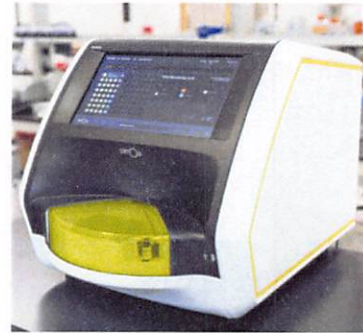


Product Number:	103400
Lot Number:	503470
Expiration:	3-May-2023
Platform(s):	SR-X

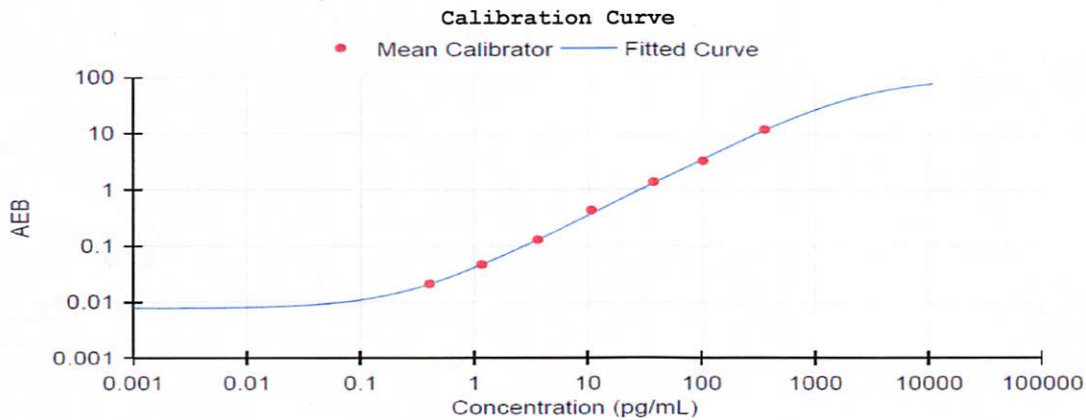


Data below represents results generated on the Quanterix SR-X® Ultra-Sensitive Biomarker Detection System

Component	Part Number	Lot Number
Bead Reagent	102246	214407
Detector Reagent	102248	214201
SBG Reagent	102250	214202
Sample Diluent	103399	215701
Calibrators	102255	213603
Control 1	102256	213604
Control 2	102257	213604
RGP Reagent	103159	N/A ¹

¹ RGP 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	213604	3.47	2.17-4.36	
Control 2	213604	109	103-155	
Panel 1	134208	7.82	3.95-9.43	
Panel 2	134208	5.42	1.86-7.15	
Panel 3	212511	90.7	73.5-116	
Panel 4	212511	248	145-372	



Calibrator Levels (pg/mL)	
A	0.000
B	0.407
C	1.17
D	3.62
E	10.8
F	37.7
G	103
H	359

Review/Approval Matt Smart Director, QC MRA 28 Jun 2022
Name
Title
Signature/Date