

Product Number:	103400
Lot Number:	503547
Expiration:	3-Aug-2023
Platform(s):	SR-X

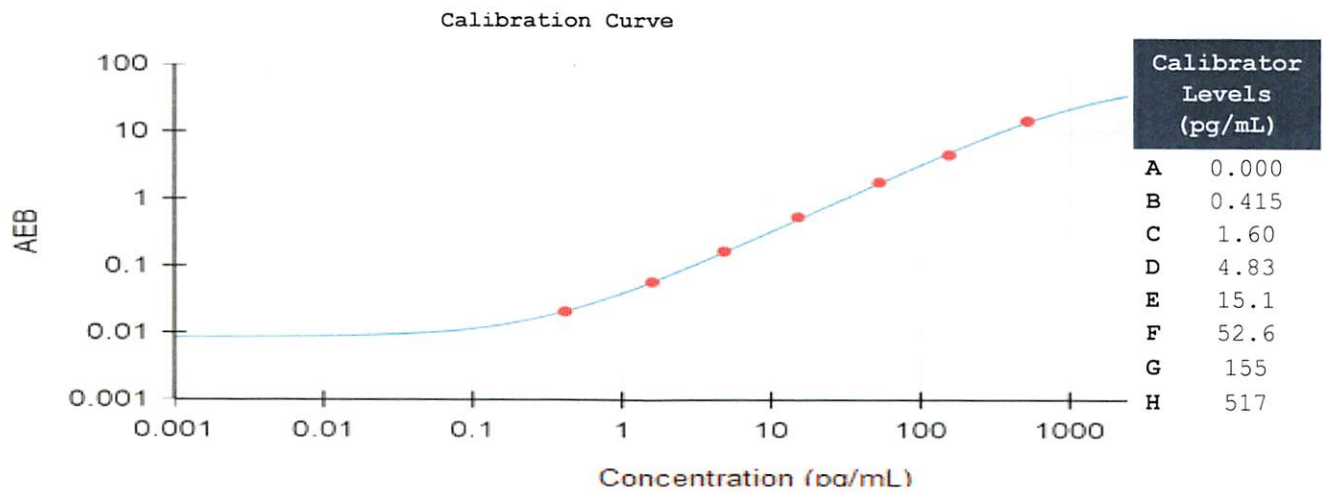


Data below represents results generated on the Quanterix SR-X® Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	222101
Detector Reagent	102248	222107
SBG Reagent	102250	222108
Sample Diluent	103399	224304
Calibrators	102255	227201
Control 1	102256	227202
Control 2	102257	227202
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP Reagents are not Kit Lot Specific

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	227202	10.1	7.61-11.5	
Control 2	227202	405	317-476	
Panel 1	224402	9.83	7.63-12.7	
Panel 2	224402	8.24	5.89-9.40	
Panel 3	224402	70.5	60.5-92.6	
Panel 4	212511	270	150-356	



Review/Approval Matt Kouloouris QC Supervisor [Signature] 20 Sept 2023

Name Title Signature/Date