

Product Number:	103400
Lot Number:	502975
Expiration:	25-May-2022
Platform(s):	SR-X

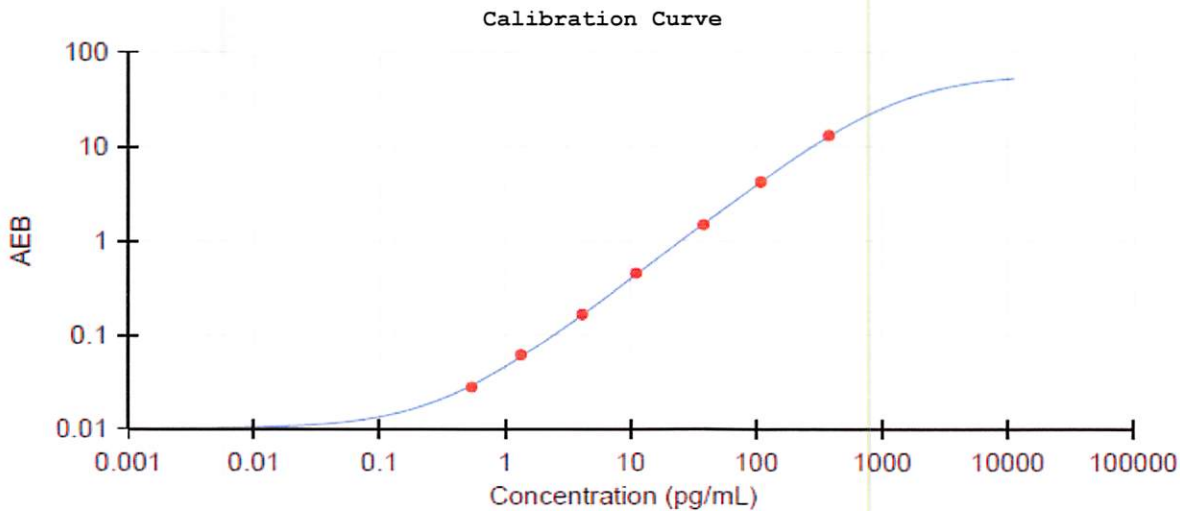


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	114511
Detector Reagent	102248	114512
SBG Reagent	102250	114513
Sample Diluent	103399	115909
Calibrators	102255	115206
Control 1	102256	115207
Control 2	102257	115207
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	115207	3.78	2.91-5.04	
Control 2	115207	236	123-395	
Panel 1	008501	12.0	8.59-20.5	
Panel 2	008501	9.07	7.39-15.1	
Panel 3	008501	76.7	43.8-91.8	
Panel 4	009401	177	128-302	



Calibrator Levels (pg/mL)	
A	0.000
B	0.539
C	1.34
D	4.09
E	11.1
F	37.8
G	108
H	376

Review/Approval: Junpil Park QC Supervisor

 02 Jul 2021