

Product Number:	102292
Lot Number:	502387
Expiration:	21-Feb-2021
Platform(s):	HD-1/HD-X and SR-X



Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

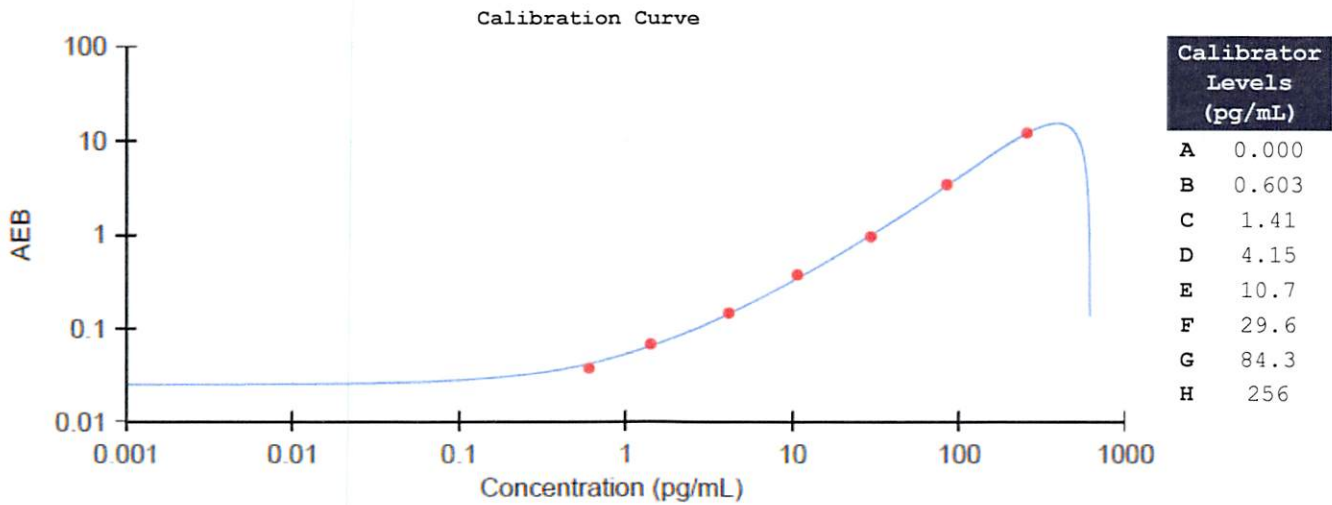
Component	Part Number	Lot Number
Bead Reagent	102280	013803
Detector Reagent	102282	013805
SBG Reagent	102284	013806
Sample Diluent	102286	013809
Calibrators	102289	014301
Controls	102290-102291	014302
SBG Diluent (SRX) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	014302	15.2	10.2-15.2
Control 2	014302	339	267-401
Panel 1	823901	24.1	15.7-31.3
Panel 2	823902	18.9	107-24.2
Panel 3	823903	57.1	44.7-71.6
Panel 4	823904	89.2	72.7-122

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



Review/Approval	Shivani Goel	QC Manager	Shivani Goel	06 09 Jun 2020
	Name	Title	Signature/Date	