

Product Number:	103400
Lot Number:	502706
Expiration:	15-Dec-2021
Platform(s):	SR-X

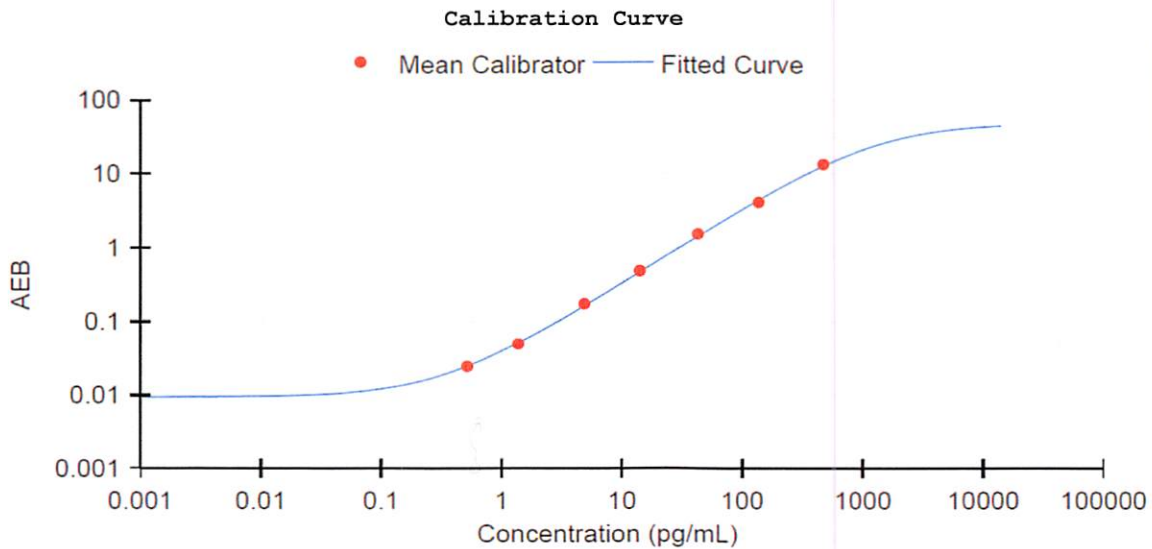


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	101301
Detector Reagent	102248	101516
SBG Reagent	102250	102301
Sample Diluent	103399	035001
Calibrators	102255	102101
Control 1	102256	102203
Control 2	102257	102203
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	102203	4.42	2.70-5.97	
Control 2	102203	190	115-207	
Panel 1	922601	21.1	14.3-21.4	
Panel 2	922602	7.42	4.97-7.45	
Panel 3	922603	41.9	28.8-43.2	
Panel 4	922604	192	133-201	



Review/Approval: Junpil Park QC Supervisor 03 Feb 2021