

Product Number:	101552
Lot Number:	502085
Expiration:	21-Oct-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	929407
Detector Reagent	101553	929508
SBG Reagent	101555	929507
Sample Diluent	101556	929506
Calibrators	101559	931701
Controls	101560-101561	919802
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

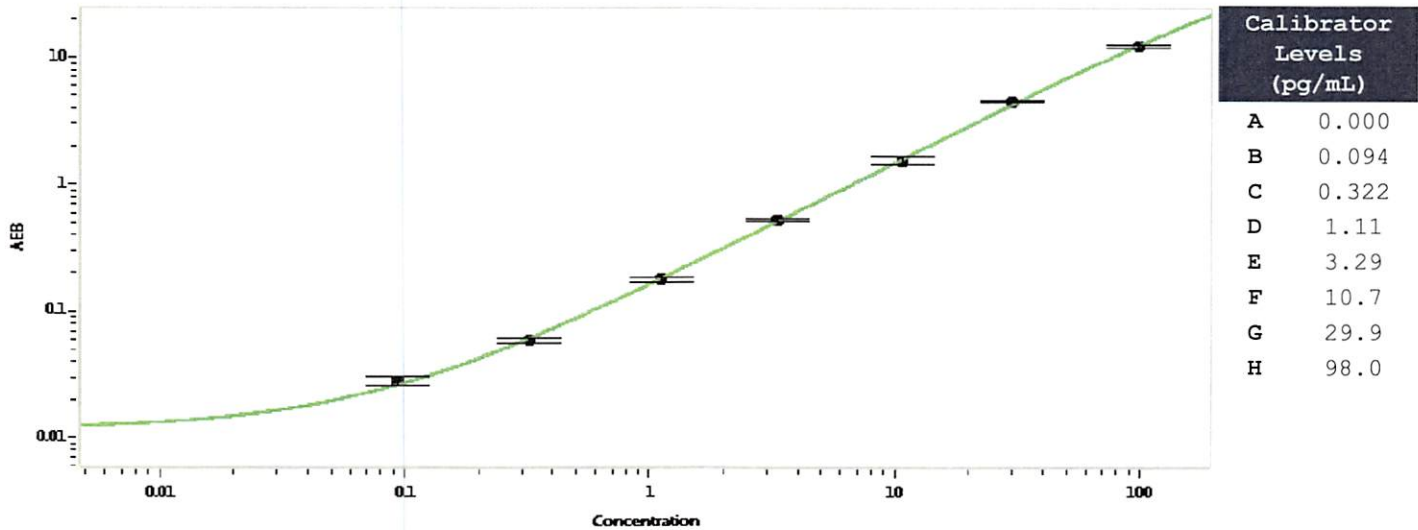
<sup>1</sup> RGP is not Kit Lot Specific

† 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	919802	0.724	0.721-1.08
Control 2	919802	40.3	35.0-52.5
Panel 1	916901	0.593	0.438-1.55
Panel 2	916902	1.19	0.822-2.11
Panel 3	916903	30.2	15.2-33.0
Panel 4	916904	92.4	61.7-98.2

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

Calibration Curve



Review/Approval

Patrizia Stadler Assoc. Director, QC  
Name Title

*P. Stadler* 21 NOV 2019  
Signature/Date