

Product Number:	101824
Lot Number:	503161
Expiration:	28-Jun-2022
Platform(s):	HD-1/ HD-X



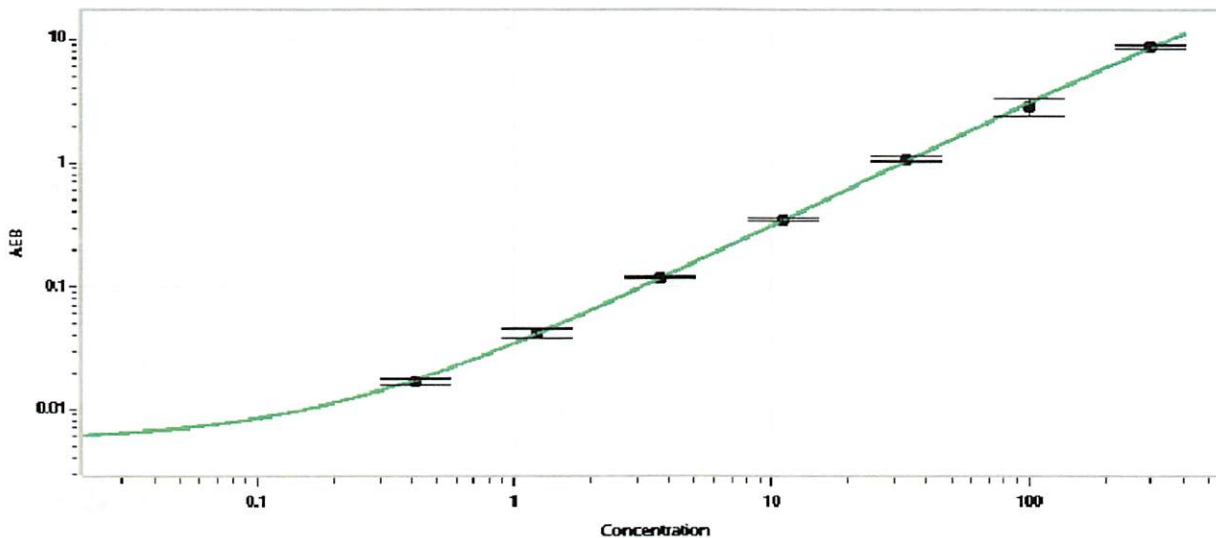
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101929	127103
Detector Stock	101930	127106
Calibrator Concentrate	101819	127003
Calibrator Diluent	101943	124911
Sample Diluent	101944	117925

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	6000 pg/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] Signature/Date, 13 Oct 2021 Date