

Product Number:	103400
Lot Number:	502626
Expiration:	24-Sep-2021
Platform(s):	SR-X

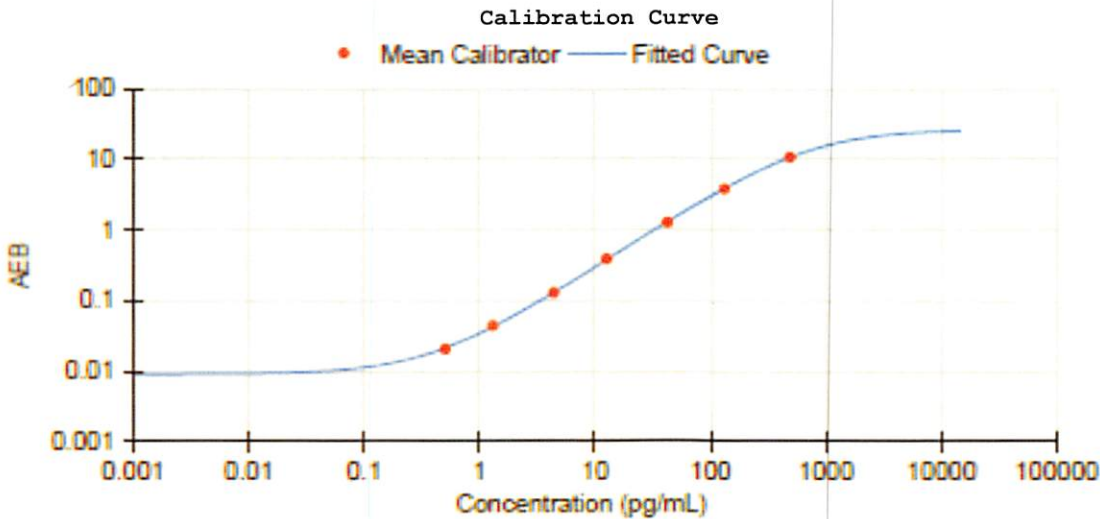


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	026804
Detector Reagent	102248	026904
SBG Reagent	102250	026908
Sample Diluent	103399	029311
Calibrators	102255	027308
Control 1	102256	027309
Control 2	102257	027309
RGP Reagent	103159	N/A ¹

¹ 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	027309	4.23	2.83-4.78	
Control 2	027309	151	120-181	
Panel 1	008501	15.5	9.91-18.2	
Panel 2	008501	10.7	8.15-13.3	
Panel 3	008501	67.9	47.2-84.0	
Panel 4	009401	202	144-253	



Calibrator Levels (pg/mL)	
A	0.000
B	0.505
C	1.31
D	4.39
E	12.7
F	42.5
G	132
H	487

Review/Approval

Junpil Park

QC Supervisor

[Signature] 04 Dec 2020