

Product Number:	103400
Lot Number:	502921
Expiration:	16-Apr-2022
Platform(s):	SR-X



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

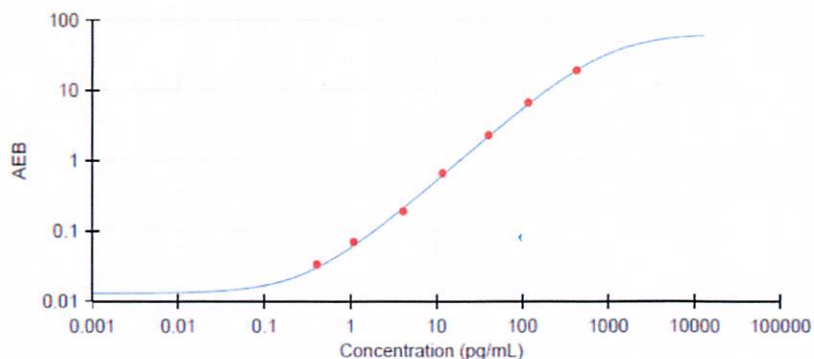
Component	Part Number	Lot Number
Bead Reagent	102246	111804
Detector Reagent	102248	111715
SBG Reagent	102250	111406
Sample Diluent	103399	110609
Calibrators	102255	112301
Control 1	102256	112302
Control 2	102257	112302
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*						* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Previously reported		Corrected **		
		Result (pg/mL)	Mean Range (pg/mL)	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	112302	4.57	2.28-5.08	3.64	1.84-4.08	
Control 2	112302	228	169-257	213	150-247	
Panel 1	008501	13.5	8.59-20.5	11.7	8.59-20.5	
Panel 2	008501	9.55	7.39-15.1	8.20	7.39-15.1	
Panel 3	008501	64.9	43.8-91.8	58.9	43.8-91.8	
Panel 4	009401	195	144-253	181	128-302	

** Values and ranges are updated as of 26Oct2021. Corrected result and calibrator levels should be used for reference.

Calibration Curve



	Calibrator Levels (pg/mL)	
	Previously reported	Corrected**
A	0.000	0.000
B	0.411	0.362
C	1.10	1.00
D	4.12	3.60
E	11.8	11.1
F	40.6	37.1
G	117	114
H	427	377

Review/Approval

Junpil Park

QC Supervisor

[Handwritten Signature] 01 Nov 2021

Name

Title

Signature/Date