

Product Number:	103400
Lot Number:	502555
Expiration:	9-Jul-2021
Platform(s):	SR-X

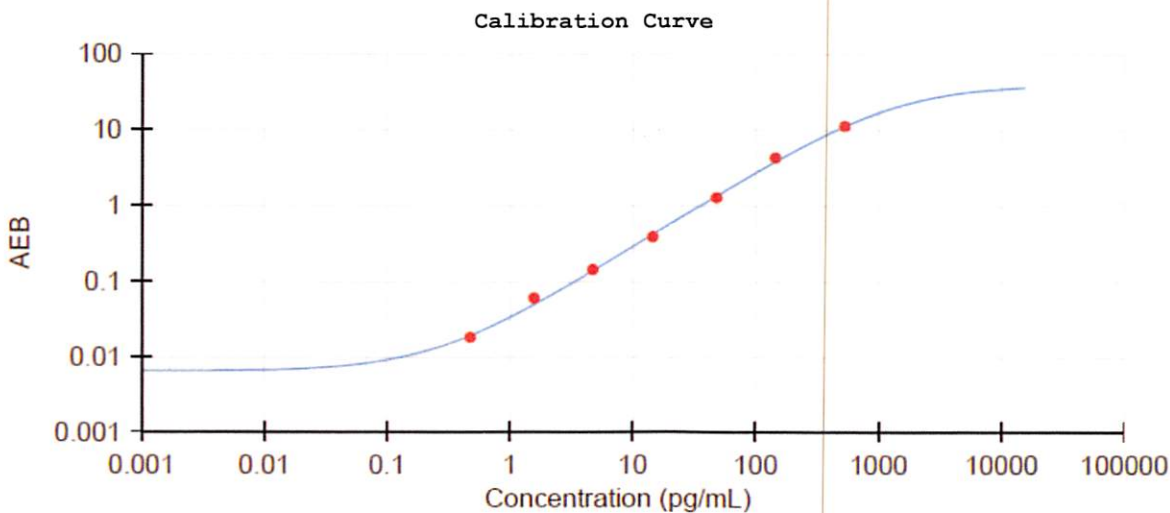


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	026804
Detector Reagent	102248	026904
SBG Reagent	102250	026908
Sample Diluent	103399	019103
Calibrators	102255	011201
Control 1	102256	011301
Control 2	102257	011301
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	011301	5.88	3.64-6.17	
Control 2	011301	205	149-224	
Panel 1	008501	14.6	10.5-15.7	
Panel 2	008501	12.0	8.22-12.3	
Panel 3	008501	68.6	49.1-73.6	
Panel 4	009401	181	151-234	



Calibrator Levels (pg/mL)	
A	0.000
B	0.475
C	1.56
D	4.71
E	14.4
F	47.4
G	143
H	521

Review/Approval: Patrizia Stadler Assoc. Director, QC

P. Stadler

06 OCT 2020