

Product Number:	103400
Lot Number:	502370
Expiration:	24-Mar-2021
Platform(s):	SR-X

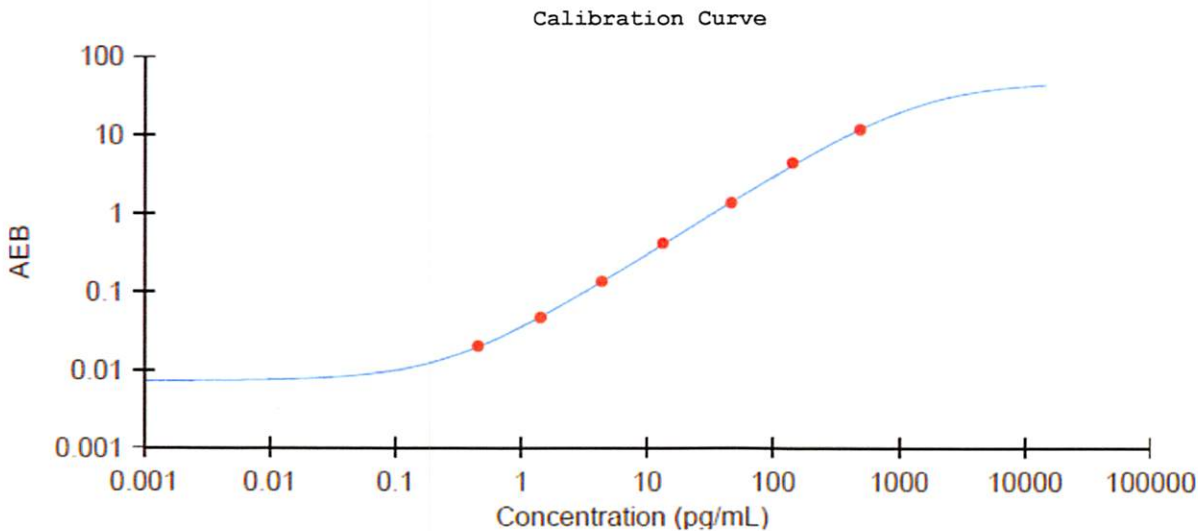


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	010802
Detector Reagent	102248	010803
SBG Reagent	102250	010902
Sample Diluent	103399	011101
Calibrators	102255	007001
Controls	102256/102257	007002
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	007002	4.20	2.52-5.10	
Control 2	007002	157	124-186	
Panel 1	922601	18.9	14.1-21.1	
Panel 2	922602	6.13	4.94-7.41	
Panel 3	922603	38.6	28.4-42.6	
Panel 4	922604	142	129-204	



Calibrator Levels (pg/mL)
A 0.000
B 0.474
C 1.36
D 4.58
E 14.1
F 51.4
G 152
H 526

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 13 May 2020
	Name	Title	Signature/Date