

<b>Product Number:</b>	103400
<b>Lot Number:</b>	502648
<b>Expiration:</b>	19-Oct-2020
<b>Platform(s):</b>	SR-X

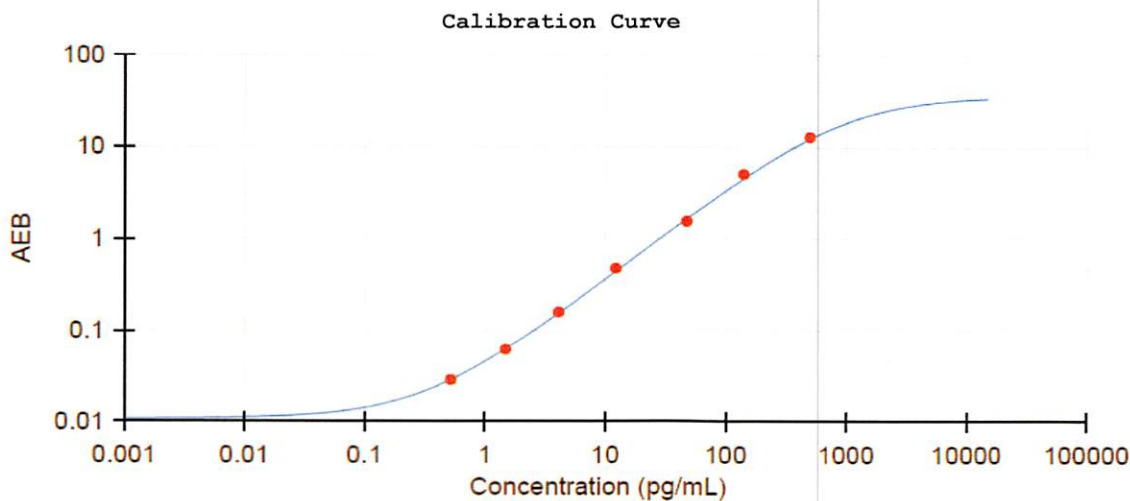


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	032812
Detector Reagent	102248	033502
SBG Reagent	102250	033501
Sample Diluent	103399	029311
Calibrators	102255	033513
Control 1	102256	033602
Control 2	102257	033602
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	033602	3.86	2.97-4.46	
Control 2	033602	170	126-189	
Panel 1	008501	14.9	9.92-18.2	
Panel 2	008501	10.5	8.16-13.3	
Panel 3	008501	69.4	47.2-84.0	
Panel 4	009401	211	144-253	



Review/Approval \_\_\_\_\_ Junpil Park QC Supervisor

*[Signature]* 14 Dec 2020