

Product Number:	103186
Lot Number:	502955
Expiration:	25-May-2022
Platform(s):	HD-1/ HD-X



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

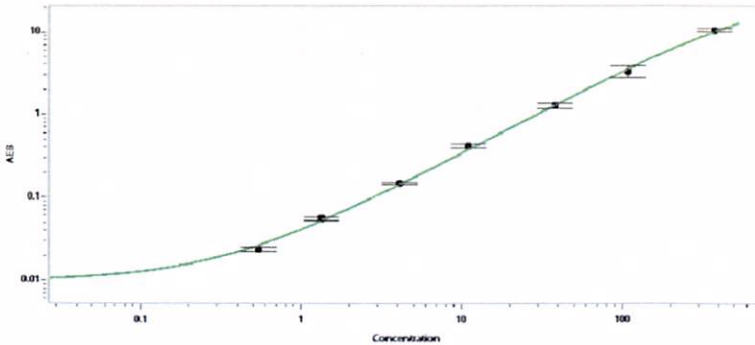
Component	Part Number	Lot Number
Bead Reagent	102246	114511
Detector Reagent	102248	114512
SBG Reagent	102250	114513
Sample Diluent	102252	114514
Calibrators	102255	115206
Control 1	102256	115207
Control 2	102257	115207
RGP Reagent	103159	N/A ¹

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	115207	4.03	3.22-5.02	3.18	2.08-4.77	
Control 2	115207	224	158-315	206	147-382	
Panel 1	008501	15.3	8.59-20.5	12.8	8.59-20.5	
Panel 2	008501	12.0	7.39-15.1	9.91	7.39-15.1	
Panel 3	008501	73.6	43.8-91.8	66.0	43.8-91.8	
Panel 4	009401	213	144-253	196	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	9.9
F	37.8	33.4
G	108	101
H	376	307

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 01 Nov 2021
Signature/Date