

Product Number:	101863
Lot Number:	502449
Expiration:	31-Mar-2021
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	101904	018205
Detector Stock	101905	018101
Calibrator Concentrate	101858	018202
Calibrator Diluent	101990	009105
Sample Diluent	101991	018106

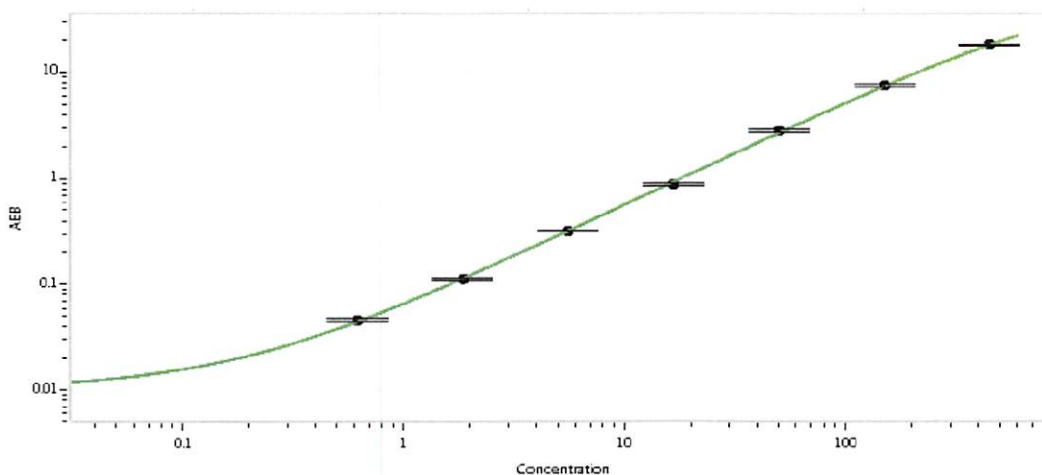
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

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

Shivani Goel

Quality Control Manager

Shivani Goel 15/01/2020

Name

Title

Signature/Date