	23-Mar-2020
Platform(s):	HD-1 and SR-X



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

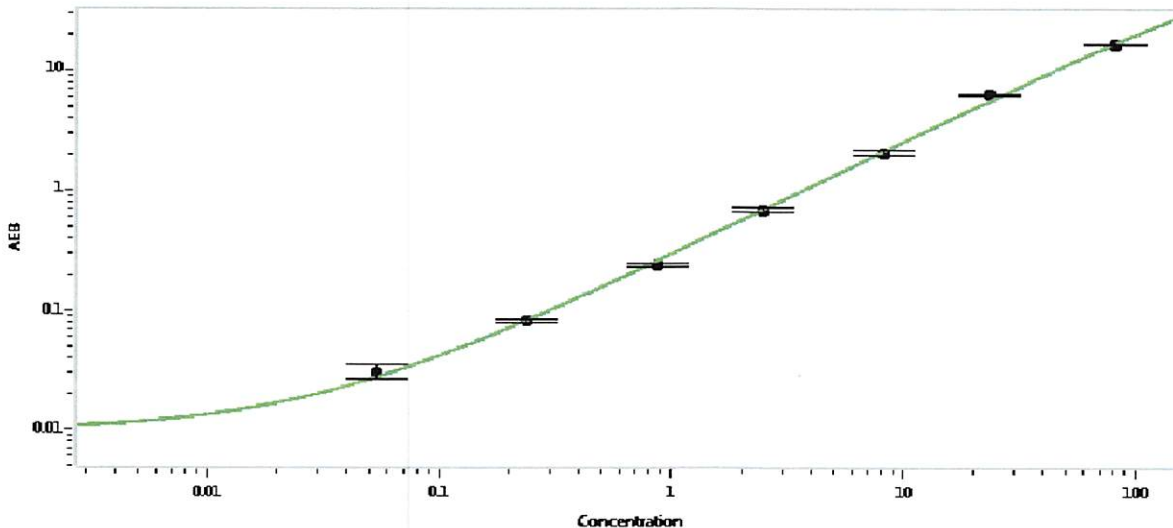
Component	Part Number	Lot Number
Bead Reagent	101554	908805
Detector Reagent	101553	908509
SBG Reagent	101555	908508
Sample Diluent	101556	908204
Calibrators	101559	912902
Controls	101560-101561	912903
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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	912903	0.766	0.680-1.28	
Control 2	912903	53.5	42.5-63.8	
Panel 1	MF-013017-02	0.942	0.251-1.23	
Panel 2	MF-050317-13	1.36	0.326-2.04	
Panel 3	MF-050417-01	69.3	38.2-115	
Panel 4	MF-050417-02	165	99.8-299	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.054
C	0.238
D	0.865
E	2.46
F	8.11
G	23.3
H	81.2

Review/Approval

Junpil Park*
Name

QC Analyst II
Title

[Signature] 23 May 2019
Signature/Date