

Product Number:	103400
Lot Number:	502500
Expiration:	9-Jul-2021
Platform(s):	SR-X

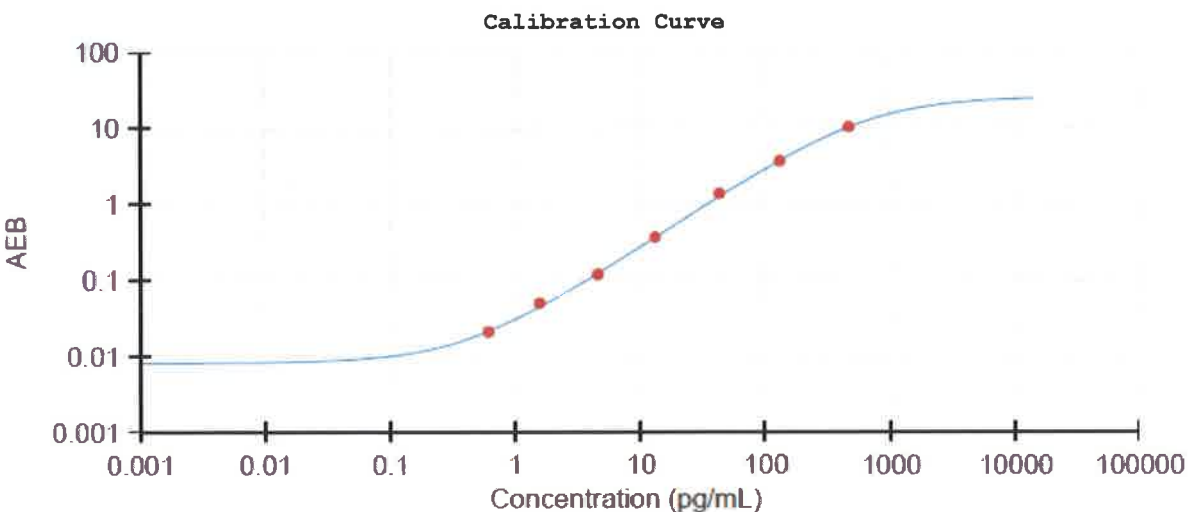


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	022302
Detector Reagent	102248	022404
SBG Reagent	102250	022403
Sample Diluent	103399	019103
Calibrators	102255	022401
Control 1	102256	022402
Control 2	102257	022402
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	022402	5.07	3.41-5.21	
Control 2	022402	150	132-205	
Panel 1	922601	20.8	14.1-21.1	
Panel 2	922602	6.64	4.94-7.41	
Panel 3	922603	41.7	28.4-42.6	
Panel 4	922604	155	129-204	



Calibrator Levels (pg/mL)

A	0.000
B	0.609
C	1.56
D	4.56
E	13.1
F	43.1
G	131
H	468

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 08 Sep 2020