

Product Number:	103186
Lot Number:	502976
Expiration:	17-May-2022
Platform(s):	HD-1/ HD-X



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

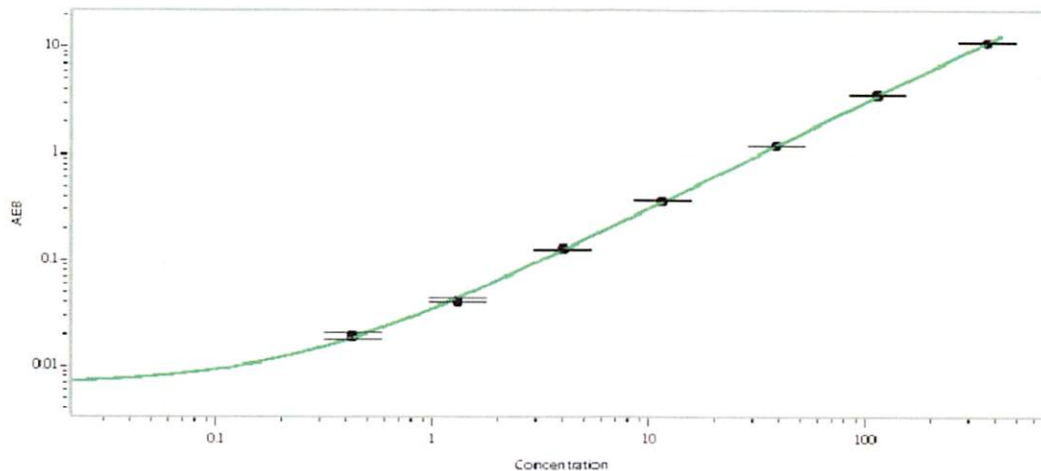
Component	Part Number	Lot Number
Bead Reagent	102246	115708
Detector Reagent	102248	115809
SBG Reagent	102250	115810
Sample Diluent	102252	115914
Calibrators	102255	115906
Control 1	102256	115907
Control 2	102257	115907
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	115907	3.87	3.06-5.03
Control 2	115907	241	140-314
Panel 1	008501	15.7	8.59-20.5
Panel 2	008501	12.1	7.39-15.1
Panel 3	008501	78.6	43.8-91.8
Panel 4	009401	275	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	0.425
C	1.30
D	3.97
E	11.4
F	38.9
G	113
H	364

Review/Approval Junpil Park QC Supervisor [Signature] 17 Jun 2021
 Name Title Signature/Date