

Product Number:	101552
Lot Number:	501843
Expiration:	26-Mar-2020
Platform(s):	HD-1 and SR-X



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

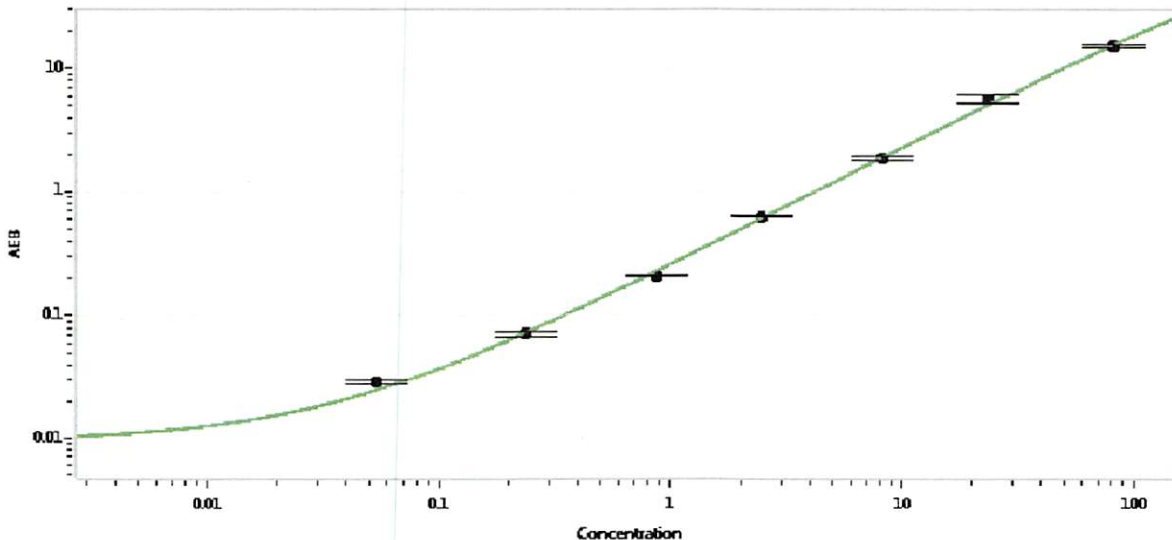
Component	Part Number	Lot Number
Bead Reagent	101554	908805
Detector Reagent	101553	908509
SBG Reagent	101555	908508
Sample Diluent	101556	908814
Calibrators	101559	912902
Controls	101560-101561	912903
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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	0.833	0.680-1.28	
Control 2	912903	55.7	42.5-63.8	
Panel 1	MF-013017-02	0.727	0.251-1.23	
Panel 2	MF-050317-13	1.03	0.326-2.04	
Panel 3	MF-050417-01	70.3	38.2-115	
Panel 4	MF-050417-02	173	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: Patrizia Stadler Associate Director, QC
 Name: Patrizia Stadler Title: QC Signature/Date: P. Stadler 21 May 2019