

Product Number:	103400
Lot Number:	502866
Expiration:	1-Mar-2022
Platform(s):	SR-X



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

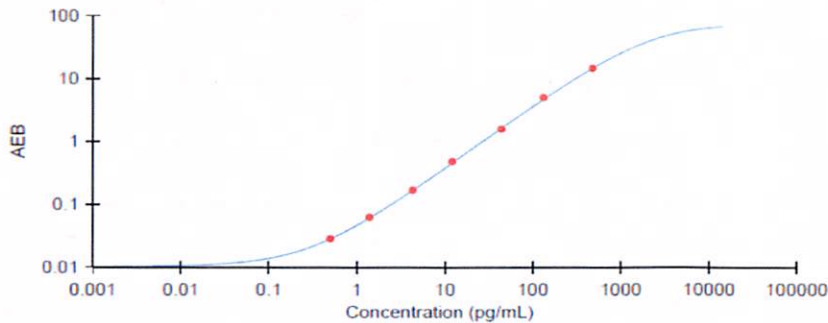
Component	Part Number	Lot Number
Bead Reagent	102246	109016
Detector Reagent	102248	108913
SBG Reagent	102250	108909
Sample Diluent	103399	106015
Calibrators	102255	109806
Control 1	102256	109807
Control 2	102257	109807
RGP Reagent	103159	N/A ¹

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	109807	5.84	3.16-6.41	5.13	3.60-5.40	
Control 2	109807	288	168-319	270	183-274	
Panel 1	008501	12.8	8.59-20.5	12.3	8.59-20.5	
Panel 2	008501	9.25	7.39-15.1	8.87	7.39-15.1	
Panel 3	008501	63.8	43.8-91.8	60.3	43.8-91.8	
Panel 4	009401	200	144-253	189	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	0.506	0.464
C	1.41	1.38
D	4.33	4.35
E	12.2	11.8
F	44.2	41.5
G	132	126
H	477	507

Review/Approval

Junpil Park

QC Supervisor

 01/NOV/2021

Name

Title

Signature/Date