

Product Number:	102462
Lot Number:	503147
Expiration:	17-Jul-2022
Platform(s):	HD-1/ HD-X



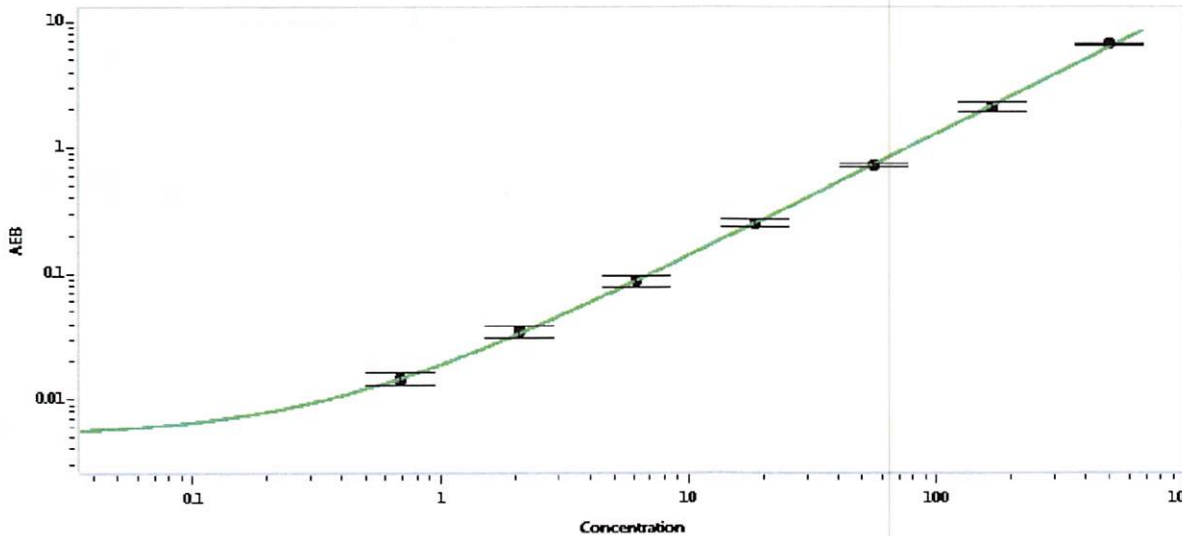
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102455	119803
Detector Stock	102456	126404
Calibrator Concentrate	102459	120304
Calibrator Diluent	102457	120010
Sample Diluent	102458	120026

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	10 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.686
C	2.06
D	6.17
E	18.5
F	55.6
G	167
H	500

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 20220924
Signature/Date