

Product Number:	102336
Lot Number:	502801
Expiration:	14-Nov-2021
Platform(s):	HD-1/HD-X



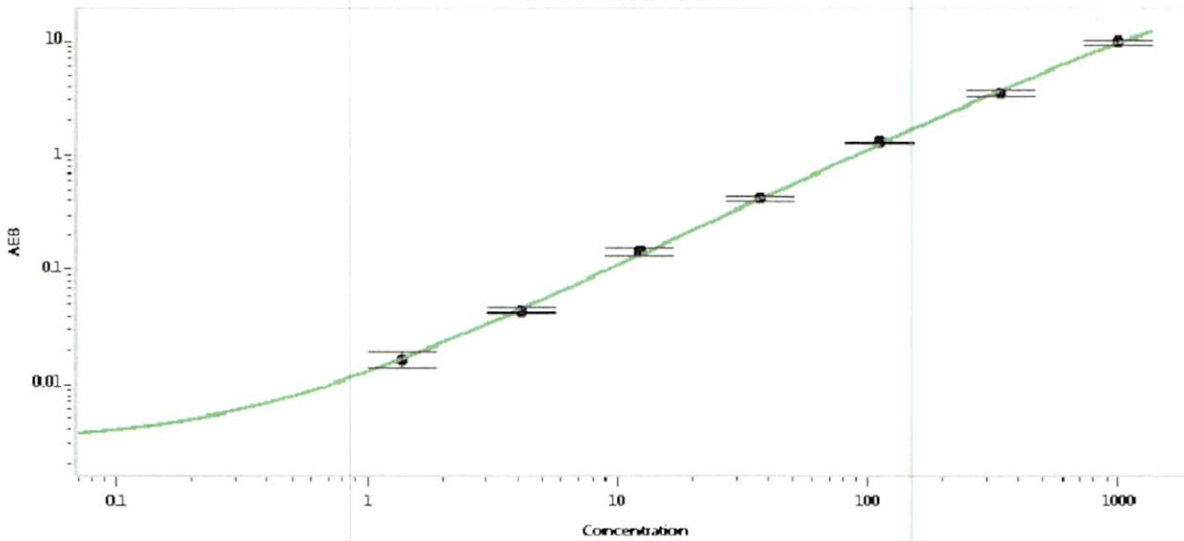
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	102908
Detector Stock	102331	102916
Calibrator Concentrate	102335	106815
Calibrator Diluent	102332	102920
Sample Diluent	102333	102917

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	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date 18 Mar 2021