

Product Number:	103400
Lot Number:	502308
Expiration:	26-Feb-2021
Platform(s):	SR-X

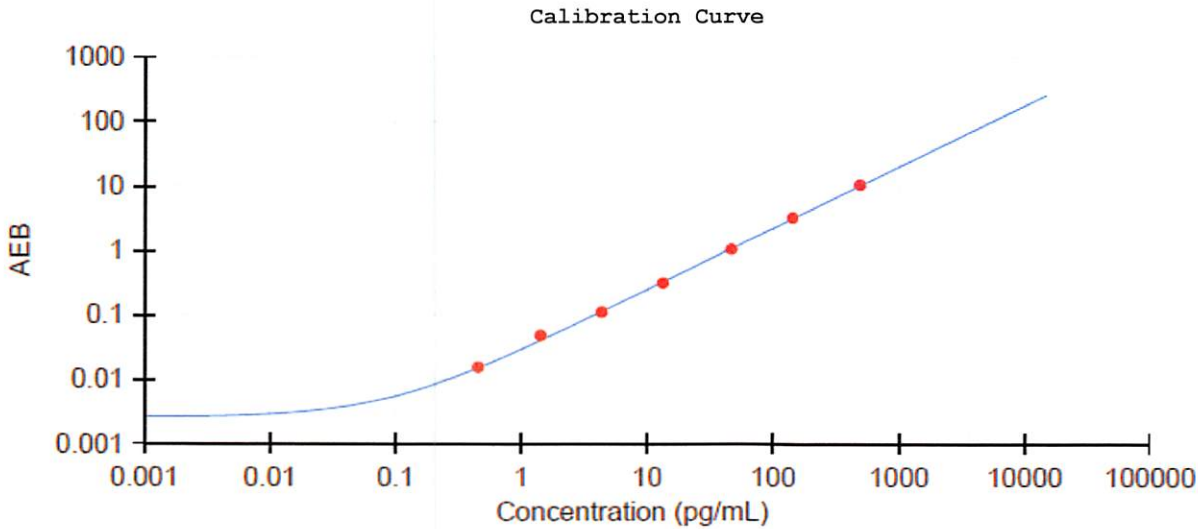


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	007811
Detector Reagent	102248	007905
SBG Reagent	102250	007907
Sample Diluent	103399	006311
Calibrators	102255	008001
Controls	102256/102257	008002
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	008002	3.72	2.43-4.30	
Control 2	008002	156	123-184	
Panel 1	922601	18.6	14.1-21.1	
Panel 2	922602	6.04	4.94-7.41	
Panel 3	922603	37.7	28.4-42.6	
Panel 4	922604	166	129-204	



Review/Approval	Patrizia Stadler Name	Assoc. Director, QC Title	<i>P. Stadler</i> 31 Nov 2020 Signature/Date
-----------------	--------------------------	------------------------------	---