

Product Number:	103400
Lot Number:	502114
Expiration:	14-Oct-2020
Platform(s):	HD-1/HD-X and SR-X



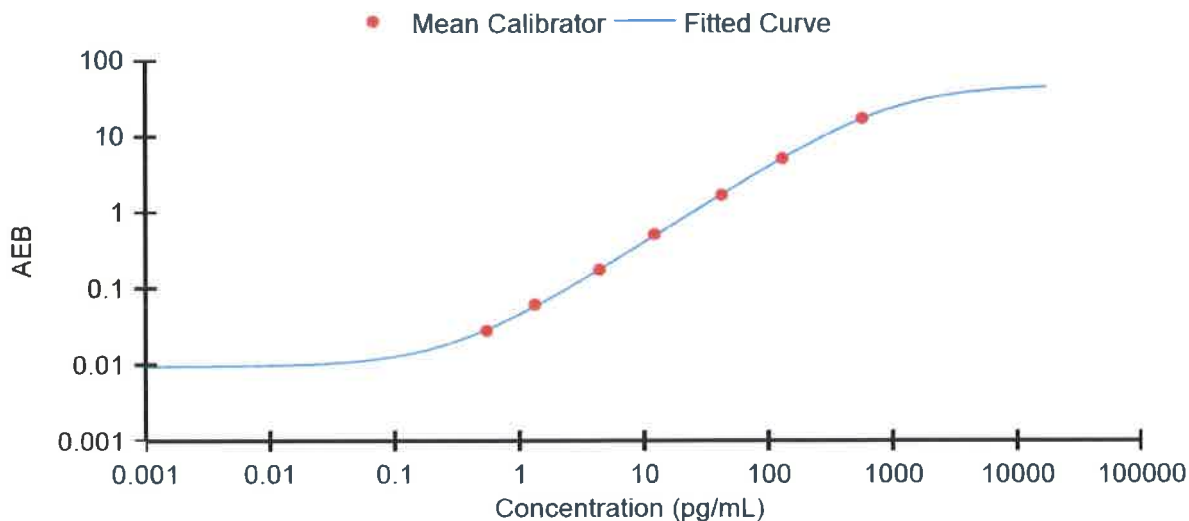
Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	928701
Detector Reagent	102248	928805
SBG Reagent	102250	928704
Sample Diluent	103399	931110
Calibrators	102255	925601
Controls	102256/102257	925901
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	925901	3.62	2.58-4.81	
Control 2	925901	142	99-173	
Panel 1	922601	17.2	14.1-21.1	
Panel 2	922602	5.74	4.94-7.41	
Panel 3	922603	38.8	28.4-42.6	
Panel 4	922604	153	129-204	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.542
C	1.31
D	4.34
E	12.0
F	41.8
G	129
H	567

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