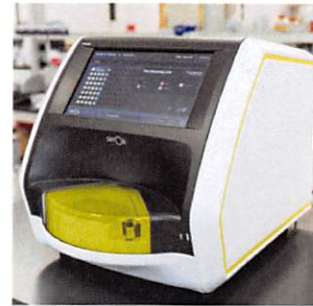


Product Number:	102292
Lot Number:	501677
Expiration:	23-Oct-2019
Platform(s):	HD-1 & SR-X



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

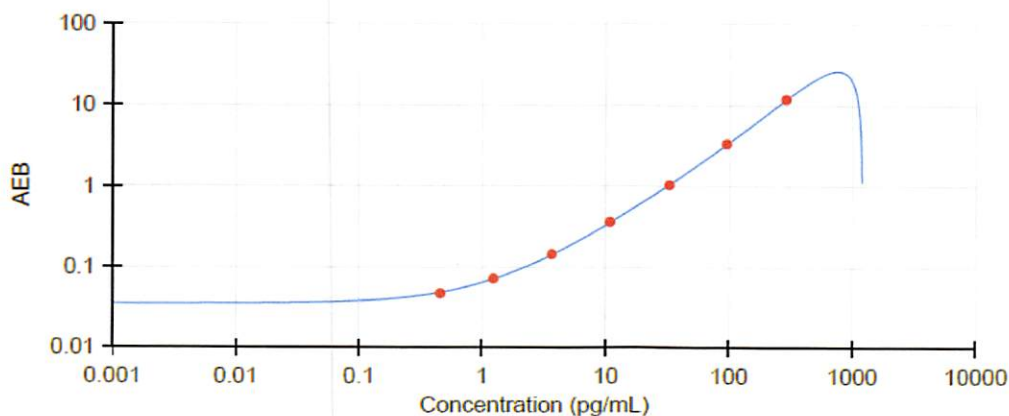
Component	Part Number	Lot Number
Bead Reagent	102280	829917
Detector Reagent	102282	829610
SBG Reagent	102284	829603
Sample Diluent	102286	829611
Calibrators	102289	825601
Controls	102290-102291	826405-825602
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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	823903	15.5	11.9-19.4	
Panel 3	823904	60.8	44.7-67.0	
Panel 3	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	287.0

Review/Approval

Patrizia Stadler

QC assoc director

Name

Title

Signature/Date

p. Stadler 27 Feb 2019