

Product Number:	103400
Lot Number:	502190
Expiration:	14-Oct-2020
Platform(s):	SR-X

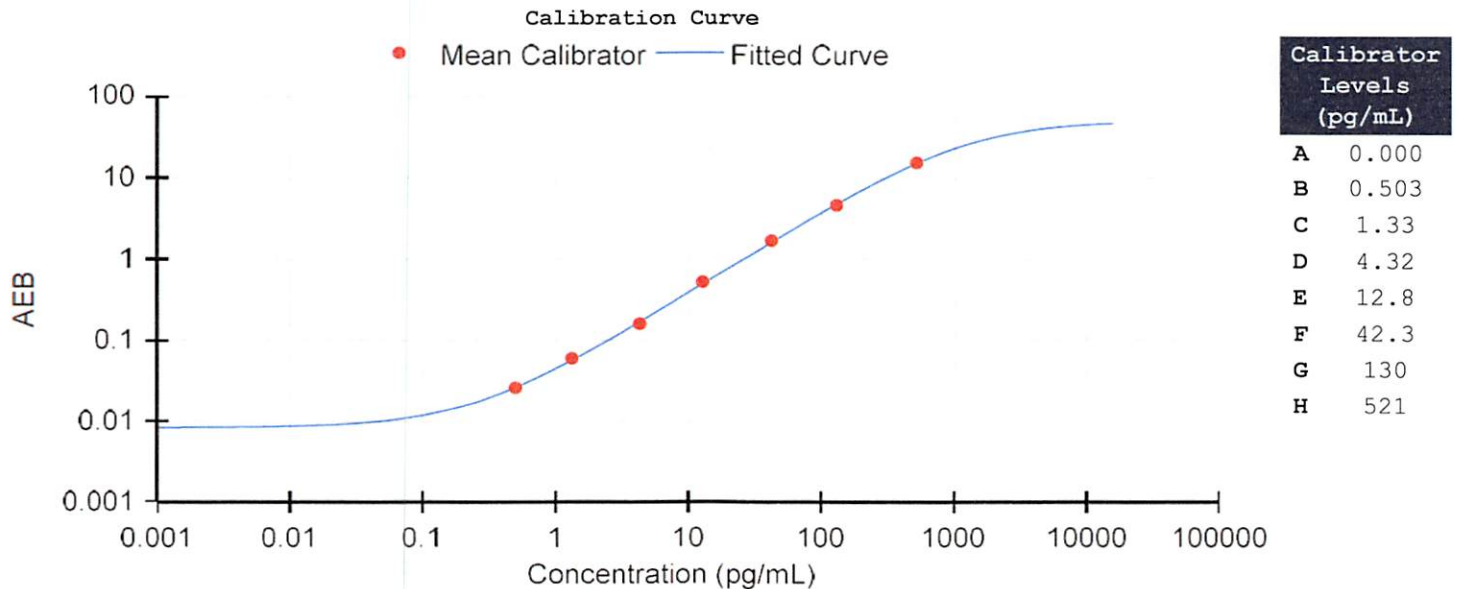


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	928701
Detector Reagent	102248	928805
SBG Reagent	102250	928704
Sample Diluent	103399	931110
Calibrators	102255	924701
Controls	102256/102257	924803
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	924803	3.54	2.50-4.49	
Control 2	924803	136	116-173	
Panel 1	922601	17.4	14.1-21.1	
Panel 2	922602	5.74	4.94-7.41	
Panel 3	922603	38.6	28.4-42.6	
Panel 4	922604	139	129-204	



<b>Review/Approval</b>	Patrizia Stadler	Assoc. Director, QC	<i>p. Stadler</i> 31 Jan 2020
	Name	Title	Signature/Date