

Product Number:	103400
Lot Number:	502424
Expiration:	17-Apr-2021
Platform(s):	SR-X

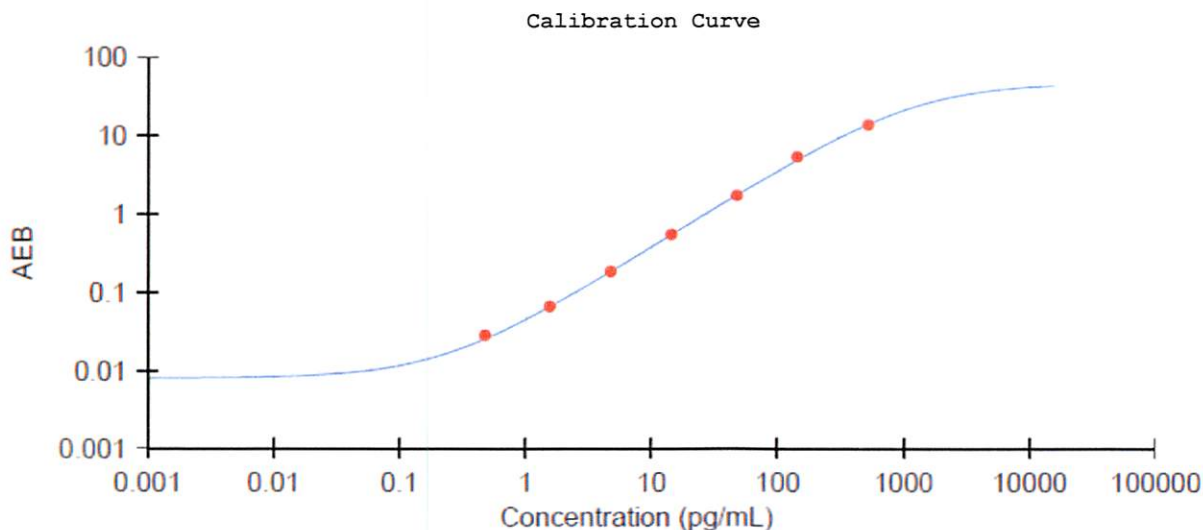


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	010802
Detector Reagent	102248	010803
SBG Reagent	102250	010902
Sample Diluent	103399	011101
Calibrators	102255	011201
Controls	102256/102257	011301
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	011301	5.92	3.64-6.17	
Control 2	011301	212	149-224	
Panel 1	922601	21.1	14.1-21.1	
Panel 2	922602	6.79	4.94-7.41	
Panel 3	922603	42.0	28.4-42.6	
Panel 4	922604	143	129-204	



**Calibrator Levels (pg/mL)**

A	0.000
B	0.475
C	1.56
D	4.71
E	14.4
F	47.4
G	143
H	521

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 17Jun2020  
Signature/Date