

Product Number:	102292
Lot Number:	503032
Expiration:	10-Jun-2022
Platform(s):	HD-1 & SR-X



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

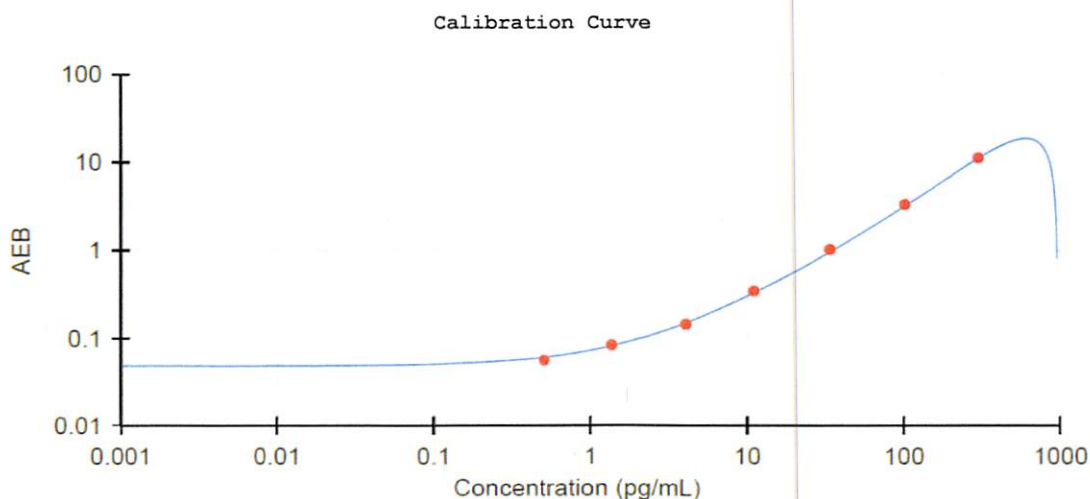
Component	Part Number	Lot Number
Bead Reagent	102280	121040
Detector Reagent	102282	118913
SBG Reagent	102284	118914
Sample Diluent	102286	118917
Reference Calibrator	102289	118911
Control 1	102290	118912
Control 2	102291	118912
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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).

Release Materials			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	118912	14.8	8.56-25.3
Control 2	118912	393	338-507
Panel 1	111816	33.1	23.7-42.4
Panel 2	111816	28.5	22.0-40.2
Panel 3	111816	15.0	4.41-16.9
Panel 4	111816	6.65	3.77-10.6

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



Review/Approval

Junpil Park
Name

QC Supervisor
Title


Signature/Date

12 Aug 2021