

<b>Product Number:</b>	103400
<b>Lot Number:</b>	501702
<b>Expiration:</b>	15-Jan-2020
<b>Platform(s):</b>	HD-1 and SR-X



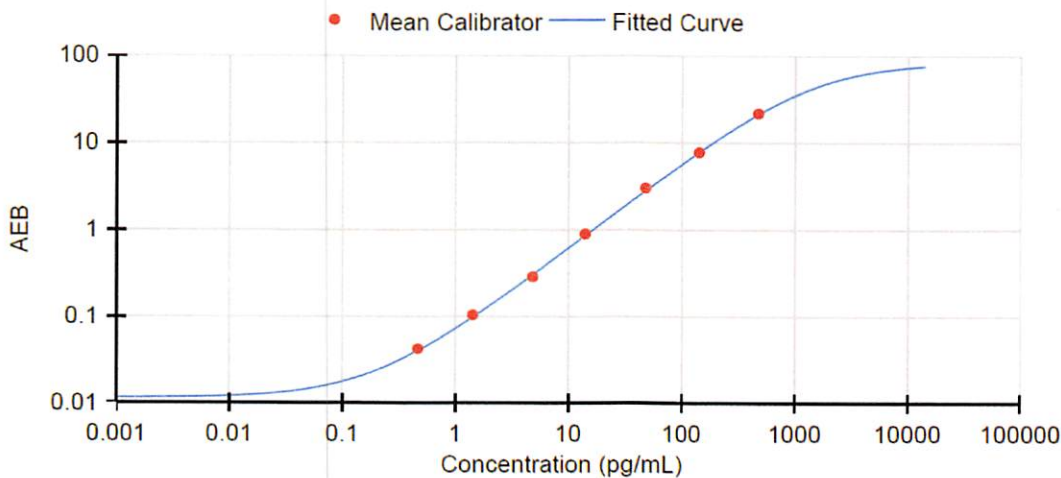
Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	901504
Detector Reagent	102248	901604
SBG Reagent	102250	901603
Sample Diluent	103399	905601
Calibrators	102255	906401
Controls	102256-102257	906402
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	906402	3.70	3.06-4.94	
Control 2	906402	162	129-193	
Panel 1	823301	4.51	4.08-6.68	
Panel 2	823302	6.75	4.90-7.61	
Panel 3	823601	89.5	59.3-101	
Panel 4	823601	352	236-402	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.466
C	1.42
D	4.79
E	14.0
F	47.6
G	142
H	472

Review/Approval	Assoc.	
	Patrizia Stadler	Director, QC
	<i>P. Stadler</i>	11 Mar 2019
	<b>Name</b>	<b>Title</b>
		<b>Signature/Date</b>