

Product Number:	102064
Lot Number:	502876
Expiration:	26-Jan-2022
Platform(s):	HD-1/HD-X



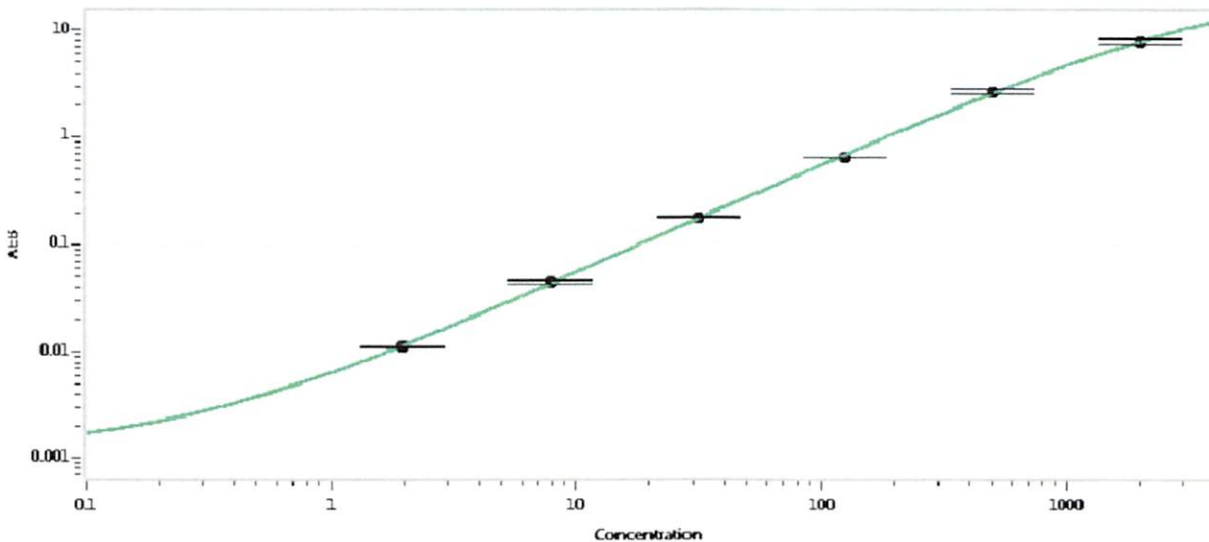
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102056	110108
Detector Stock	102057	110216
Calibrator Concentrate	102060	110408
Calibrator Diluent	102058	110214
Sample Diluent	102059	110213

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	40 ng/mL
--	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.95
C	7.81
D	31.3
E	125
F	500
G	2000

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 26 Apr 2021
Signature/Date