

Product Number:	102318
Lot Number:	503418
Expiration:	12-Dec-2022
Platform(s):	HD-1/HDX



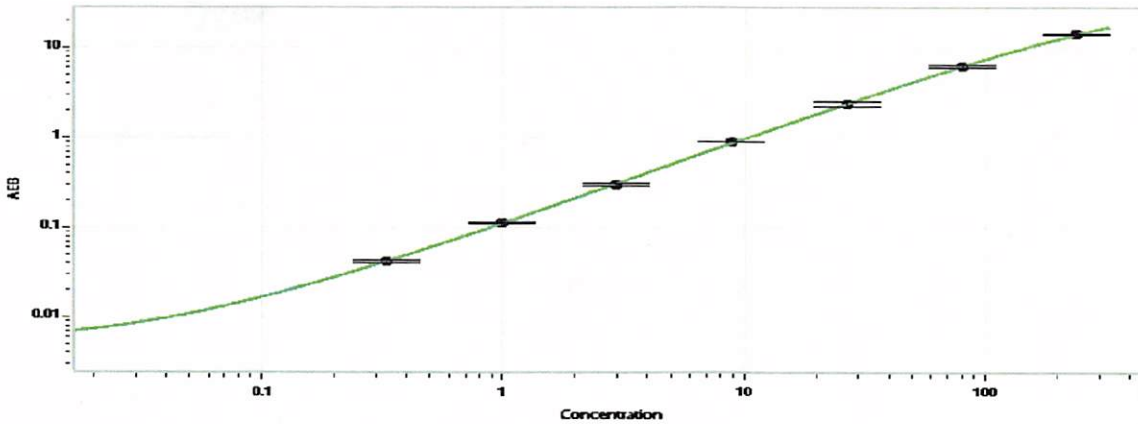
Component	Part Number	Lot Number
Bead Stock	102309	209604
Detector Stock	102310	210112
Calibrator Concentrate	102314	119507
Calibrator Diluent	102311	209713
Sample Diluent	102312	209712

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 / HD-X Analyzer.

Calibrator Concentrate Target Value	4800 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

Review/Approval

Richa Shrestha

QC Supervisor

[Signature] 03 May 2022

Name

Title

Signature/Date