

Product Number:	102794
Lot Number:	502743
Expiration:	14-Nov-2021
Platform(s):	HD-1/HD-X



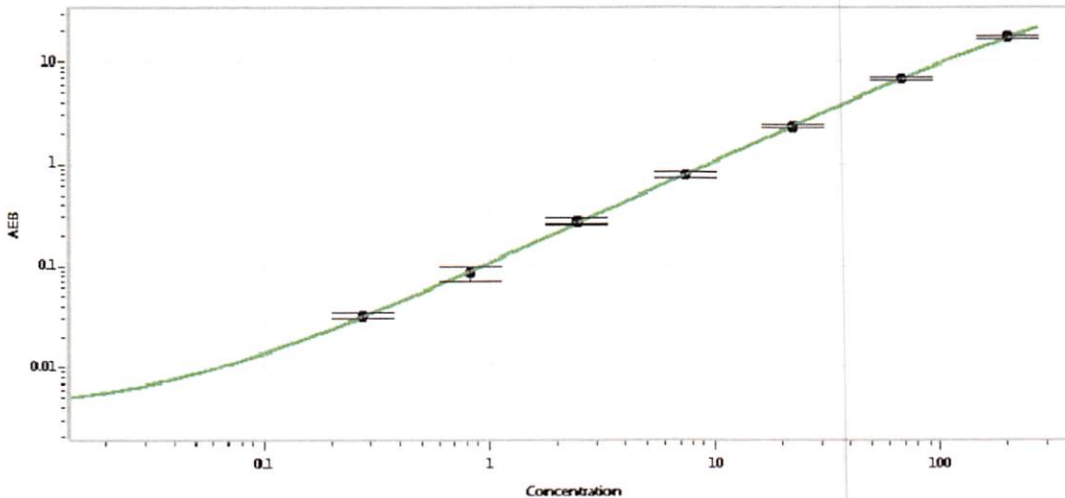
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102787	104205
Detector Stock	102788	104330
Calibrator Concentrate	102791	936007
Calibrator Diluent	102789	102921
Sample Diluent	102790	104213

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	4 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Junpil Park  
Name

QC Supervisor  
Title

*[Signature]* 04 Mar 2021  
Signature/Date