

Product Number:	103400
Lot Number:	502975
Expiration:	25-May-2022
Platform(s):	SR-X



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

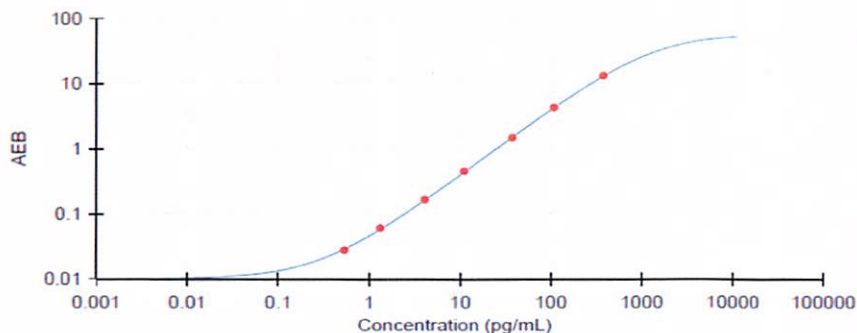
Component	Part Number	Lot Number
Bead Reagent	102246	114511
Detector Reagent	102248	114512
SBG Reagent	102250	114513
Sample Diluent	103399	115909
Calibrators	102255	115206
Control 1	102256	115207
Control 2	102257	115207
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	115207	3.78	2.91-5.04	2.98	2.08-4.77	
Control 2	115207	236	123-395	215	147-383	
Panel 1	008501	12.0	8.59-20.5	9.93	8.59-20.5	
Panel 2	008501	9.07	7.39-15.1	8.38	7.39-15.1	
Panel 3	008501	76.7	43.8-91.8	68.2	43.8-91.8	
Panel 4	009401	177	128-302	161	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	0.539	0.393
C	1.34	1.14
D	4.09	3.44
E	11.1	9.90
F	37.8	33.4
G	108	101
H	376	307

Review/Approval

Junpil Park  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date 01 Nov 2021