

Product Number:	103400
Lot Number:	502467
Expiration:	18-Jun-2021
Platform(s):	SR-X



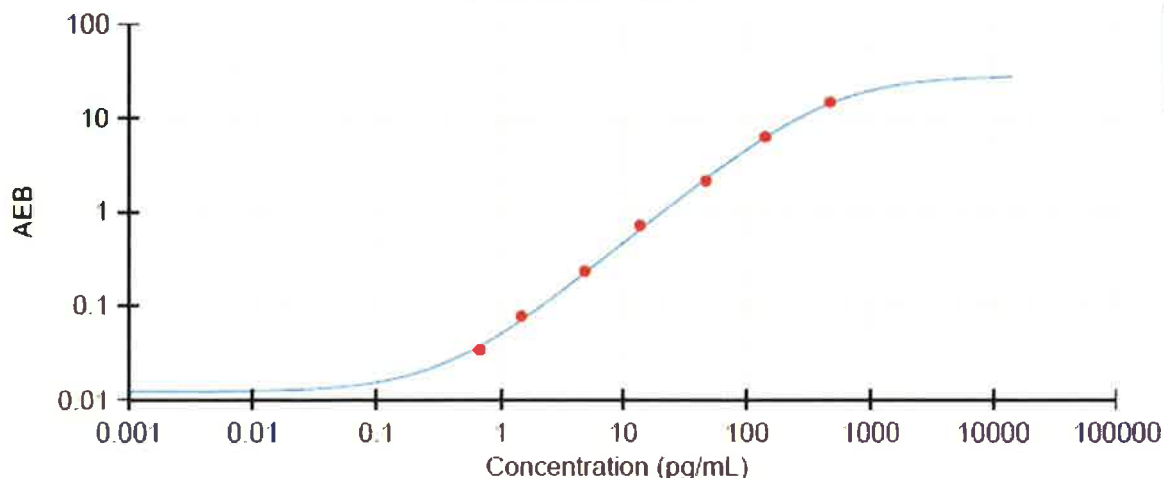
Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	017009
Detector Reagent	102248	017004
SBG Reagent	102250	017007
Sample Diluent	103399	019103
Calibrators	102255	018301
Control 1	102256	018401
Control 2	102257	018401
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	018401	4.65	3.30-4.95	
Control 2	018401	154	117-176	
Panel 1	922601	19.9	14.1-21.1	
Panel 2	922602	7.09	4.94-7.41	
Panel 3	922603	42.6	28.4-42.6	
Panel 4	922604	167	129-204	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.673
C	1.48
D	4.94
E	13.9
F	47.3
G	142
H	472

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 23 Jul 2020