

Product Number:	103186
Lot Number:	503019
Expiration:	27-Jun-2022
Platform(s):	HD-1/ HD-X



Data below represents results generated on the Simoa HD-X® Analyzer.

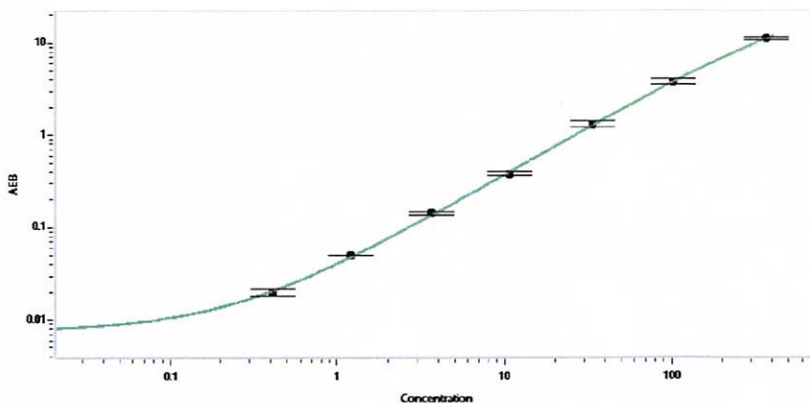
Component	Part Number	Lot Number
Bead Reagent	102246	117801
Detector Reagent	102248	117803
SBG Reagent	102250	117808
Sample Diluent	102252	117914
Calibrators	102255	118706
Control 1	102256	118707
Control 2	102257	118707
RGP Reagent	103159	N/A ¹

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	118707	3.43	2.62-3.93	2.80	2.07-3.74	
Control 2	118707	186	138-214	173	147-220	
Panel 1	008501	14.4	8.59-20.5	12.4	8.59-20.5	
Panel 2	008501	11.4	7.39-15.1	9.76	7.39-15.1	
Panel 3	008501	61.4	43.8-91.8	55.5	43.8-91.8	
Panel 4	009401	189	128-302	176	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	0.410	0.337
C	1.21	1.01
D	3.67	3.08
E	10.6	9.15
F	33.6	32.0
G	102	97.7
H	369	339

Review/Approval	Junpil Park	QC Supervisor		01/Nov/2021
	Name	Title	Signature/Date	