

Product Number:	102700
Lot Number:	501698
Expiration:	22-Jan-2020
Platform(s):	HD-1



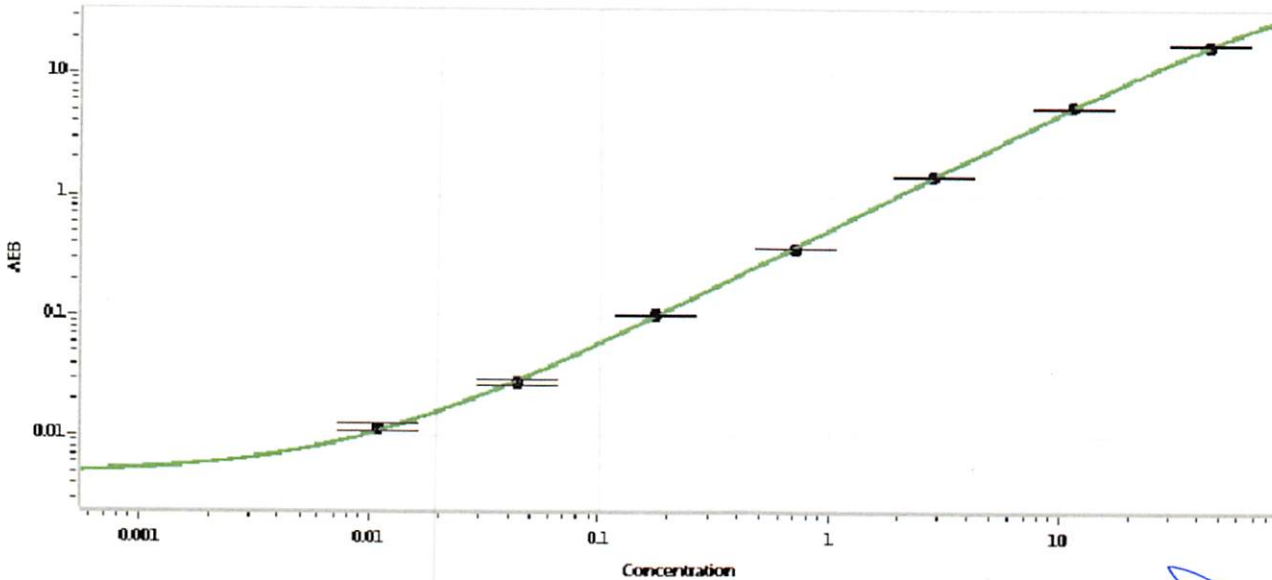
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102693	903205
Detector Stock	102694	903107
Calibrator Concentrate	102697	MF-060418-01
Calibrator Diluent	102695	902922
Sample Diluent	102696	902921

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.011
C	0.044
D	0.176
E	0.703
F	2.81
G	11.3
H	45.0

Review/Approval

*Junpil Park
Name

QC Analyst II
Title

[Signature] 06 Mar 2019
Signature/Date

* On behalf of Patrizia Stadler Assoc. Director, QC