

Product Number:	102109
Lot Number:	501678
Expiration:	5-Jan-2020
Platform(s):	HD-1



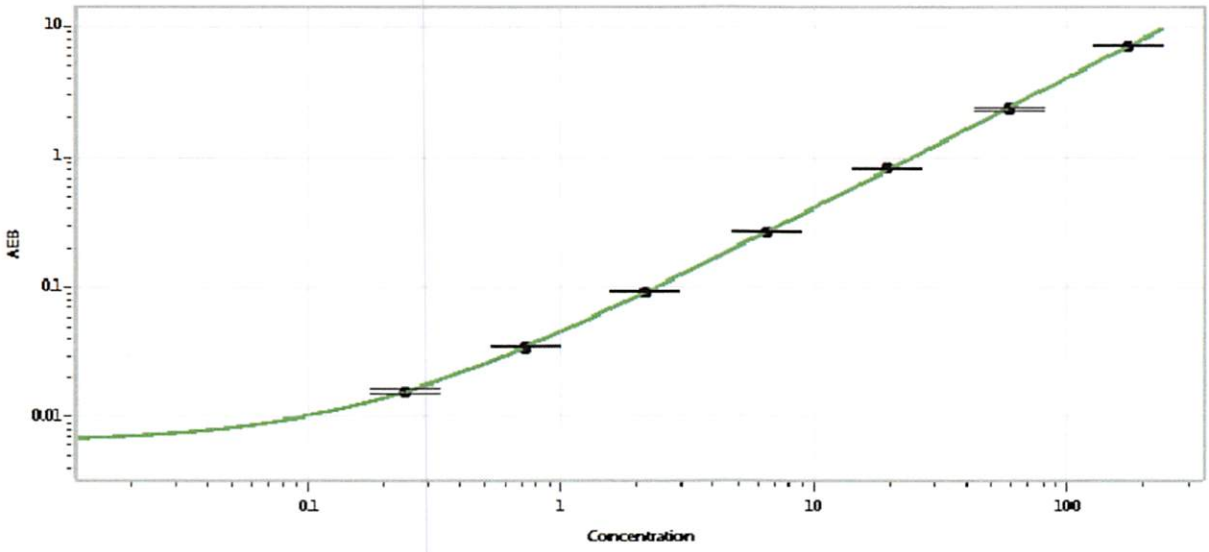
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102101	904407
Detector Stock	102102	904207
Calibrator Concentrate	102105	MF-010518-05
Calibrator Diluent	102103	901016
Sample Diluent	102104	902923

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	3500 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.240
C	0.720
D	2.16
E	6.48
F	19.4
G	58.3
H	175

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P.S 27 Feb 2019
Signature/Date