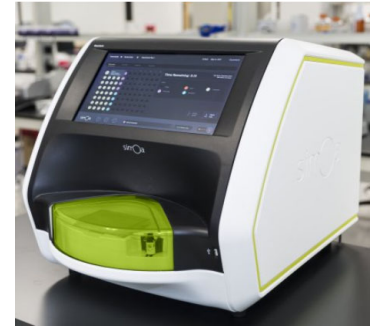


<b>Product Number:</b>	103400
<b>Lot Number:</b>	502114
<b>Expiration:</b>	14-Oct-2020
<b>Platform(s):</b>	HD-1/HD-X and SR-X



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

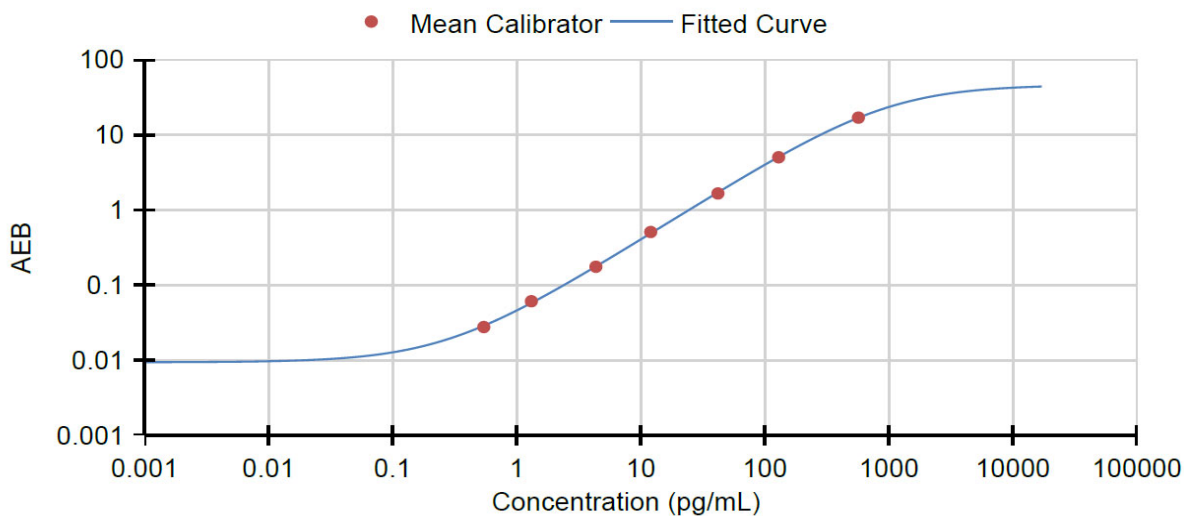
Component	Part Number	Lot Number
Bead Reagent	102246	928701
Detector Reagent	102248	928805
SBG Reagent	102250	928704
Sample Diluent	103399	931110
Calibrators	102255	925601
Controls	102256/102257	925901
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific.

Release Materials			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	925901	3.62	2.58-4.81
Control 2	925901	142	99-173
Panel 1	922601	17.2	14.1-21.1
Panel 2	922602	5.74	4.94-7.41
Panel 3	922603	38.8	28.4-42.6
Panel 4	922604	153	129-204

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.542
C	1.31
D	4.34
E	12.0
F	41.8
G	129
H	567

<b>Review/Approval</b>	Patrizia Stadler	Assoc. Director, QC	
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>