

Product Number:	103400
Lot Number:	503333
Expiration:	7-Feb-2023
Platform(s):	SR-X

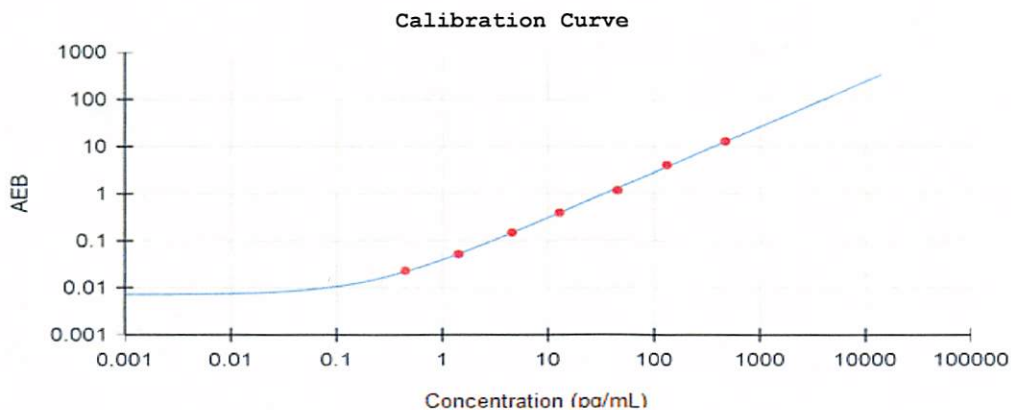


Data below represents results generated on the Quanterix SR-X® Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	203806
Detector Reagent	102248	203910
SBG Reagent	102250	203911
Sample Diluent	103399	204301
Calibrators	102255	203104
Control 1	102256	203105
Control 2	102257	203105
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)	Mean Range (pg/mL)	
Control 1	203105	3.73	2.91-4.38	
Control 2	203105	257	160-285	
Panel 1	134208	6.28	3.95-9.43	
Panel 2	134208	3.36	1.86-7.15	
Panel 3	134208	60.6	49.9-74.8	
Panel 4	135801	381	274-411	



Calibrator Levels (pg/mL)	
A	0.000
B	0.448
C	1.42
D	4.53
E	12.8
F	45.1
G	133
H	472

Review/Approval Shivani Goel QC Manager Shivani Goel 31 Mar 2022  
 Name Title Signature/Date