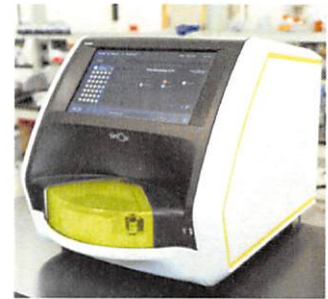


Product Number:	102292
Lot Number:	501996
Expiration:	13-Jun-2020
Platform(s):	HD-1/HD-X and SR-X



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

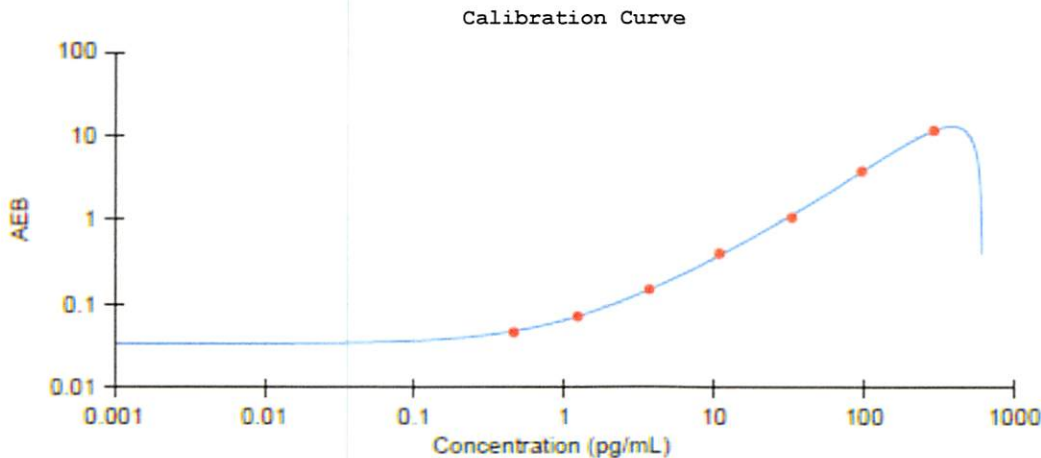
Component	Part Number	Lot Number
Bead Reagent	102280	923410
Detector Reagent	102282	923107
SBG Reagent	102284	923108
Sample Diluent	102286	923203
Calibrators	102289	825601
Controls	102290-102291	826405   825602
SBG Diluent (SRX) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<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			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	826405	14.8	7.83-17.7
Control 2	825602	387	282-501
Panel 1	823901	23.5	17.5-26.3
Panel 2	823902	15.7	11.9-19.4
Panel 3	823903	55.9	44.7-67.0
Panel 4	823904	89.3	75.1-113

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



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

Review/Approval	Patrizia Stadler Name	Assoc. Director, QC Title	<i>P Stadler</i> Signature/Date	30 Aug 2019 Date
-----------------	--------------------------	------------------------------	------------------------------------	---------------------