

Product Number:	102475
Lot Number:	503157
Expiration:	14-Feb-2022
Platform(s):	HD-1/HD-X



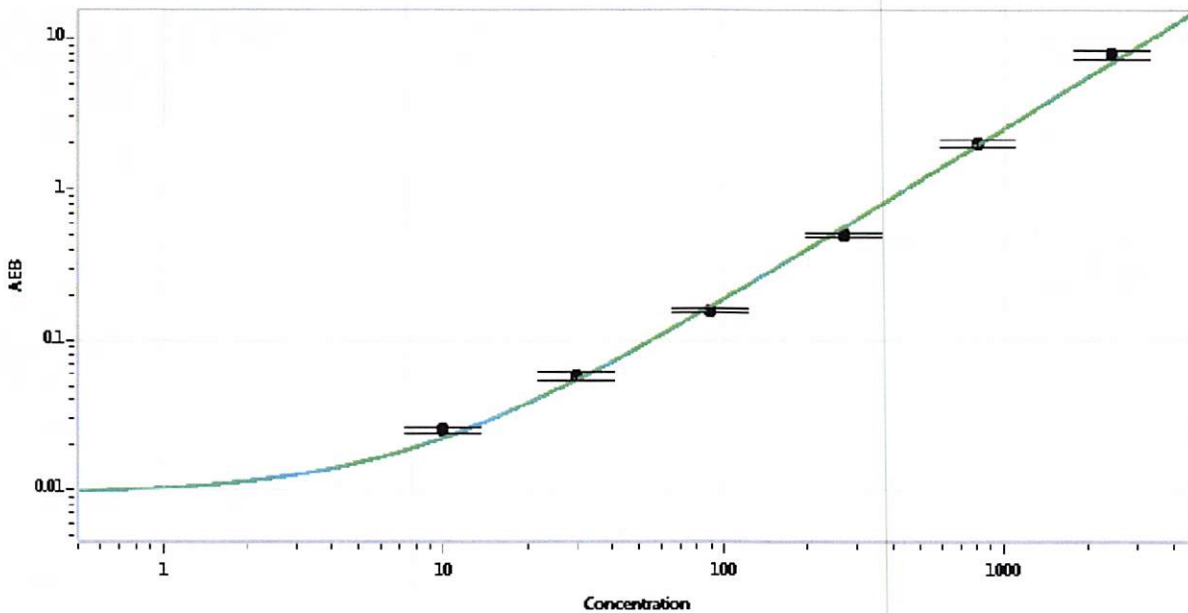
Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102468	127001
Detector Stock	102469	127008
Calibrator Concentrate	102472	004511
Calibrator Diluent	102470	126808
Sample Diluent	102471	121114

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	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.88
C	29.6
D	88.9
E	267
F	800
G	2400

Review/Approval

Jungil Park
Name

QC supervisor
Title

[Signature] 08 Oct 2021
Signature/Date