

Product Number:	103400
Lot Number:	502262
Expiration:	29-Jan-2021
Platform(s):	SR-X

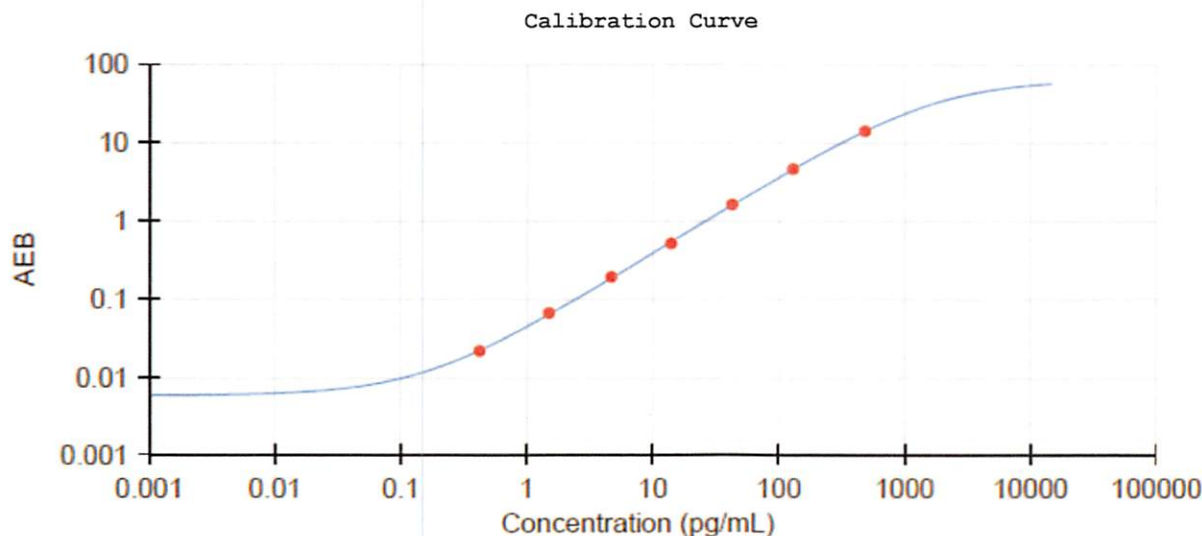


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	005504
Detector Reagent	102248	005005
SBG Reagent	102250	005001
Sample Diluent	103399	004501
Calibrators	102255	003501
Controls	102256/102257	003601
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	003601	4.53	3.33-4.99	
Control 2	003601	181	114-234	
Panel 1	922601	18.3	14.1-21.1	
Panel 2	922602	5.76	4.94-7.41	
Panel 3	922603	34.6	28.4-42.6	
Panel 4	922604	153	129-204	



Calibrator Levels (pg/mL)	
A	0.000
B	0.419
C	1.50
D	4.66
E	14.0
F	42.5
G	130
H	484

Review/Approval: Patrizia Stadler Assoc. Director, QC P. Stadler 04 Feb 2020
 Name Title Signature/Date