

Product Number:	102233
Lot Number:	502980
Expiration:	21-Oct-2021
Platform(s):	HD-1/HD-X, SRX



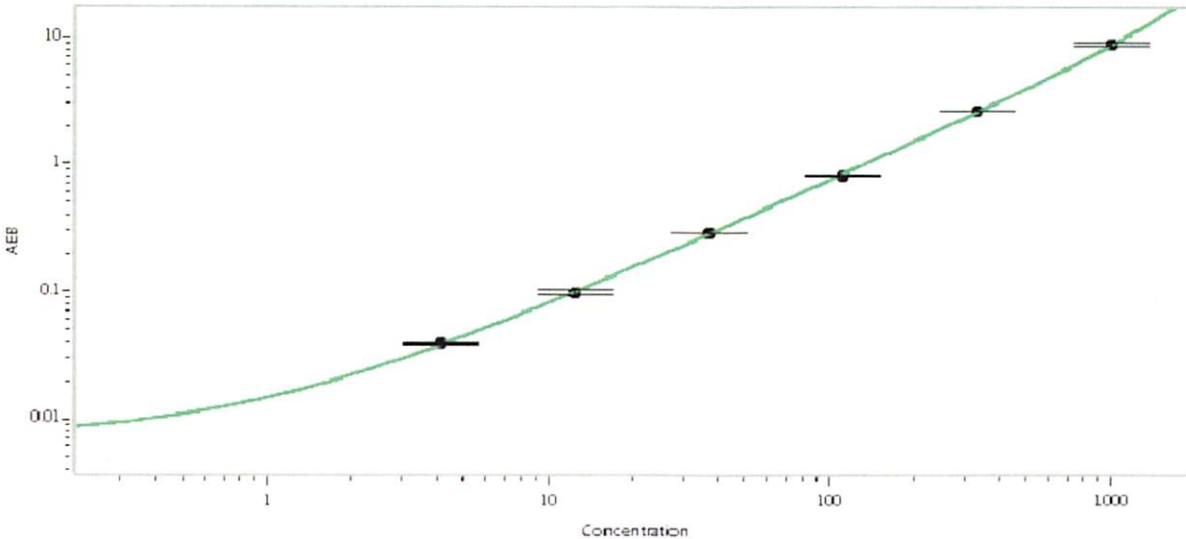
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	111105
Detector Stock	102224	115815
Calibrator Concentrate	102229	027601
Calibrator Diluent	102225	103310
Sample Diluent	102226	103309

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	4.12
C	12.3
D	37.0
E	111
F	333
G	1000

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 11 Jun 2021
Signature/Date