

Product Number:	102343
Lot Number:	502825
Expiration:	18-Nov-2021
Platform(s):	HD-1/HD-X



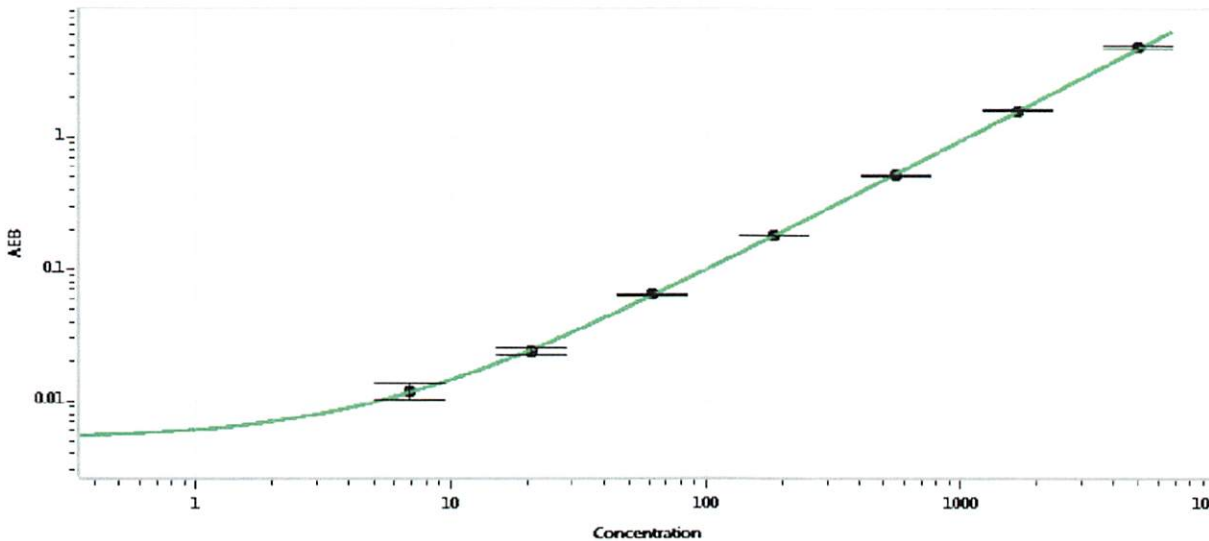
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102337	107903
Detector Stock	102338	108304
Calibrator Concentrate	102342	024404
Calibrator Diluent	102339	107521
Sample Diluent	102340	107909

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	100 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.00
B	6.86
C	20.6
D	61.7
E	185
F	556
G	1667
H	5000

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 07Apr2021
Signature/Date