

Product Number:	102635
Lot Number:	502920
Expiration:	4-Mar-2022
Platform(s):	HD-1/ HD-X



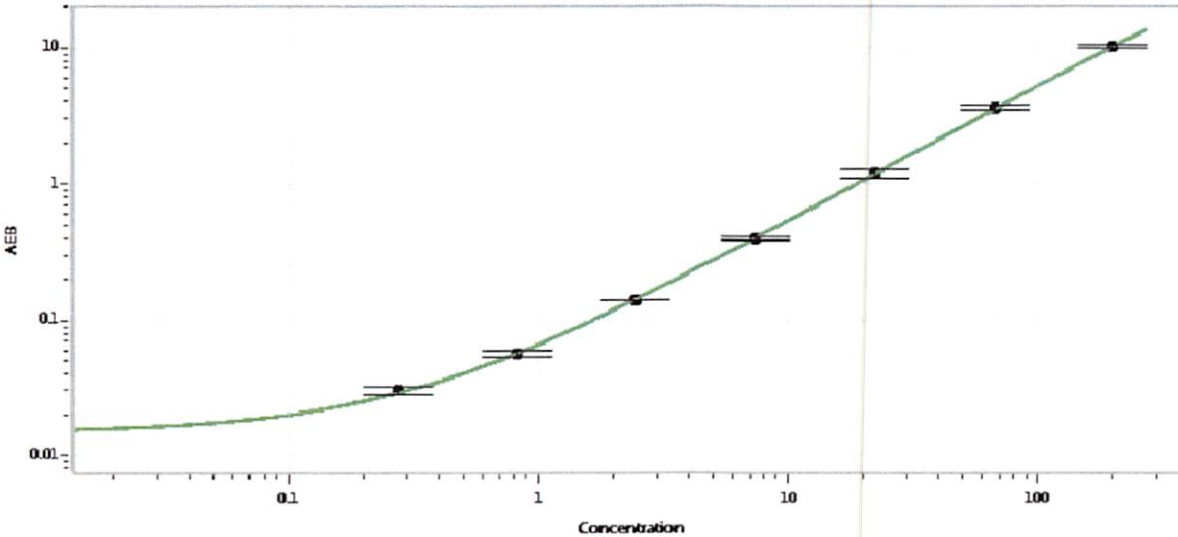
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102628	108602
Detector Stock	102629	113012
Calibrator Concentrate	102632	013308
Calibrator Diluent	102630	106319
Sample Diluent	102631	108707

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	4 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 19 May 2021
Signature/Date