

Product Number:	102100
Lot Number:	503003
Expiration:	10-Apr-2022
Platform(s):	HD-1 / HD-X



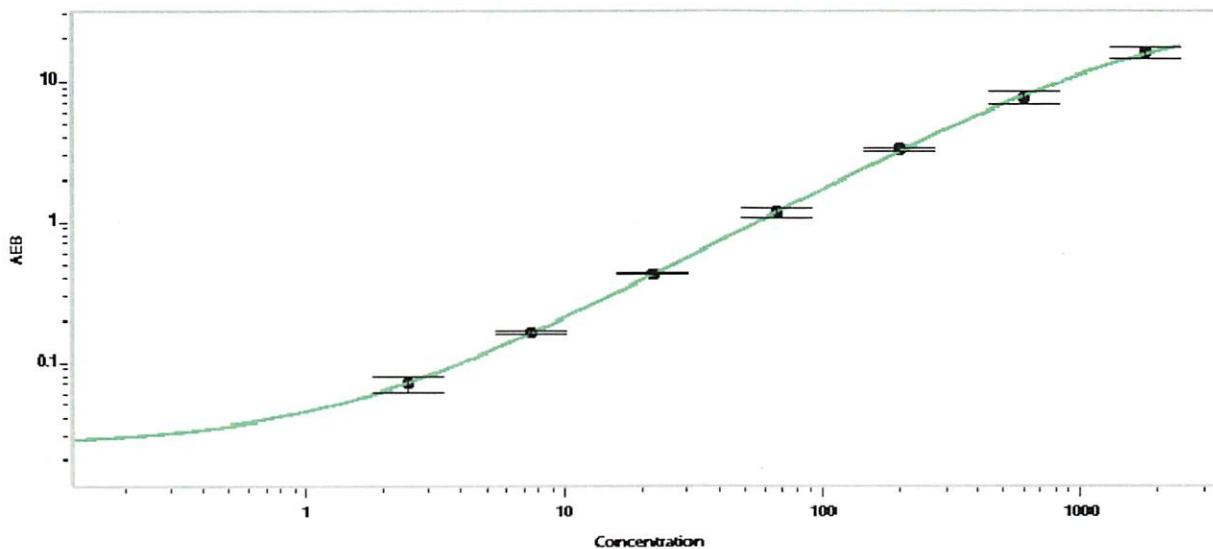
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102092	116510
Detector Stock	102093	116811
Calibrator Concentrate	102096	117407
Calibrator Diluent	102094	112721
Sample Diluent	102095	112722

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	36 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.47
C	7.41
D	22.2
E	66.7
F	200
G	600
H	1800

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 28 Jun 2021
Signature/Date