

Product Number:	103400
Lot Number:	502254
Expiration:	14-Oct-2020
Platform(s):	SR-X

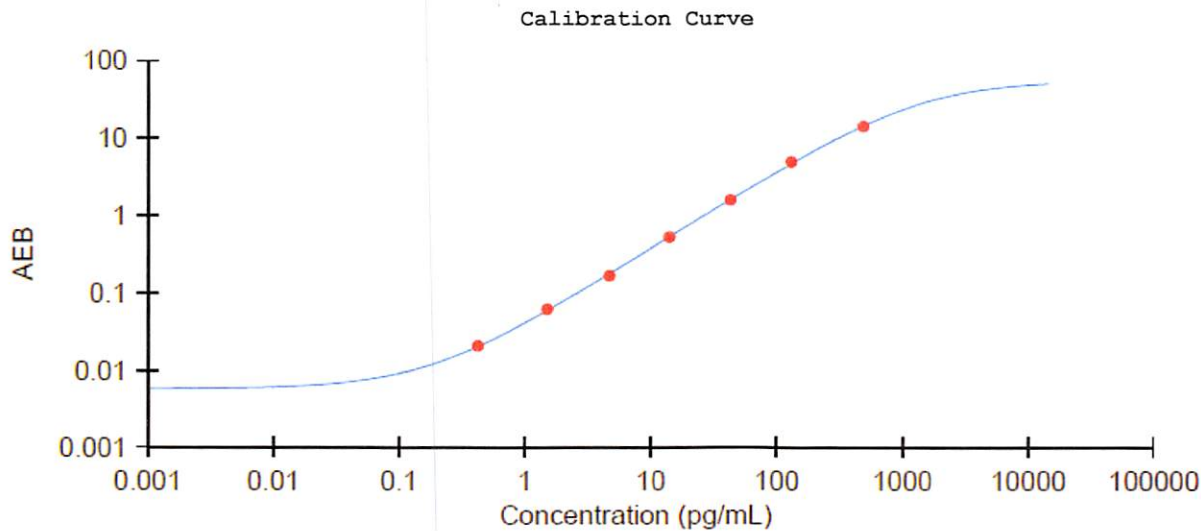


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	002705
Detector Reagent	102248	928805
SBG Reagent	102250	928704
Sample Diluent	103399	004501
Calibrators	102255	003501
Controls	102256/102257	003601
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	003601	4.68	3.33-4.99	
Control 2	003601	178	114-234	
Panel 1	922601	18.2	14.1-21.1	
Panel 2	922602	5.35	4.94-7.41	
Panel 3	922603	36.5	28.4-42.6	
Panel 4	922604	160	129-204	



Calibrator Levels (pg/mL)	
A	0.000
B	0.419
C	1.50
D	4.66
E	14.0
F	42.5
G	130
H	484

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>
	Name	Title	Signature/Date