

Product Number:	102100
Lot Number:	502621
Expiration:	16-Apr-2021
Platform(s):	HD-1 / HD-X



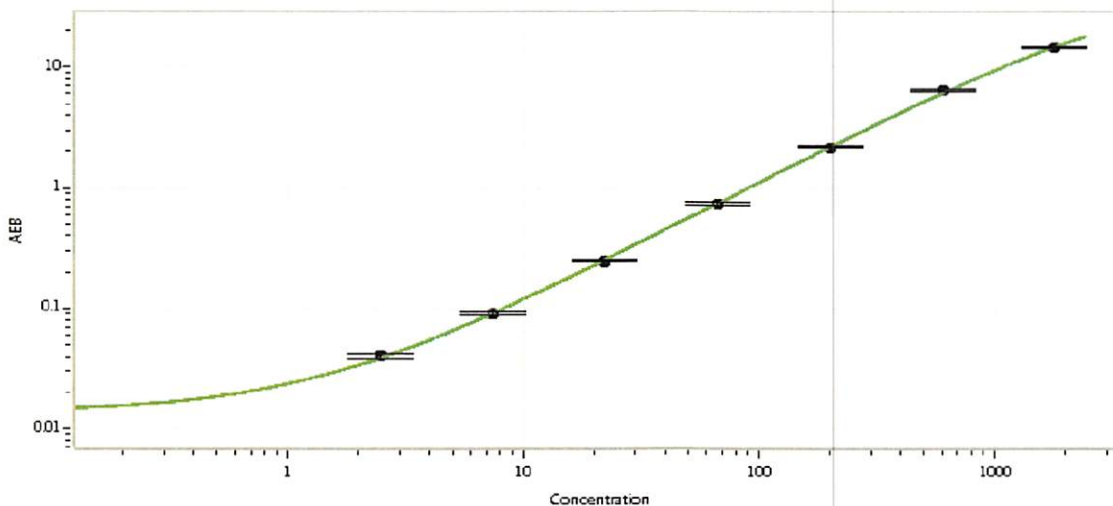
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102092	011703
Detector Stock	102093	011807
Calibrator Concentrate	102096	919111
Calibrator Diluent	102094	010713
Sample Diluent	102095	010714

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	36 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.47
C	7.41
D	22.2
E	66.7
F	200
G	600
H	1800

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 04 Dec 2020
Signature/Date