

Product Number:	103186
Lot Number:	502994
Expiration:	28-Jul-2022
Platform(s):	HD-1/ HD-X



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

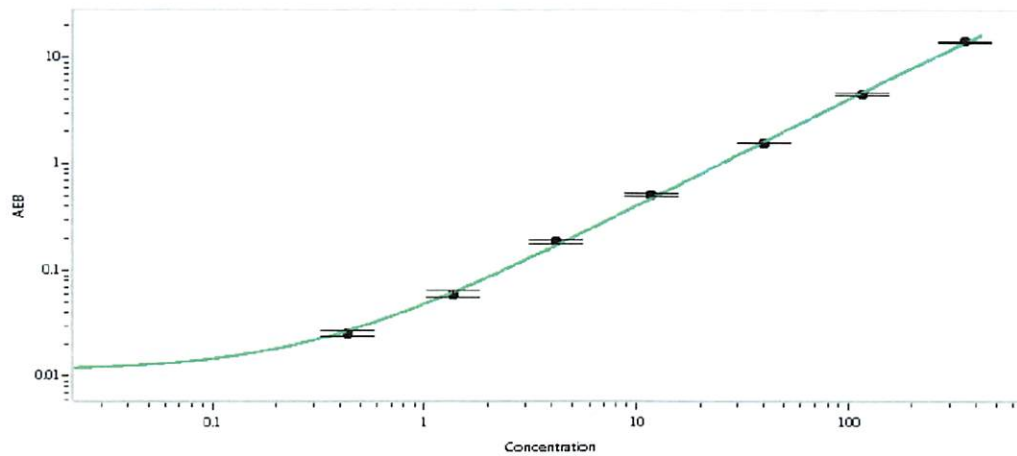
Component	Part Number	Lot Number
Bead Reagent	102246	121015
Detector Reagent	102248	121028
SBG Reagent	102250	121030
Sample Diluent	102252	121031
Calibrators	102255	121411
Control 1	102256	117908
Control 2	102257	117908
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	117908	4.63	3.00-5.12
Control 2	117908	261	198-297
Panel 1	008501	17.1	8.59-20.5
Panel 2	008501	13.4	7.39-15.1
Panel 3	008501	78.4	43.8-91.8
Panel 4	009401	249	128-302

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.431
C	1.36
D	4.18
E	11.7
F	39.7
G	117
H	354

Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

12 Aug 2021