

Product Number:	102292
Lot Number:	503386
Expiration:	1-Apr-2023
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Reagent	102280	209105
Detector Reagent	102282	209201
SBG Reagent	102284	209202
Sample Diluent	102286	209405
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.

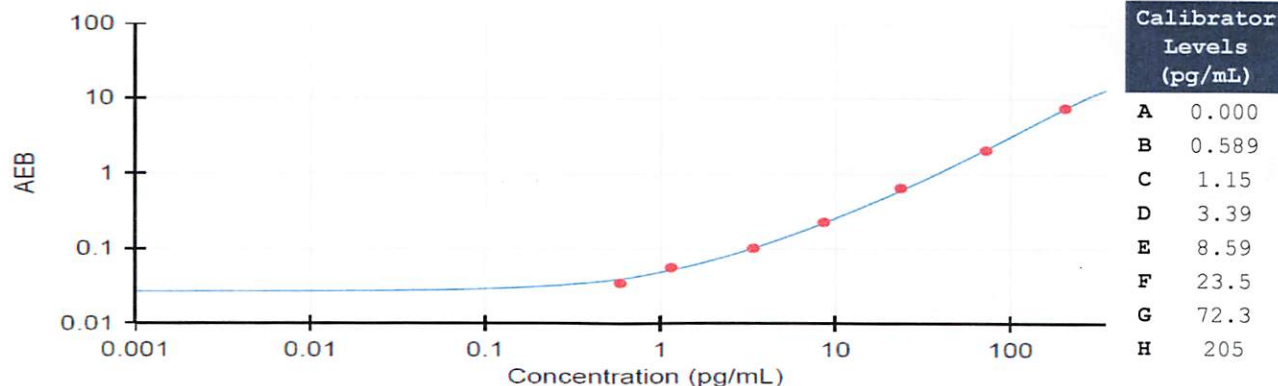
† Specific for use on the SR-X platform - refer to the Simoa™ pTau-231 SR-X Kit Instructions (KI-0110).



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	218106	14.5	7.83-17.1	
Control 2	218106	311	201-352	
Panel 1	111816	31.3	23.7-42.4	
Panel 2	111816	31.1	22.0-40.2	
Panel 3	111816	5.81	4.41-16.9	
Panel 4	111816	5.56	3.77-10.6	

Calibration Curve



Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 04 AUG 2022
Signature/Date