

Product Number:	102100
Lot Number:	501633
Expiration:	8-Jun-2019
Platform(s):	HD-1



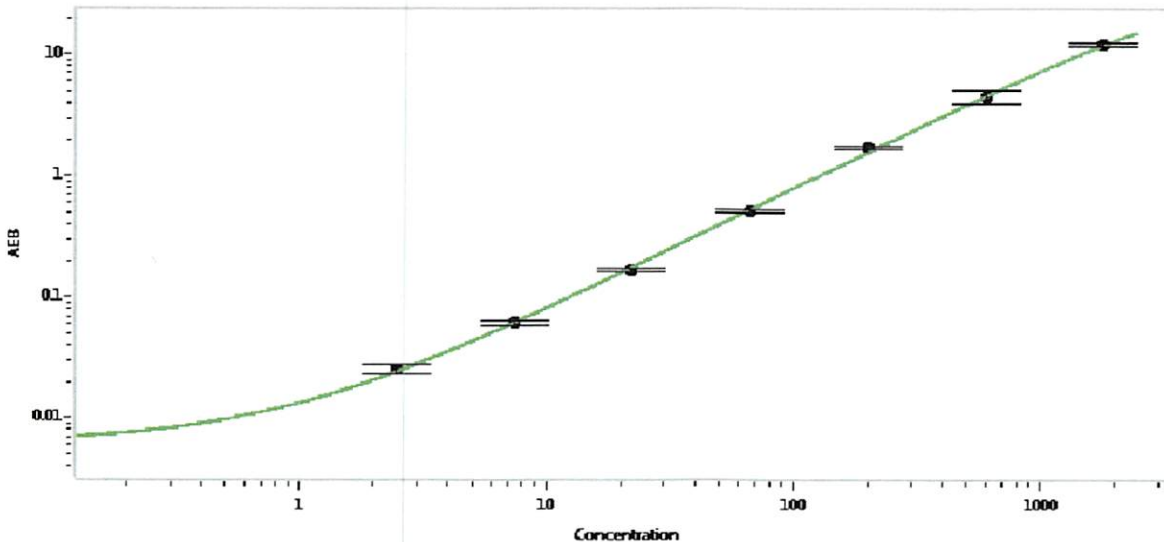
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102092	902804
Detector Stock	102093	903302
Calibrator Concentrate	102096	MF-060817-03
Calibrator Diluent	102094	901913
Sample Diluent	102095	901912

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.47
C	7.41
D	22.2
E	66.6
F	200
G	600
H	1800

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 13 Feb 2019
	Name	Title	Signature/Date