

Product Number:	102570
Lot Number:	502755
Expiration:	23-Nov-2021
Platform(s):	HD-1/HD-X

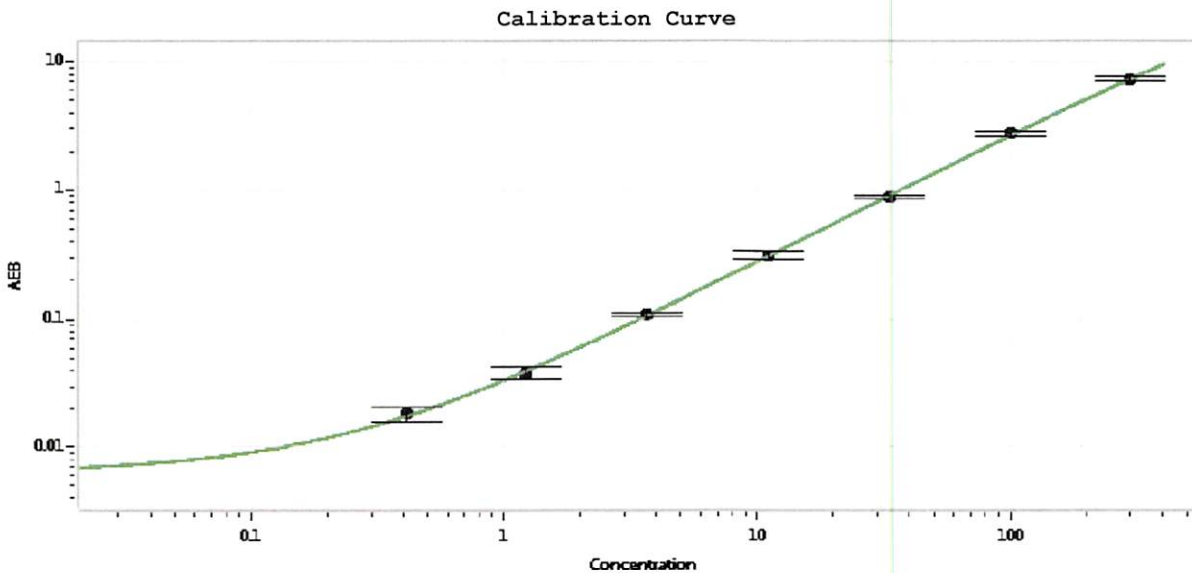


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102563	104402
Detector Stock	102564	105401
Detector Diluent	102590	104407
Calibrator Concentrate	102567	104326
Calibrator Diluent	102565	036322
Sample Diluent	102566	036321

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
-------------------------------------	---------



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

Junpil Park
Name

QC Supervisor
Title

[Signature] 07 Apr 2021
Signature/Date