

Product Number:	102039
Lot Number:	503047
Expiration:	23-Apr-2022
Platform(s):	HD-1/HD-X & SR-X



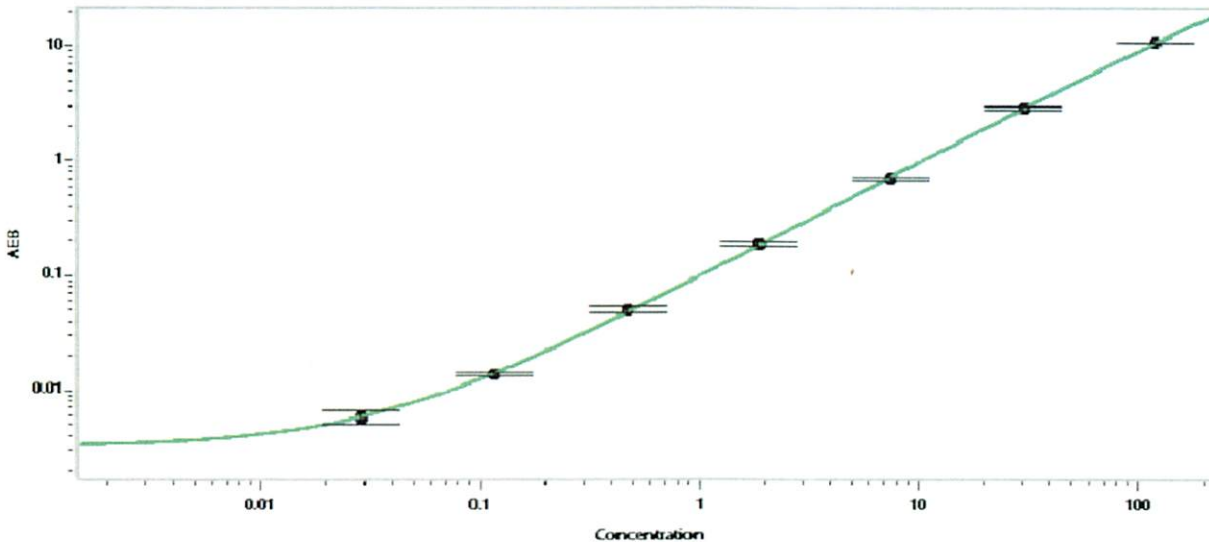
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102028	118915
Detector Stock	102029	119203
Calibrator Concentrate	102032	119607
Calibrator Diluent	102030	118023
Sample Diluent	103402	111918

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

Junpil Park

QC Supervisor

 26 Jul 2021

Name

Title

Signature/Date