

Product Number:	102556
Lot Number:	502882
Expiration:	26-Jan-2022
Platform(s):	HD-1 / HD-X



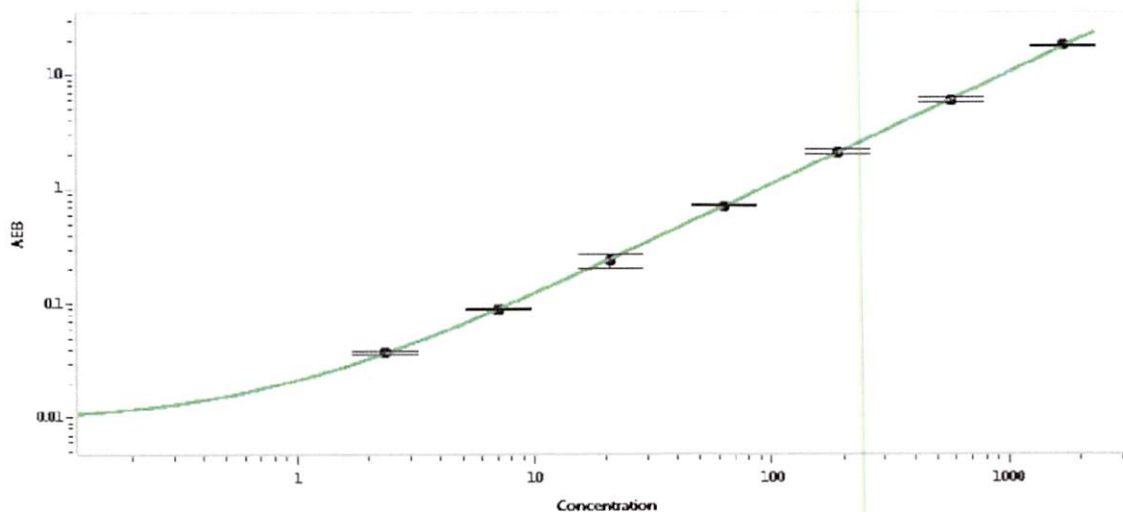
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102549	110305
Detector Stock	102550	110309
Calibrator Concentrate	102553	109012
Calibrator Diluent	102551	106321
Sample Diluent	102552	110011

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	40 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.33
C	7.00
D	21.0
E	63.0
F	189
G	567
H	1700

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 26 Apr 2021
Signature/Date