

Product Number:	102336
Lot Number:	503138
Expiration:	1-Jun-2022
Platform(s):	HD-1 / HD-X

Component	Part Number	Lot Number
Bead Stock	102330	115508
Detector Stock	102331	115822
Calibrator Concentrate	102335	125608
Calibrator Diluent	102332	115220
Sample Diluent	102333	115821

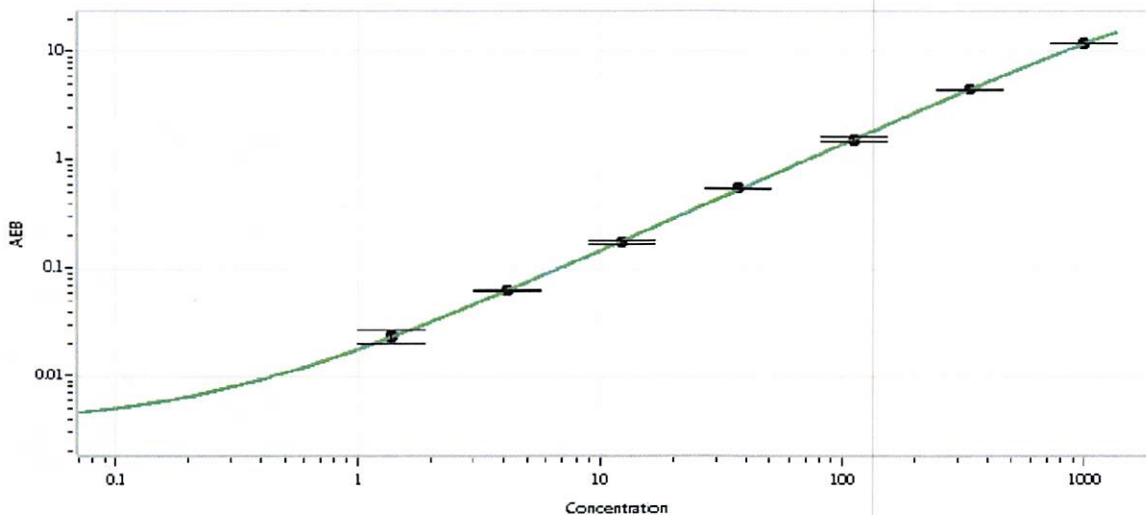
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent	101954	N/A
Discovery Detector Diluent	101955	N/A
Discovery SBG Diluent	101956	N/A
Discovery SBG Concentrate	103297	N/A
RGP Reagent	103159	N/A



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

Calibrator Concentrate Target Value	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Jungil Park
Name

QC Supervisor
Title

[Signature] 21 Sep 2021
Signature/Date