

<b>Product Number:</b>	101552
<b>Lot Number:</b>	501764
<b>Expiration:</b>	18-Dec-2019
<b>Platform(s):</b>	HD-1 and SR-X



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

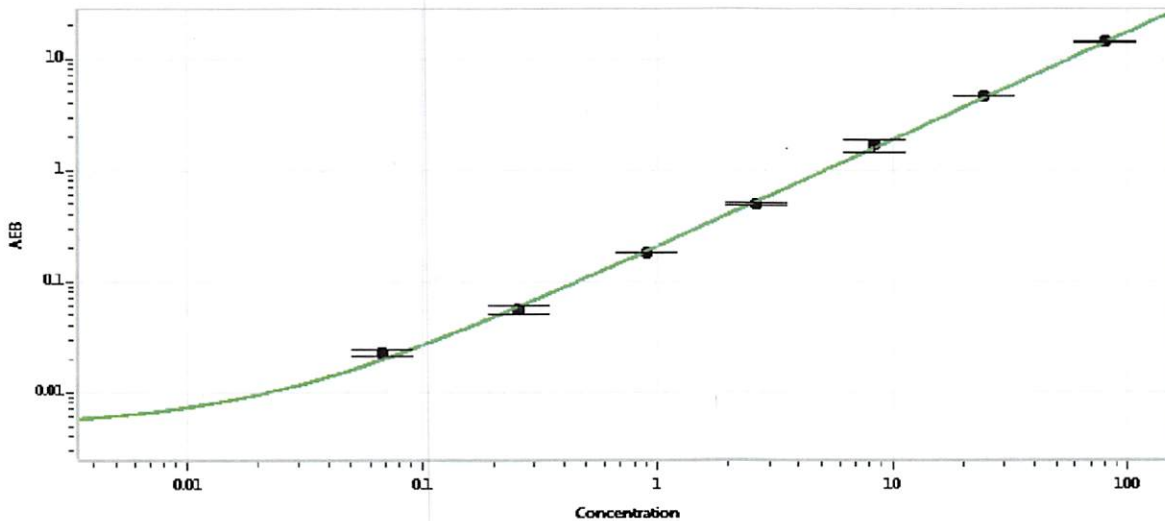
Component	Part Number	Lot Number
Bead Reagent	101554	836101
Detector Reagent	101553	835204
SBG Reagent	101555	835404
Sample Diluent	101556	908201
Calibrators	101559	900801
Controls	101560-101561	900905
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<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	900905	1.17	1.02-1.54	
Control 2	900905	60.6	47.6-71.4	
Panel 1	MF-013017-02	0.837	0.251-1.23	
Panel 2	MF-050317-13	1.30	0.326-2.04	
Panel 3	MF-050417-01	69.0	38.2-115	
Panel 4	MF-050417-02	156	99.8-299	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.068
C	0.256
D	0.902
E	2.64
F	8.44
G	24.5
H	80.8

<b>Review/Approval</b>	Associate Director, QC Patrizia Stadler	P. Stadler	11 Apr 2019
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>