

Product Number:	102570
Lot Number:	503242
Expiration:	14-Jun-2022
Platform(s):	HD-1/HD-X



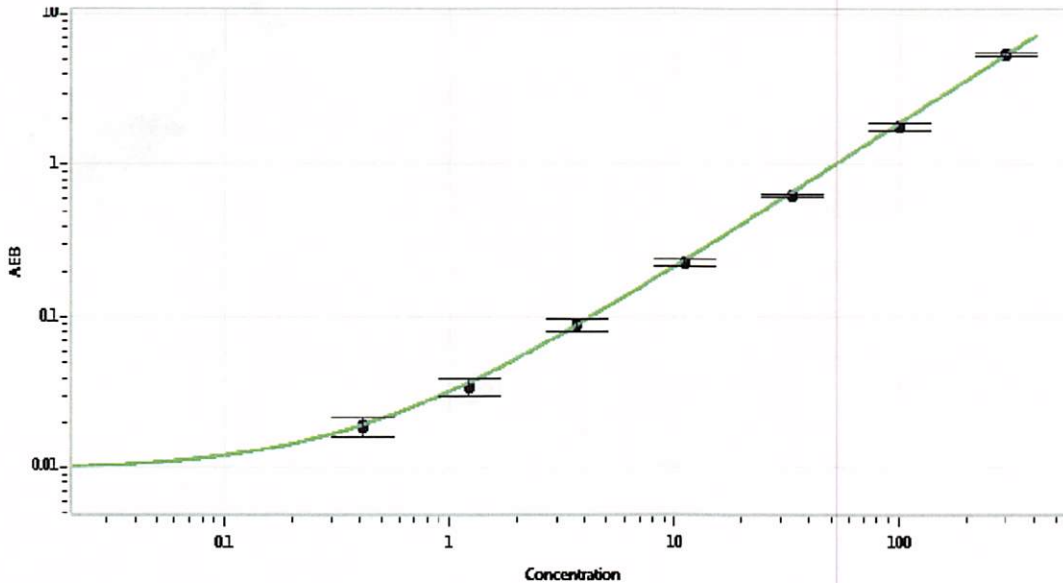
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102563	133205
Detector Stock	102564	133710
Detector Diluent	102590	132510
Calibrator Concentrate	102567	133306
Calibrator Diluent	102565	121119
Sample Diluent	102566	121120

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] 21 Dec 2021
Signature/Date