

Product Number:	103400
Lot Number:	502770
Expiration:	15-Dec-2021
Platform(s):	SR-X

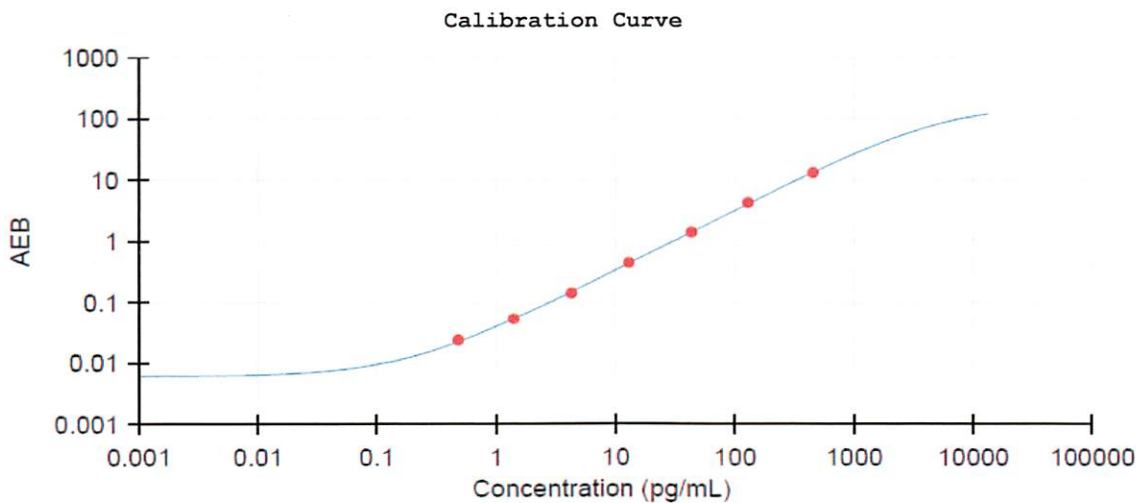


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	104109
Detector Reagent	102248	104101
SBG Reagent	102250	104015
Sample Diluent	103399	035001
Calibrators	102255	104002
Control 1	102256	104003
Control 2	102257	105001
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	104003	3.43	2.52-4.09	
Control 2	105001	162	136-204	
Panel 1	008501	14.9	9.92-18.2	
Panel 2	008501	11.2	8.15-13.3	
Panel 3	008501	66.6	47.2-84.0	
Panel 4	009401	215	144-253	



Review/Approval _____

Junpil Park

QC Supervisor

01 Mar 2021