

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

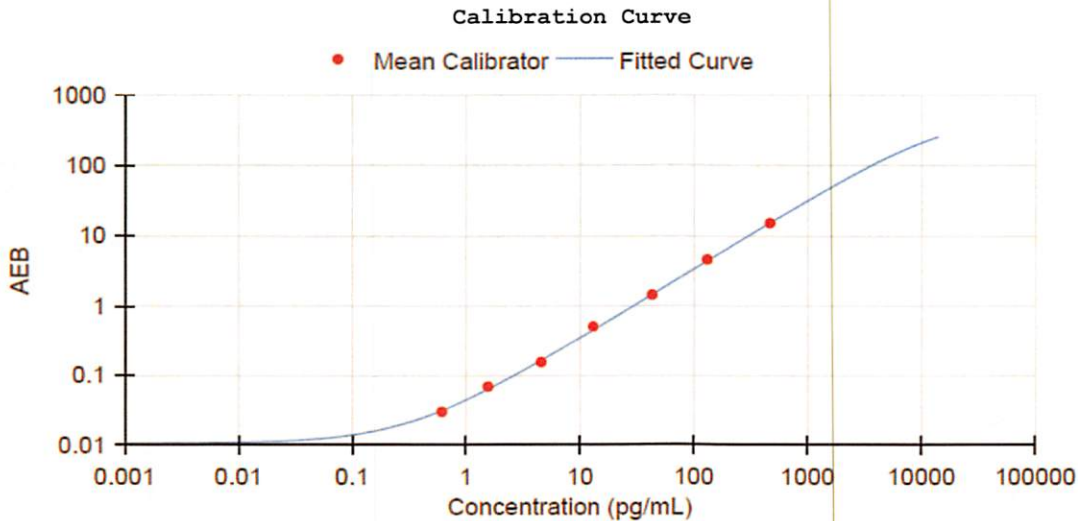


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

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 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	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	



Review/Approval: Patrizia Stadler Assoc. Director, QC

p. Stadler 18 Nov 2020