

Product Number:	102292
Lot Number:	503935
Expiration:	5-Aug-2024
Platform(s):	HD-1/HD-X and SR-X



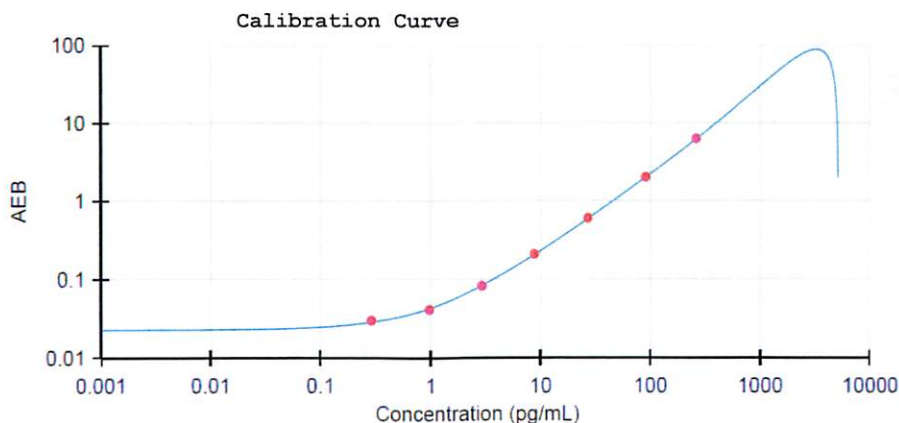
Component	Part Number	Lot Number
Bead Reagent	102280	326302
Detector Reagent	102282	326204
SBG Reagent	102284	325905
Sample Diluent	102286	326206
Reference Calibrator	102289	326202
Control 1	102290	326405
Control 2	102291	326405
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific.

† Specific 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	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	



Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 16 Oct 2023
Signature/Date