

Product Number:	101552
Lot Number:	501987
Expiration:	23-Mar-2020
Platform(s):	HD-1/HD-x & 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 <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

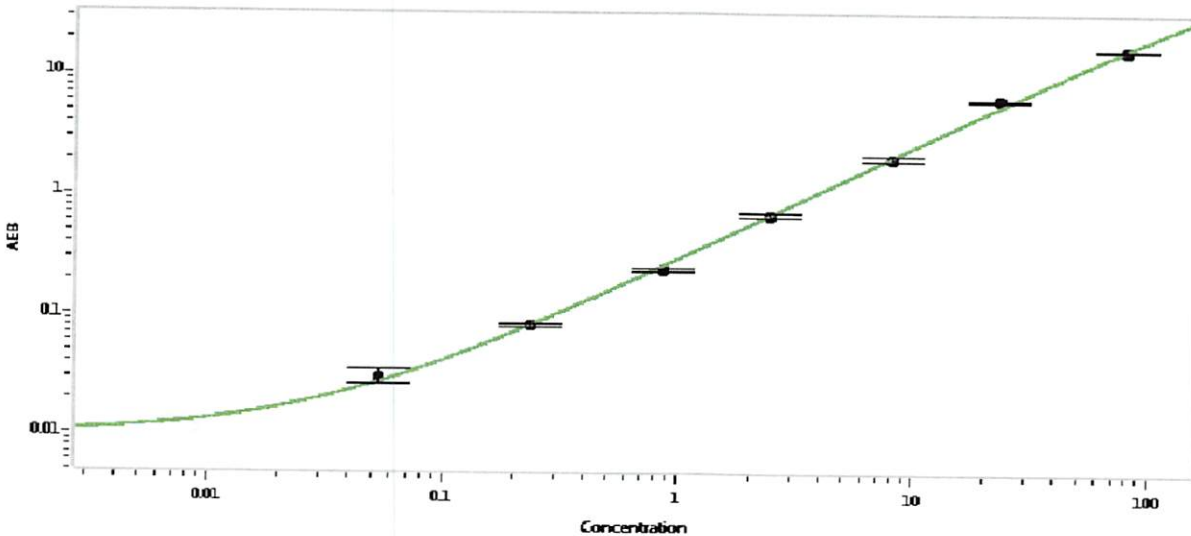
<sup>1</sup> RGP is not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Human Total Tau Advantage SR-X Kit Instructions (KI-0034).

Release Materials*			
	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

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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

Daniel Ritchie Senior QC Analyst

Name

Title

Signature/Date

*[Handwritten Signature]* / 22 Aug 2019

\*Signing on behalf of Patrizia Stadler