

Product Number:	103186
Lot Number:	502700
Expiration:	15-Dec-2021
Platform(s):	HD-1/ HD-X



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

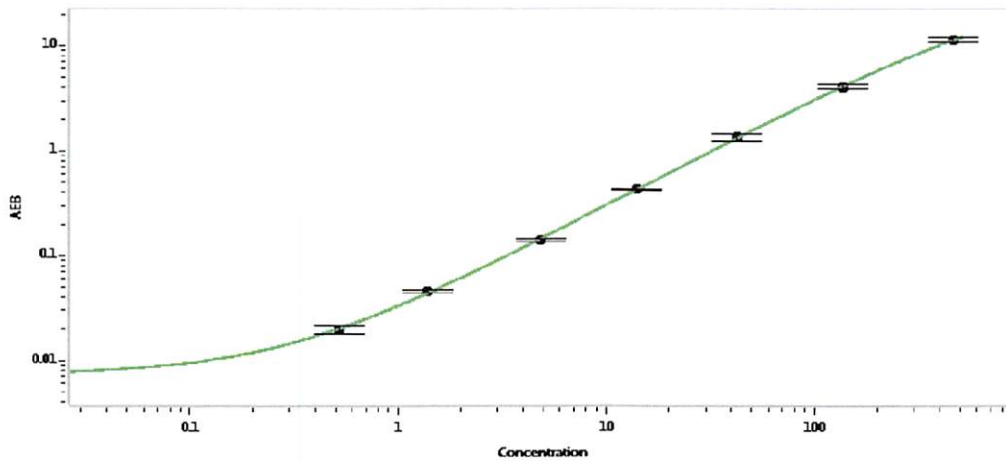
Component	Part Number	Lot Number
Bead Reagent	102246	101301
Detector Reagent	102248	101516
SBG Reagent	102250	102301
Sample Diluent	102252	102004
Calibrators	102255	102101
Control 1	102256	102203
Control 2	102257	102203
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	102203	4.22	2.62 - 6.02
Control 2	102203	144	123 - 184
Panel 1	922601	18.3	14.3 - 21.4
Panel 2	922602	7.08	4.97 - 7.45
Panel 3	922603	41.2	28.8 - 43.2
Panel 4	922604	182	133 - 201

* 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	0.519
C	1.37
D	4.87
E	14.0
F	42.3
G	136
H	466

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 28 Jan 2021
Signature/Date