

Product Number:	102233
Lot Number:	503111
Expiration:	18-Feb-2022
Platform(s):	HD-1/HD-X, SRX



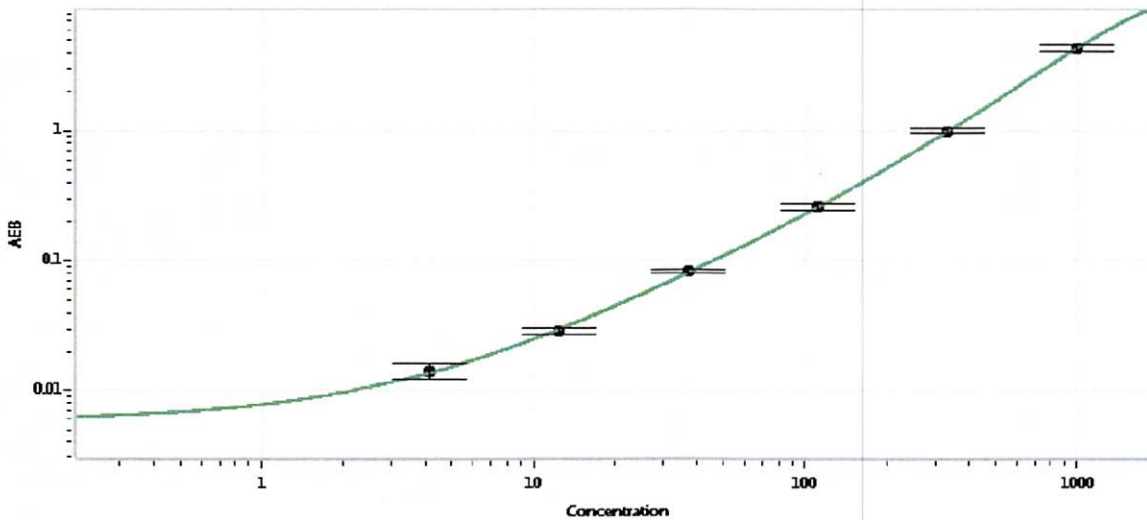
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	123207
Detector Stock	102224	123008
Calibrator Concentrate	102229	120404
Calibrator Diluent	102225	123209
Sample Diluent	102226	123208

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.11
C	12.3
D	37.0
E	111
F	333
G	1000

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 4 Sep 2021
Signature/Date