

Product Number:	103400
Lot Number:	502610
Expiration:	14-Oct-2021
Platform(s):	SR-X



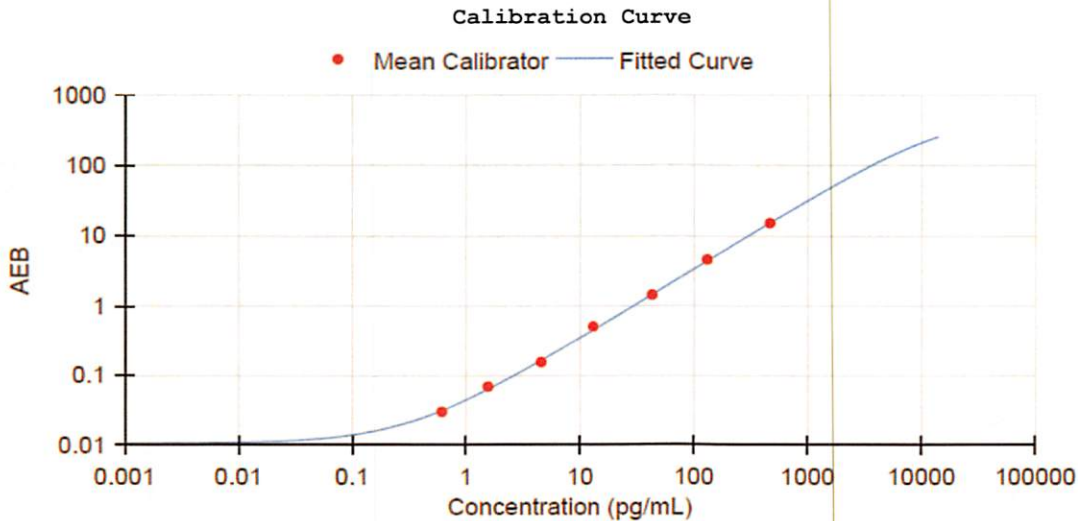
Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	028804
Detector Reagent	102248	029003
SBG Reagent	102250	031118
Sample Diluent	103399	029311
Calibrators	102255	022401
Control 1	102256	022402
Control 2	102257	022402
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

Release Materials			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	022402	4.84	3.36-5.68
Control 2	022402	204	127-226
Panel 1	008501	16.5	9.58-18.0
Panel 2	008501	12.6	7.93-13.5
Panel 3	008501	80.2	45.5-84.8
Panel 4	009401	216	139-256

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.



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: Patrizia Stadler Assoc. Director, QC

*p. Stadler* 18 Nov 2020