

Product Number:	102184
Lot Number:	503013
Expiration:	24-Feb-2022
Platform(s):	HD-1/ HD-X



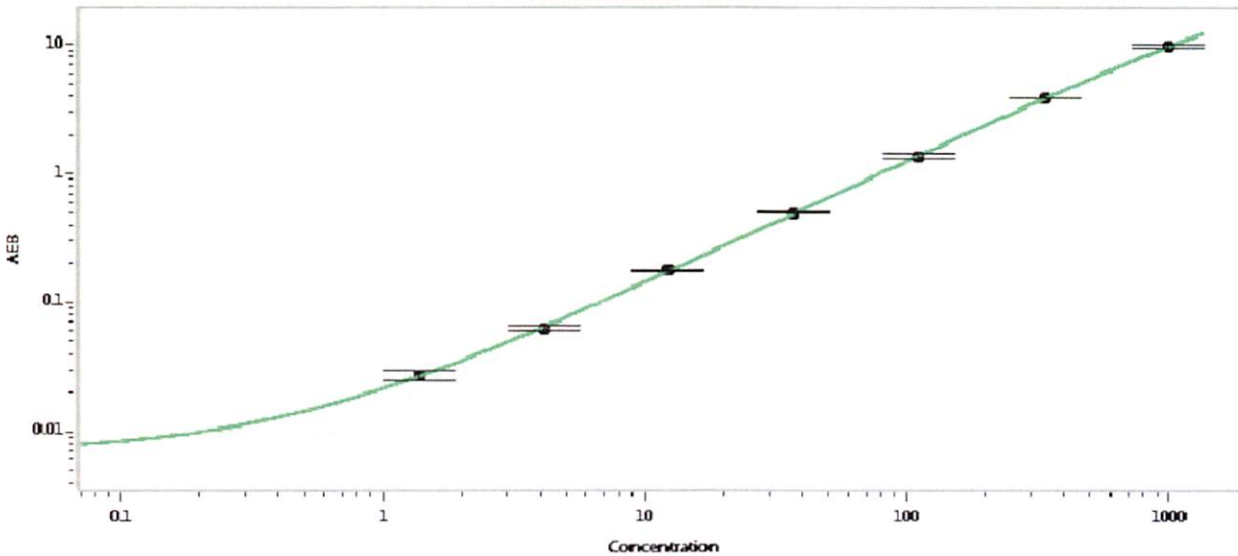
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102176	117305
Detector Stock	102177	117606
Calibrator Concentrate	102180	005605
Calibrator Diluent	102178	112914
Sample Diluent	102179	109025

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	40 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 02 Jul 2021
Signature/Date