

Product Number:	102292
Lot Number:	504212
Expiration Date:	19-Sep-25
Platform(s):	SR-X and HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	102280	410205
Detector Reagent	102282	410210
SBG Reagent	102284	410211
Sample Diluent	102286	410302
Calibrators	102289	326202
Control 1	102290	326405
Control 2	102291	326405
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

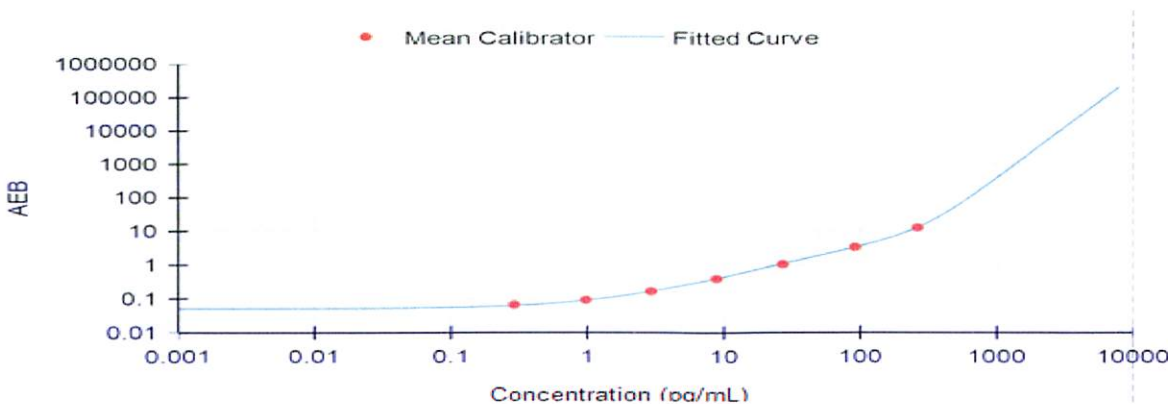
¹ Reagents are not Kit Lot Specific

† Specific to use on the SRX platform.
Refer to the pTau-231 Kit Instructions - SR-X (KI-0110).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	326405	13.7	10.4 - 16.8
Control 2	326405	436	242 - 488
Panel 1	N/A	N/A	N/A
Panel 2	N/A	N/A	N/A
Panel 3	313804	22.4	20.6 - 31.5
Panel 4	313804	74.8	66.2 - 99.3

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.290
C	0.975
D	2.93
E	8.79
F	27.0
G	91.0
H	262

Review/Approval

Cody Carlson

QC Manager

Name

Title

[Signature] 18 APR 2024

Signature/Date