

Product Number:	102292
Lot Number:	502166
Expiration:	13-Sep-2020
Platform(s):	HD-1/HD-X and SR-X



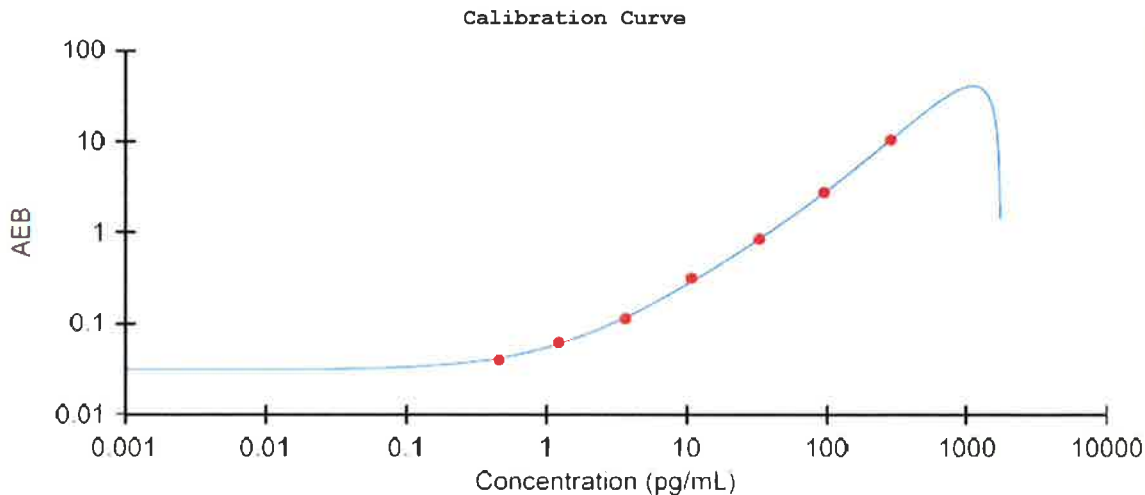
Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102280	930404
Detector Reagent	102282	930213
SBG Reagent	102284	934310
Sample Diluent	102286	930211
Calibrators	102289	825601
Controls	102290-102291	826405   825602
SBG Diluent (SRX) †	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™ pTau-231 SR-X Kit Instructions (KI-0110).

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	826405	14.2	8.14-18.5	
Control 2	825602	401	281-519	
Panel 1	823901	29.7	15.3-31.5	
Panel 2	823902	22.8	10.1-24.3	
Panel 3	823903	65.1	43.9-72.8	
Panel 4	823904	112	73.9-123	



Calibrator Levels (pg/mL)	
A	0.000
B	0.458
C	1.23
D	3.66
E	10.8
F	32.6
G	95.3
H	287

<b>Review/Approval</b>	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	16 Dec 2019
	Name	Title	Signature/Date	