

Product Number:	103186
Lot Number:	502907
Expiration:	23-Apr-2022
Platform(s):	HD-1/ HD-X



Data below represents results generated on the Simoa HD-1® Analyzer.

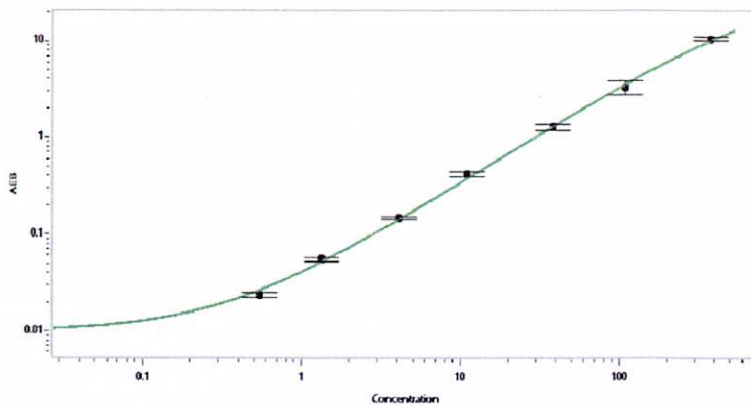
Component	Part Number	Lot Number
Bead Reagent	102246	111804
Detector Reagent	102248	111715
SBG Reagent	102250	111406
Sample Diluent	102252	111405
Calibrators	102255	112301
Control 1	102256	112302
Control 2	102257	112302
RGP Reagent	103159	N/A ¹

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	112302	3.69	2.77-4.15	3.27	1.84 - 4.07	
Control 2	112302	203	167-251	190	150 - 247	
Panel 1	008501	13.3	8.59-20.5	12.1	8.59 - 20.5	
Panel 2	008501	10.4	7.39-15.1	9.37	7.39 - 15.1	
Panel 3	008501	58.8	43.8-91.8	54.5	43.8 - 91.8	
Panel 4	009401	191	144-253	179	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	0.411	0.362
C	1.10	1.00
D	4.12	3.60
E	11.8	11.1
F	40.6	37.1
G	117	114
H	427	377

Review/Approval

Junpil Park
Name

QC Supervisor
Title


Signature/Date
6/11/2021