

Product Number:	102504
Lot Number:	502672
Expiration:	14-Sep-2021
Platform(s):	HD-1



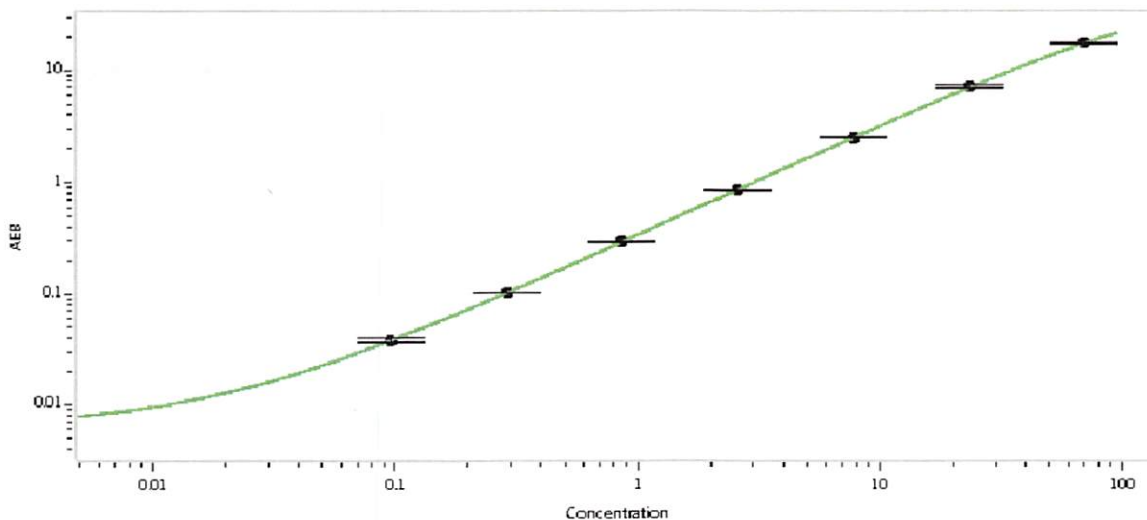
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102497	035703
Detector Stock	102498	035706
Calibrator Concentrate	102501	036602
Calibrator Diluent	102499	035804
Sample Diluent	102500	035803

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	1.4 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.096
C	0.288
D	0.864
E	2.59
F	7.77
G	23.3
H	70.0

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 15Jan2021

Name

Title

Signature/Date