

Product Number: 103400
 Lot Number: 502609
 Expiration: 14-Sep-2021
 Platform(s): SR-X



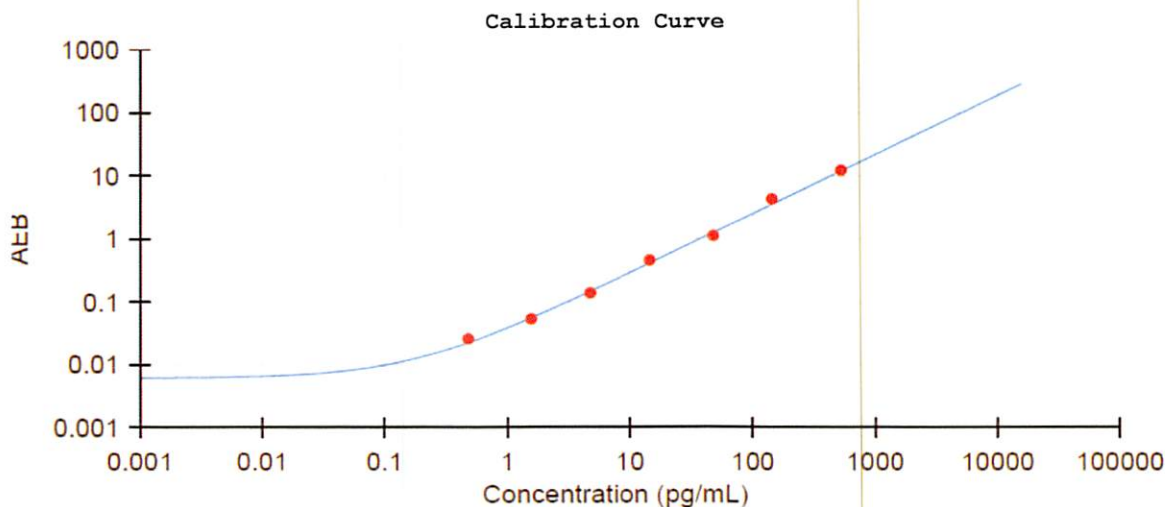
Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	028804
Detector Reagent	102248	029003
SBG Reagent	102250	031118
Sample Diluent	103399	029311
Calibrators	102255	011201
Control 1	102256	011301
Control 2	102257	011301
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	011301	6.30	3.56-7.16
Control 2	011301	264	113-287
Panel 1	922601	21.3	14.3-21.4
Panel 2	922602	7.18	4.97-7.45
Panel 3	922603	41.1	28.8-43.2
Panel 4	922604	197	133-201

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



Calibrator Levels (pg/mL)	
A	0.000
B	0.475
C	1.56
D	4.71
E	14.4
F	47.4
G	143
H	521

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 14/13 Nov 2020
 SG 13 Nov 2020