

Product Number:	103400
Lot Number:	502165
Expiration:	7-Nov-2020
Platform(s):	HD-1/HD-X and SR-X

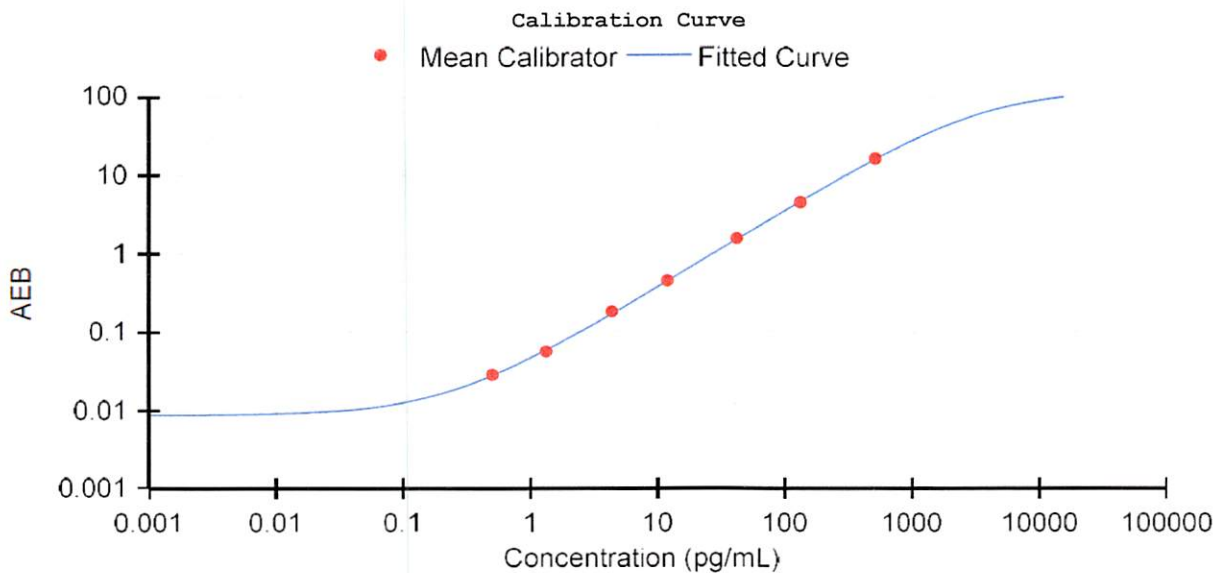


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	933607
Detector Reagent	102248	932706
SBG Reagent	102250	931201
Sample Diluent	103399	931110
Calibrators	102255	933601
Controls	102256/102257	933901
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	933901	3.43	2.43-4.80	
Control 2	933901	156	118-183	
Panel 1	922601	17.7	14.1-21.1	
Panel 2	922602	5.80	4.94-7.41	
Panel 3	922603	41.2	28.4-42.6	
Panel 4	922604	173	129-204	



Calibrator Levels (pg/mL)	
A	0.000
B	0.501
C	1.32
D	4.32
E	11.8
F	41.1
G	131
H	506

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 18 Dec 2019
	Name	Title	Signature/Date