

Product Number:	101863
Lot Number:	503735
Expiration:	25-Oct-2023
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	101904	231203
Detector Stock	101905	231412
Calibrator Concentrate	101858	231307
Calibrator Diluent	101990	231315
Sample Diluent	101991	231304

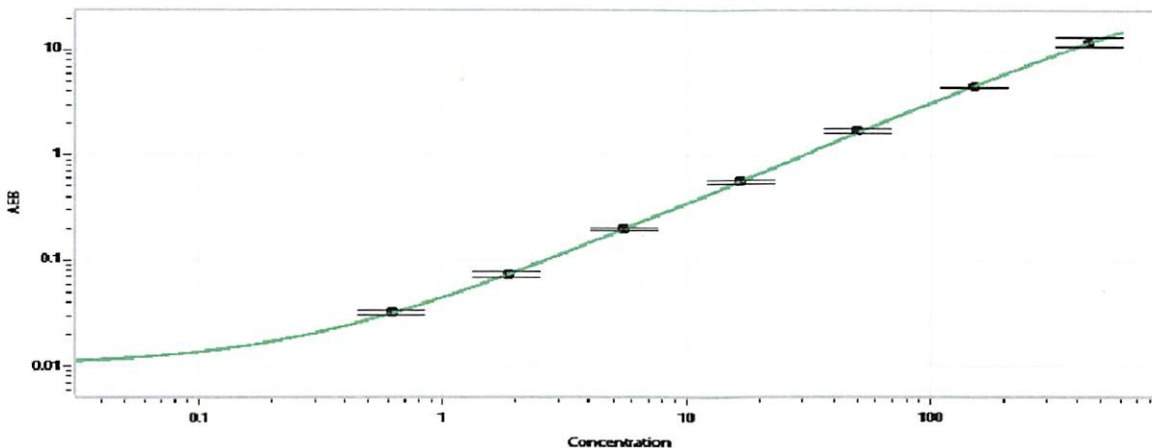


Data below represents results generated on the Simoa® HD-X Analyzer.

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 (ng/mL)	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: Tony Iem Name: QC Title: [Signature] Signature/Date: 17NOV2022