

Product Number:	102233
Lot Number:	503031
Expiration:	6-Jan-2022
Platform(s):	HD-1/HD-X, SRX

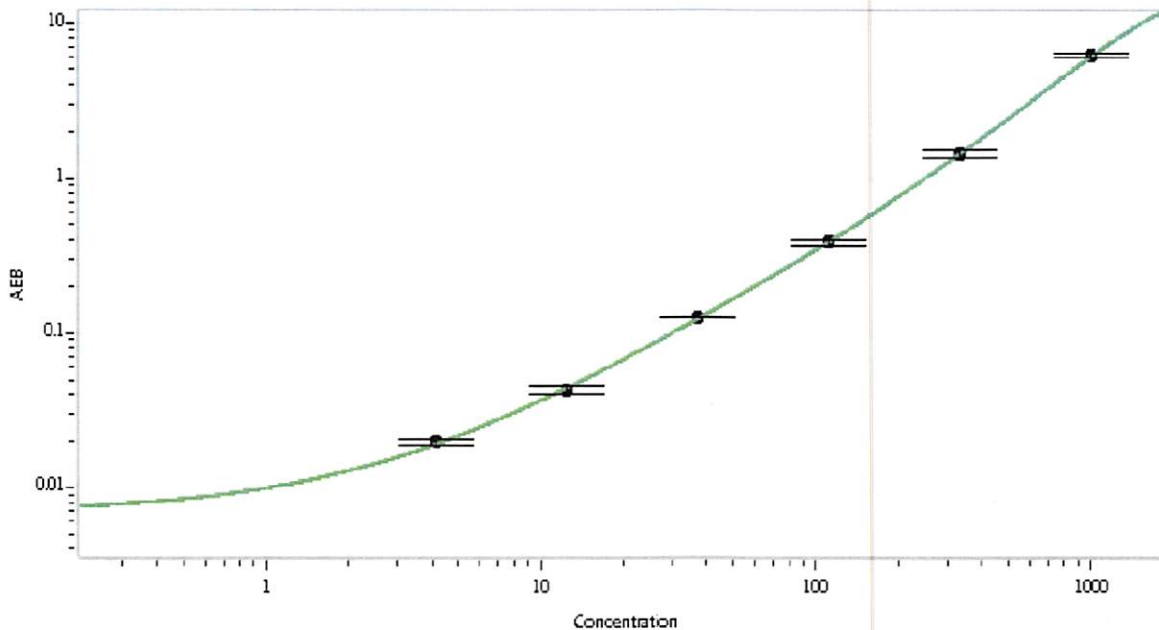


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	118704
Detector Stock	102224	119407
Calibrator Concentrate	102229	120404
Calibrator Diluent	102225	118811
Sample Diluent	102226	118809

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	20 ng/mL
-------------------------------------	----------



Calibrator Levels (pg/mL)	
A	0.000
B	4.12
C	12.3
D	37.0
E	111
F	333
G	1000

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 06 Aug 2021
Signature/Date