

Product Number:	101968
Lot Number:	502924
Expiration:	7-Jan-2022
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	101966	112101
Detector Stock	101967	112304
Calibrator Concentrate	101965	000719
Calibrator Diluent	101963	111724
Sample Diluent	101964	111408

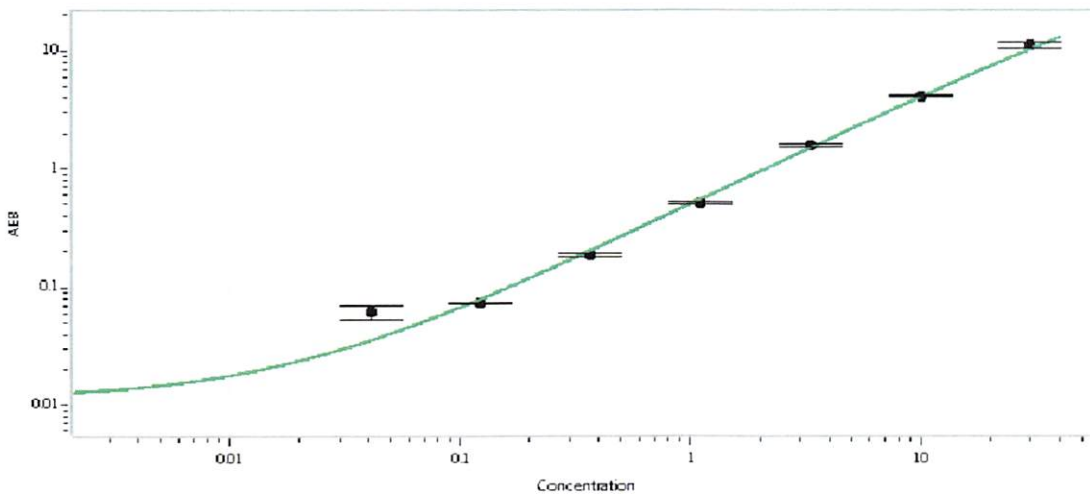


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	600 pg/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Junpil Park

QC Supervisor

02 Jun 2021

Name

Title

Signature/Date