

Product Number:	103400
Lot Number:	503260
Expiration:	8-Jun-2022
Platform(s):	SR-X

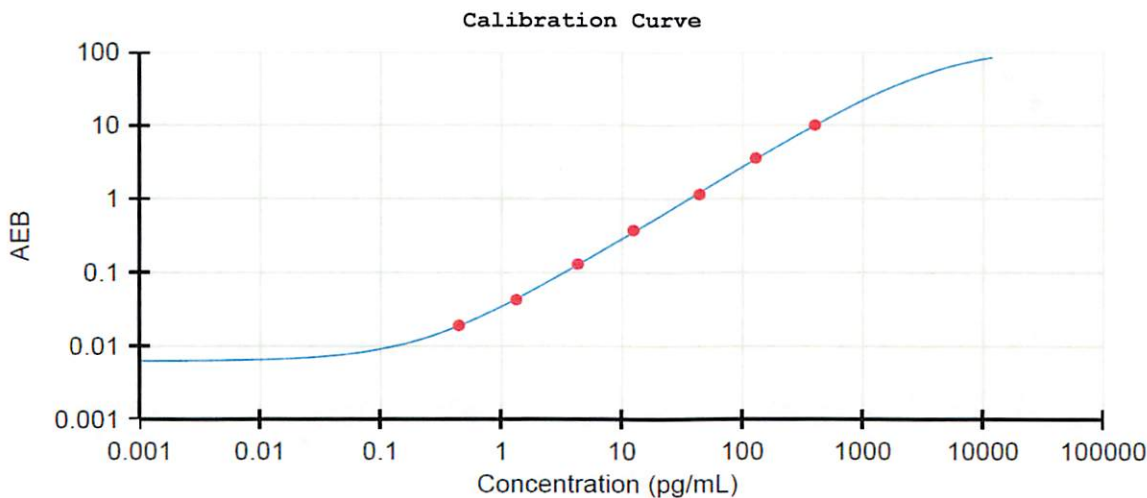


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	124505
Detector Reagent	102248	123907
SBG Reagent	102250	127405
Sample Diluent	103399	115909
Calibrators	102255	133605
Control 1	102256	133606
Control 2	102257	133606
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	133606	3.65	3.02 - 5.23	
Control 2	133606	236	200 - 300	
Panel 1	134208	5.29	3.70 - 9.61	
Panel 2	134208	4.01	2.97 - 6.06	
Panel 3	134208	63.3	49.7 - 74.6	
Panel 4	135801	351	271 - 407	



Review/Approval

Wenka Tsai

QC

[Signature]

14 Jan 2022