

Product Number:	103400
Lot Number:	502224
Expiration:	7-Nov-2020
Platform(s):	SR-X

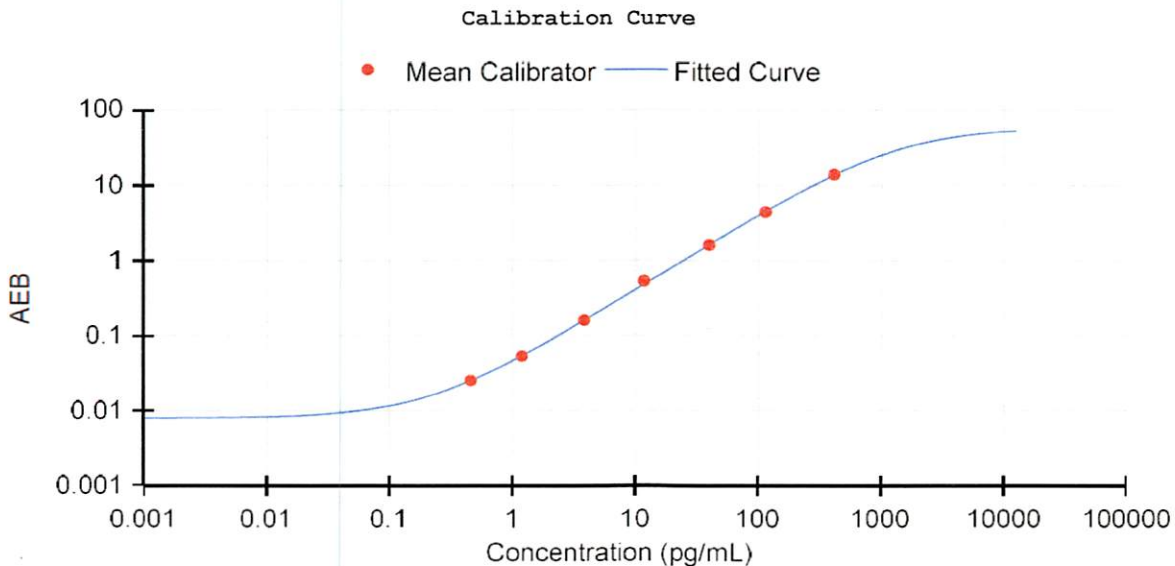


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	002705
Detector Reagent	102248	002701
SBG Reagent	102250	002711
Sample Diluent	103399	931110
Calibrators	102255	935001
Controls	102256/102257	935701
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	935701	3.72	2.99-5.24	
Control 2	935701	151	130-195	
Panel 1	922601	16.6	14.1-21.1	
Panel 2	922602	5.35	4.94-7.41	
Panel 3	922603	38.6	28.4-42.6	
Panel 4	922604	140	129-204	



Calibrator Levels (pg/mL)	
A	0.000
B	0.461
C	1.20
D	3.87
E	11.8
F	40.2
G	115
H	415

<b>Review/Approval</b>	Patrizia Stadler Name	Assoc. Director, QC <i>p Stadler</i> Title	<i>06Feb 2020</i> Signature/Date
------------------------	--------------------------	---	-------------------------------------