

Product Number:	103400
Lot Number:	501940
Expiration:	16-Feb-2020
Platform(s):	HD-1/HD-X and SR-X

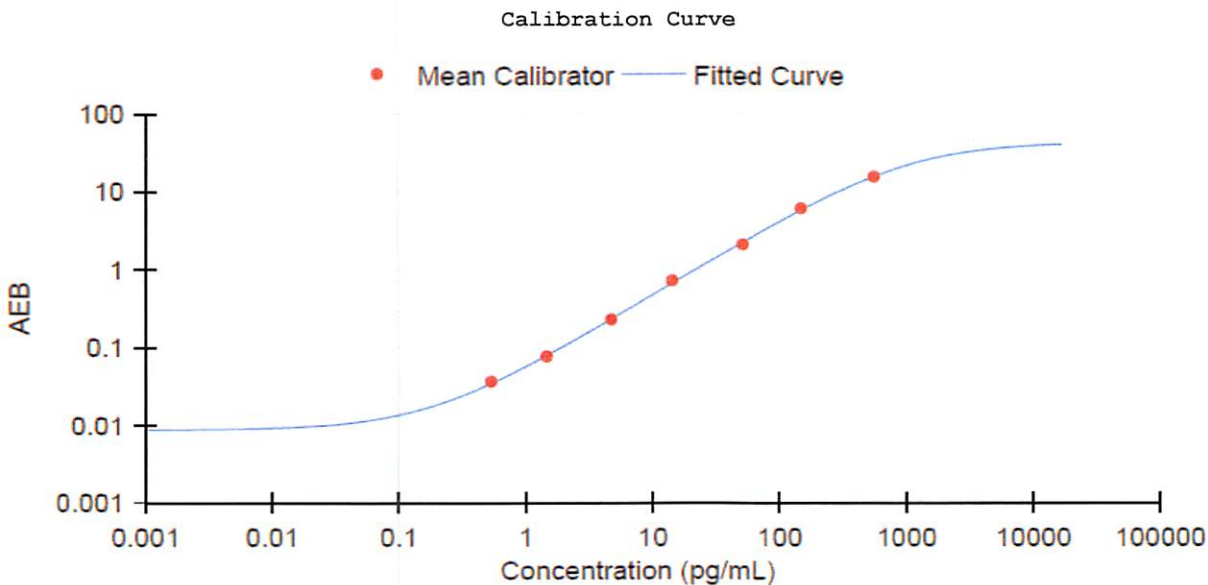


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	906503
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	103399	910505
Calibrators	102255	910701
Controls	102256/102257	910702
RGP Reagent	103159	N/A ¹

¹ 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	910702	5.21	4.38-6.78	
Control 2	910702	149	127-191	
Panel 1	823301	5.29	4.08-6.68	
Panel 2	823302	6.42	4.90-7.61	
Panel 3	823601	85.7	59.3-101	
Panel 4	823602	306	236-402	



Calibrator Levels (pg/mL)	
A	0.000
B	0.535
C	1.46
D	4.70
E	14.2
F	50.9
G	147
H	549

Review/Approval	Patrizia Stadler Name	Assoc. Director, QC Title	<i>p. stadler</i> 18 Sep 2019 Signature/Date
------------------------	--------------------------	------------------------------	---