

Product Number:	101984
Lot Number:	502886
Expiration:	26-Jan-2022
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	101910	107908
Detector Stock	101911	110608
Calibrator Concentrate	101983	108306
Calibrator Diluent	101908	108005
Sample Diluent	101909	108001

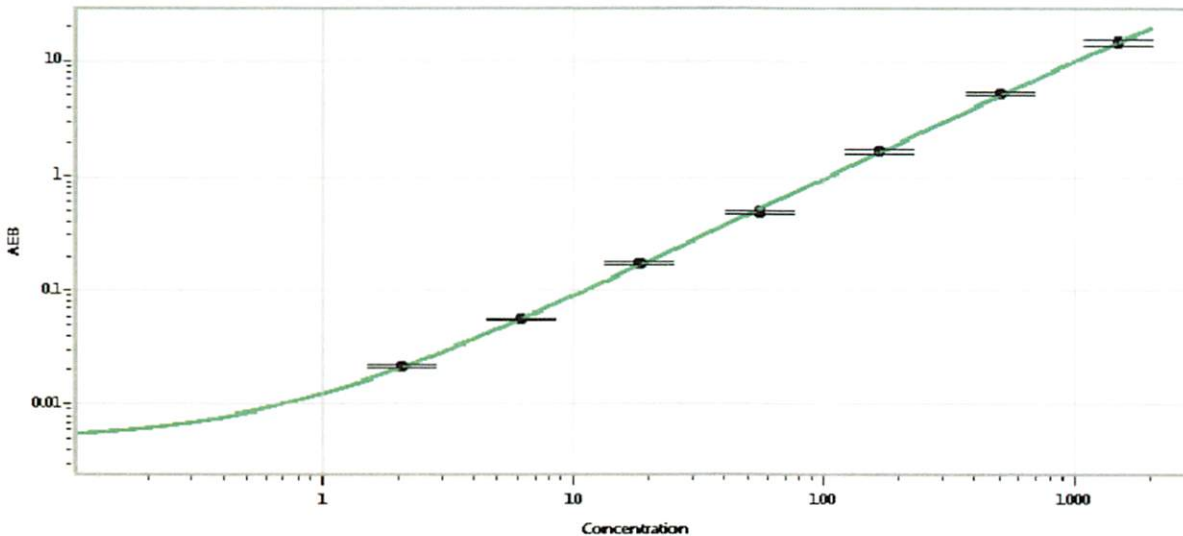


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	30 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.06
C	6.17
D	18.5
E	55.6
F	167
G	500
H	1500

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 26 Apr 2021
Signature/Date