

Product Number:	102570
Lot Number:	503242
Expiration:	13-Jul-2022
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102563	133205
Detector Stock	102564	133710
Detector Diluent	102590	132510
Calibrator Concentrate	102567	133306
Calibrator Diluent	102565	121119
Sample Diluent	102566	121120

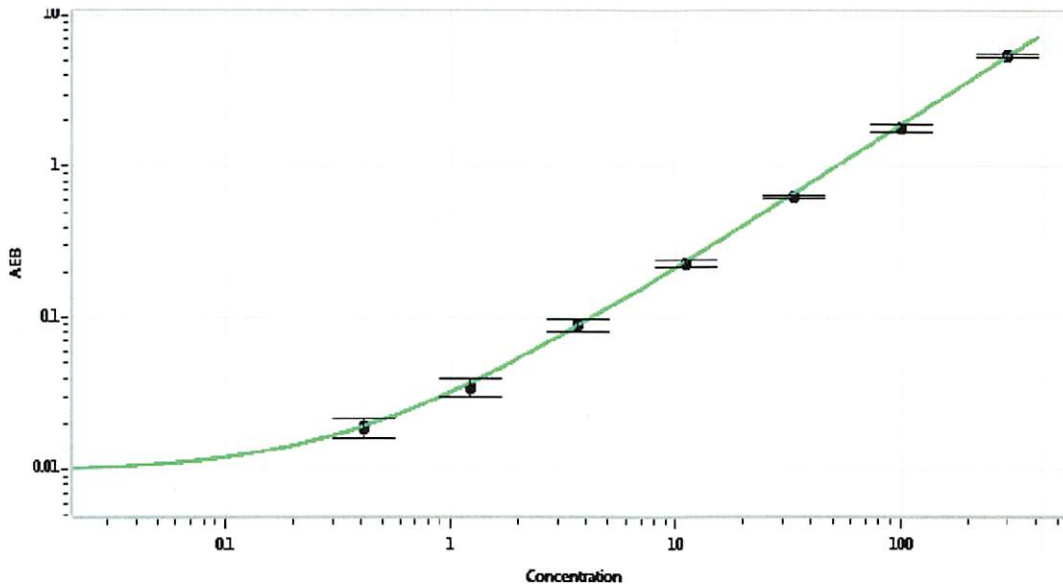


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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

Calibrator Concentrate Target Value	6 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

Wenkei Tsai QC Supervisor [Signature] 17 Dec 2021

Name Title Signature/Date