

Product Number:	103400
Lot Number:	502856
Expiration:	1-Mar-2022
Platform(s):	SR-X



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

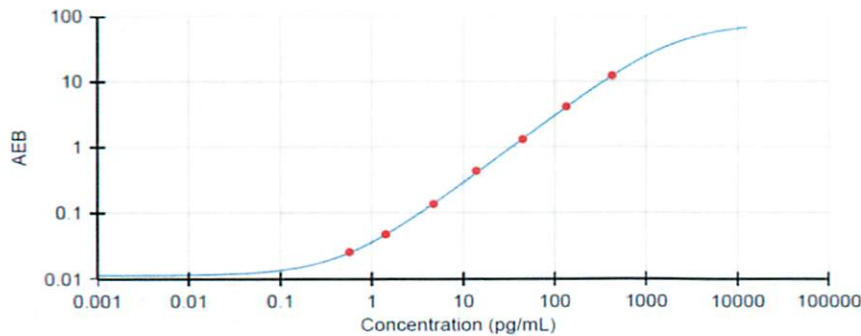
Component	Part Number	Lot Number
Bead Reagent	102246	106912
Detector Reagent	102248	106911
SBG Reagent	102250	106716
Sample Diluent	103399	106015
Calibrators	102255	106801
Control 1	102256	106818
Control 2	102257	109005
RGP Reagent	103159	N/A ¹

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	106818	5.01	2.73-5.70	3.89	2.27-4.32	
Control 2	109005	295	165-305	234	144-250	
Panel 1	008501	19.3	8.59-20.5	15.5	8.59-20.5	
Panel 2	008501	13.86	7.39-15.1	10.9	7.39-15.1	
Panel 3	008501	85.4	43.8-91.8	67.4	43.8-91.8	
Panel 4	009401	230	144-253	182	144-253	

** 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	0.576	0.479
C	1.43	1.12
D	4.75	3.77
E	13.9	10.8
F	44.8	36.4
G	134	108
H	424	350

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 22 NOV 2021
Signature/Date