

Product Number:	101863
Lot Number:	502941
Expiration:	17-Mar-2022
Platform(s):	HD-1/HD-X



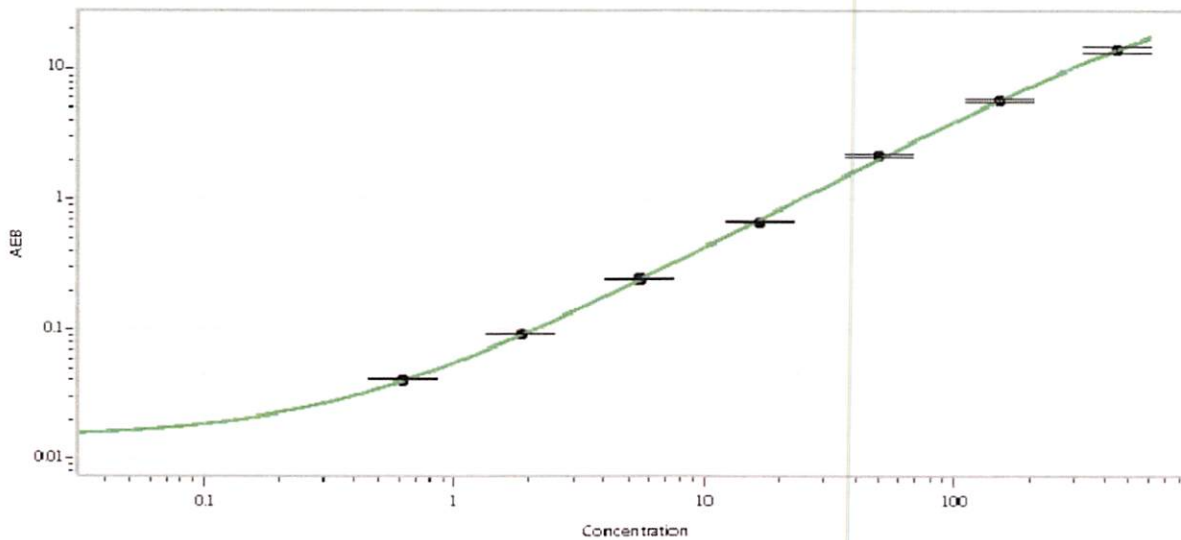
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101904	113708
Detector Stock	101905	113911
Calibrator Concentrate	101858	018202
Calibrator Diluent	101990	113312
Sample Diluent	101991	113809

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	9 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450

Review/Approval

Junpil Park

QC Supervisor

[Signature] 04 Jun 2021

Name

Title

Signature/Date