

Product Number:	102570
Lot Number:	502931
Expiration:	24-Feb-2022
Platform(s):	HD-1/HD-X



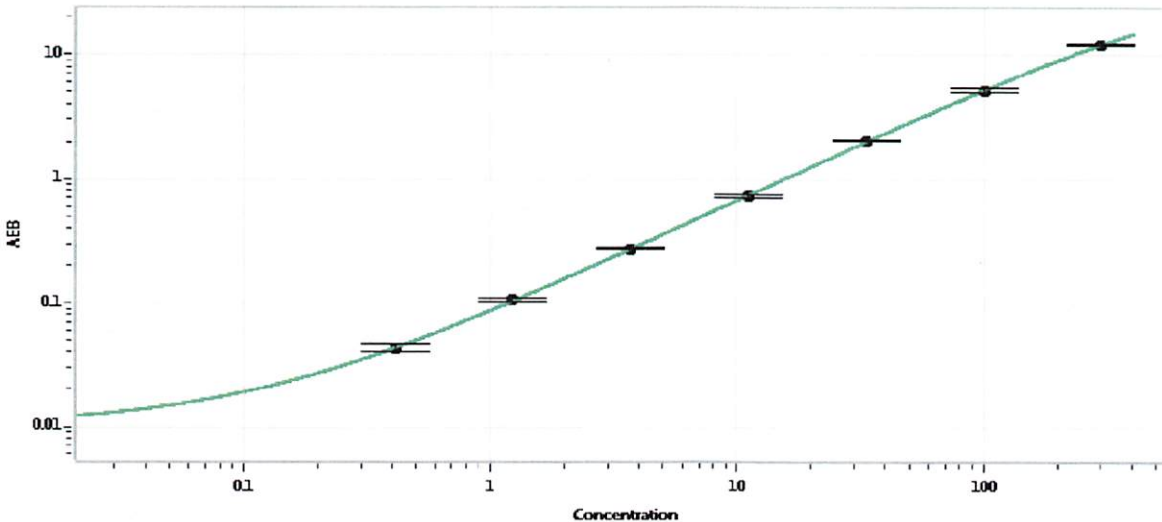
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102563	112801
Detector Stock	102564	112711
Detector Diluent	102590	113018
Calibrator Concentrate	102567	104326
Calibrator Diluent	102565	111911
Sample Diluent	102566	105521

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	6 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

Junpil Park  
Name

QC Supervisor  
Title

*[Signature]* 03 Jun 20 21  
Signature/Date