

Product Number:	102292
Lot Number:	502717
Expiration:	26-Jan-2022
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Reagent	102280	103007
Detector Reagent	102282	103603
SBG Reagent	102284	103609
Sample Diluent	102286	103608
Reference Calibrator	102289	014301
Control 1	102290	014302
Control 2	102291	014302
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

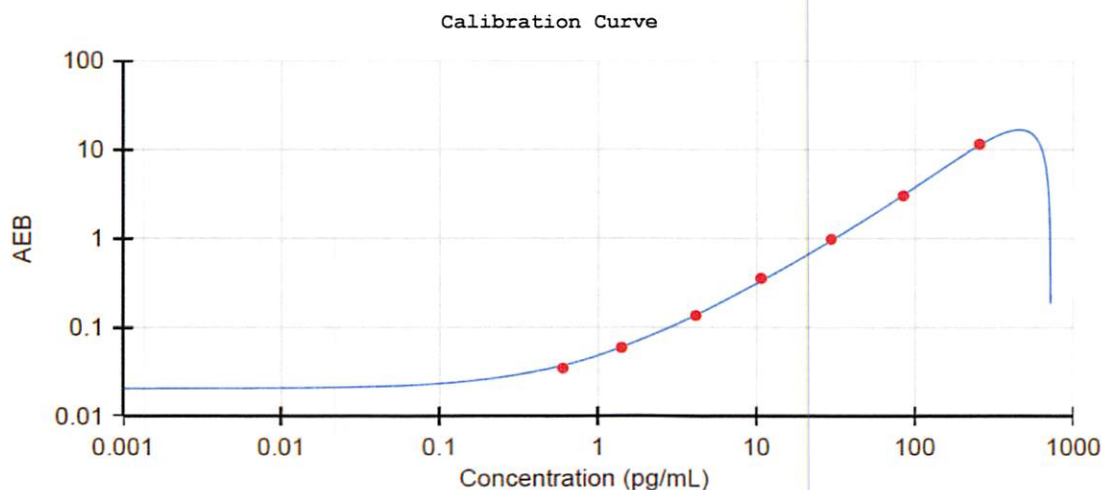
¹ Reagents are not Kit Lot Specific.

† Specific for use on the SR-X platform - refer to the Simoa™ pTau-231 SR-X Kit Instructions (KI-0110).



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

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	014302	14.5	10.2-15.2	
Control 2	014302	320	267-401	
Panel 1	111816	37.0	22.9-43.2	
Panel 2	111816	35.8	23.4-40.7	
Panel 3	111816	7.00	6.25-12.4	
Panel 4	111816	5.53	3.98-10.7	



Calibrator Levels (pg/mL)	
A	0.000
B	0.603
C	1.41
D	4.15
E	10.7
F	29.6
G	84.3
H	256

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 13 May 2021
Signature/Date