

Product Number:	102209
Lot Number:	502883
Expiration:	4-Mar-2022
Platform(s):	HD-1/HD-X & SR-X



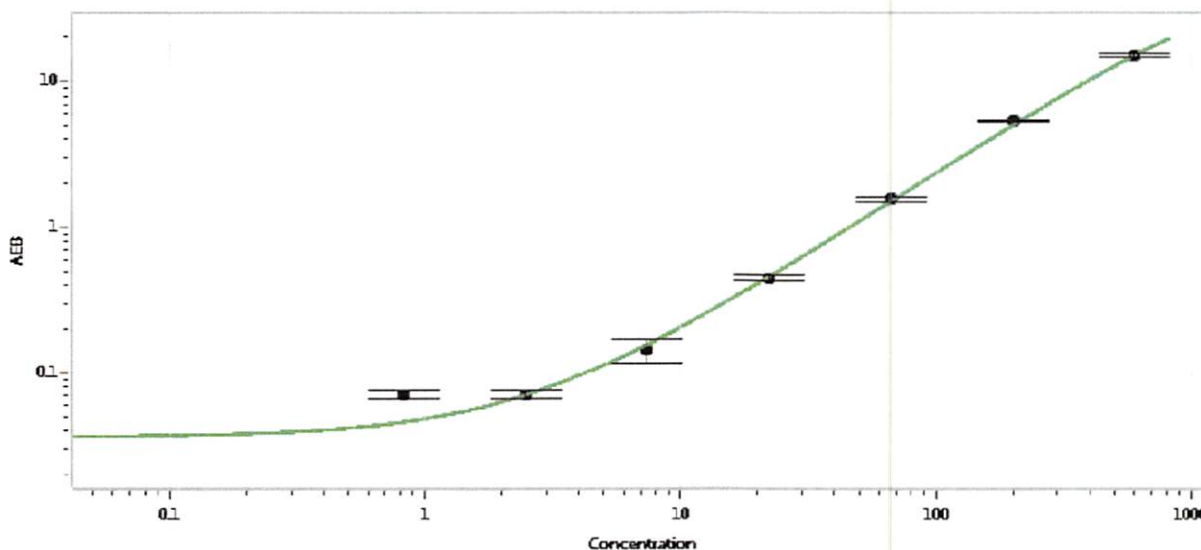
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102201	110912
Detector Stock	102202	110915
Calibrator Concentrate	102205	023103
Calibrator Diluent	102203	110918
Sample Diluent	102204	110917

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	12 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval

Junpil Park QC Supervisor *[Signature]* 18 May 2021
Name Title Signature/Date