

Product Number:	102292
Lot Number:	503704
Expiration:	7-Sep-2023
Platform(s):	HD-1 & SR-X



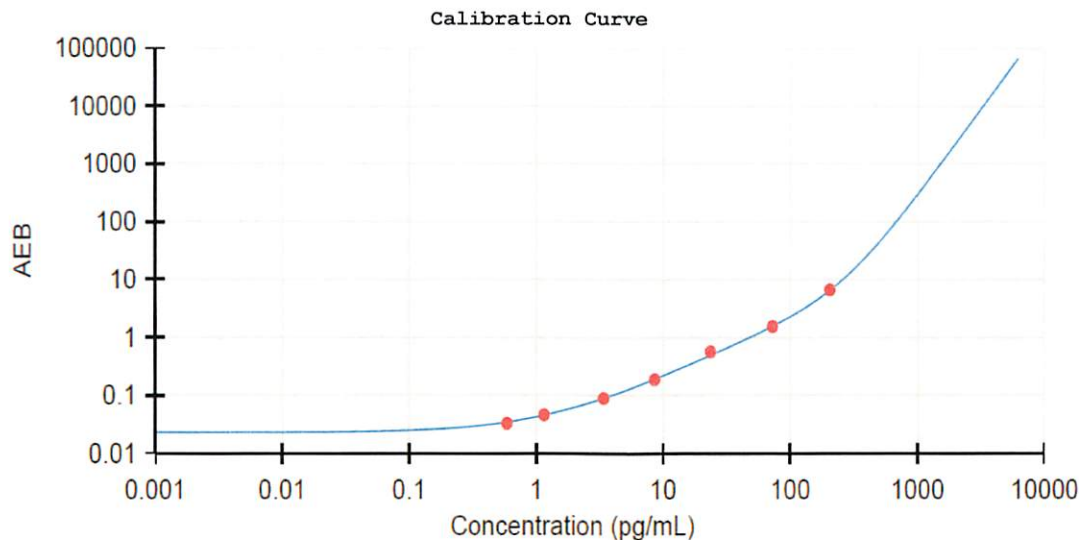
Data below represents results generated on the Simoa® SR-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102280	225703
Detector Reagent	102282	225704
SBG Reagent	102284	225705
Sample Diluent	102286	225706
Reference Calibrator	102289	218105
Control 1	102290	218106
Control 2	102291	218106
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)	Spec. Range (pg/mL)	
Control 1	218106	15.2	7.74 - 18.3	
Control 2	218106	362	178 - 410	
Panel 1	111816	37.9	23.7 - 44.3	
Panel 2	111816	37.2	22.0 - 42.7	
Panel 3	111816	6.38	2.31 - 17.0	
Panel 4	111816	6.35	3.69 - 10.3	



Calibrator Levels (pg/mL)	
A	0.000
B	0.589
C	1.15
D	3.39
E	8.59
F	23.5
G	72.3
H	205

Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Khillan / 10NOV2022
Signature/Date