

Product Number:	103400
Lot Number:	502856
Expiration:	1-Mar-2022
Platform(s):	SR-X

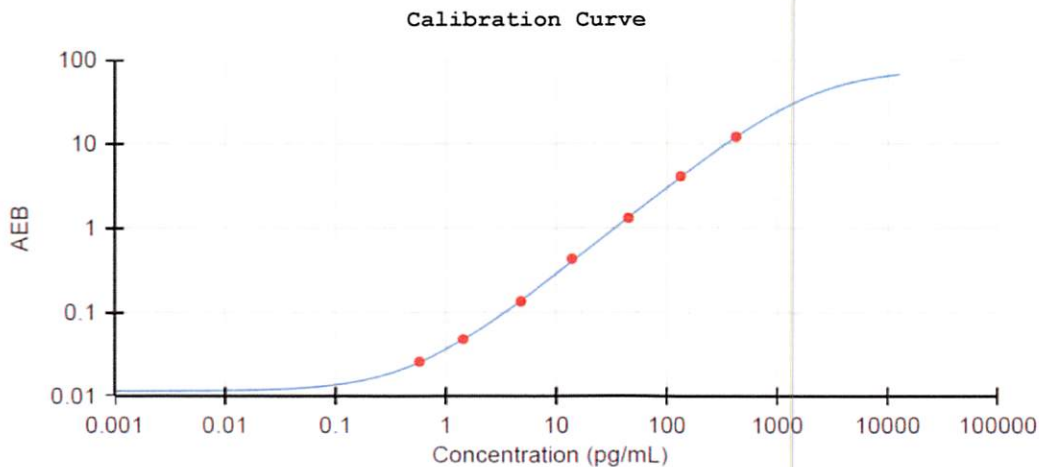


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	106912
Detector Reagent	102248	106911
SBG Reagent	102250	106716
Sample Diluent	103399	106015
Calibrators	102255	106801
Control 1	102256	106818
Control 2	102257	109005
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	106818	5.01	2.73-5.70	
Control 2	109005	295	165-305	
Panel 1	008501	19.3	8.59-20.5	
Panel 2	008501	13.9	7.39-15.1	
Panel 3	008501	85.4	43.8-91.8	
Panel 4	009401	230	144-253	



Calibrator Levels (pg/mL)

A	0.000
B	0.576
C	1.43
D	4.75
E	13.9
F	44.8
G	134
H	424

Review/Approval

Junpil Park

QC Supervisor

15 Jun 2021