

Product Number:	102292
Lot Number:	503798
Expiration:	30-Jan-2024
Platform(s):	HD-1/HD-X and SR-X



Component	Part Number	Lot Number
Bead Reagent	102280	303007
Detector Reagent	102282	303104
SBG Reagent	102284	303105
Sample Diluent	102286	303205
Reference Calibrator	102289	306702
Control 1	102290	306703
Control 2	102291	306703
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

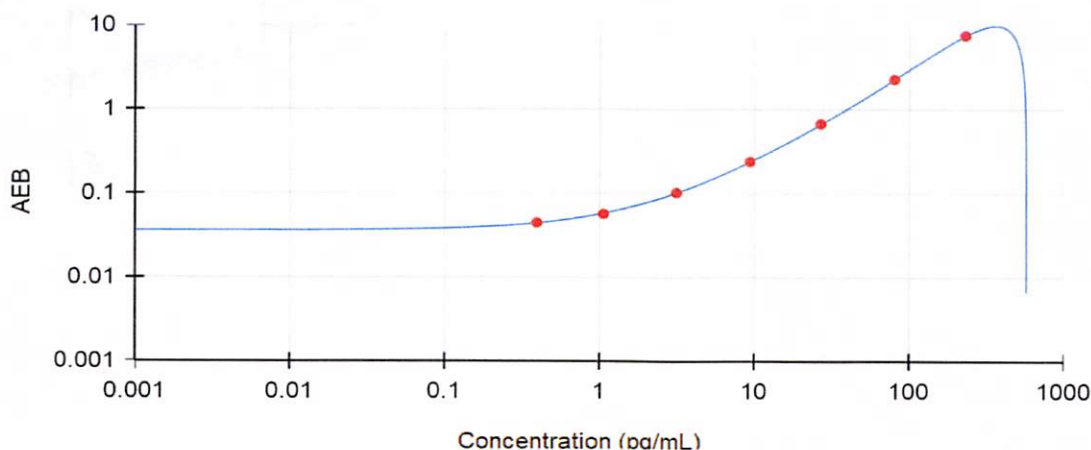
¹ Reagents are not Kit Lot Specific.

† 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 Simoa® SR-X Analyzer.

Release Materials				Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1	306703	14.7	10.8 - 16.2	
Control 2	306703	316	266 - 518	
Panel 3	313804	27.5	21.5 - 32.3	
Panel 4	313804	85.4	68 - 101	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.417
C	1.15
D	2.97
E	9.19
F	26.9
G	79.4
H	228

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 27 JUN 2023
Signature/Date