

Product Number:	102292
Lot Number:	501476
Expiration:	23-Apr-2019
Platform(s):	HD-1 & SR-X



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

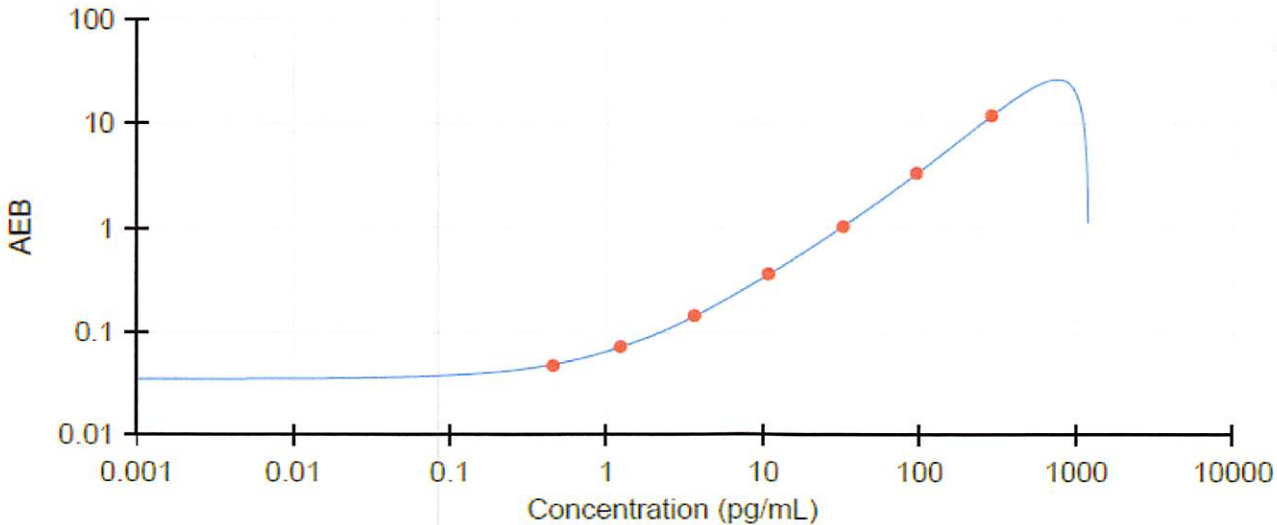
Component	Part Number	Lot Number
Bead Reagent	102280	829907
Detector Reagent	102282	829606
SBG Reagent	102284	829603
Sample Diluent	102286	829604
Reference Calibrators	102289	825601
Controls	102290-102291	826405 825602
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ 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	15.6	7.83-17.7	
Control 2	825602	375	282-501	
Panel 1	823901	24.1	17.5-26.3	
Panel 2	823902	15.5	11.9-19.4	
Panel 3	823903	60.8	44.7-67.0	
Panel 4	823904	101	75.1-113	

Calibration Curve



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

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 20DEC2018
Signature/Date