

Product Number:	102292
Lot Number:	504213
Expiration Date:	16-Mar-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	326302
Detector Reagent	102282	326204
SBG Reagent	102284	325905
Sample Diluent	102286	326206
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

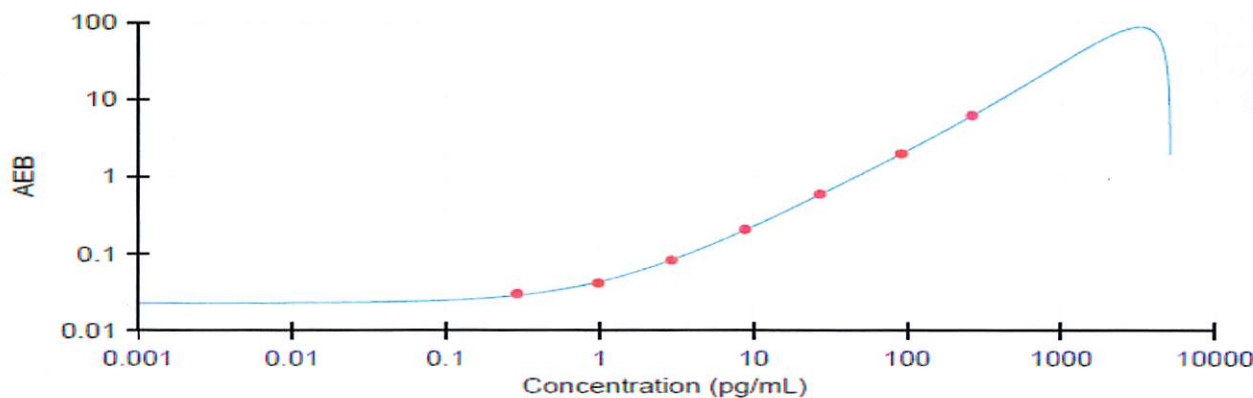
¹ Reagents are not Kit Lot Specific

† 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	10.7 - 16.6
Control 2	326405	361	273 - 410
Panel 3	313804	26.1	21.4 - 32.1
Panel 4	313804	83.6	67.5 - 101

* 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] 11 APR 2024
Signature/Date