

Product Number:	102336
Lot Number:	502862
Expiration:	4-Mar-2022
Platform(s):	HD-1/HD-X



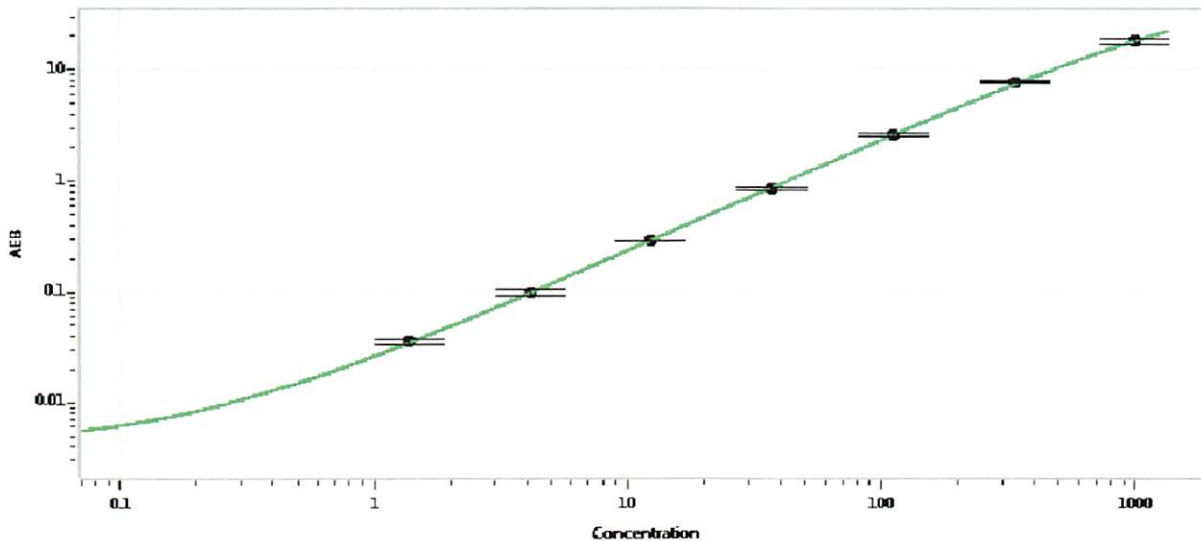
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	111401
Detector Stock	102331	111402
Calibrator Concentrate	102335	108915
Calibrator Diluent	102332	111404
Sample Diluent	102333	111403

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] Signature/Date: 06 May 2021