

Product Number:	101968
Lot Number:	501451
Expiration:	20-Feb-2019
Platform(s):	HD-1



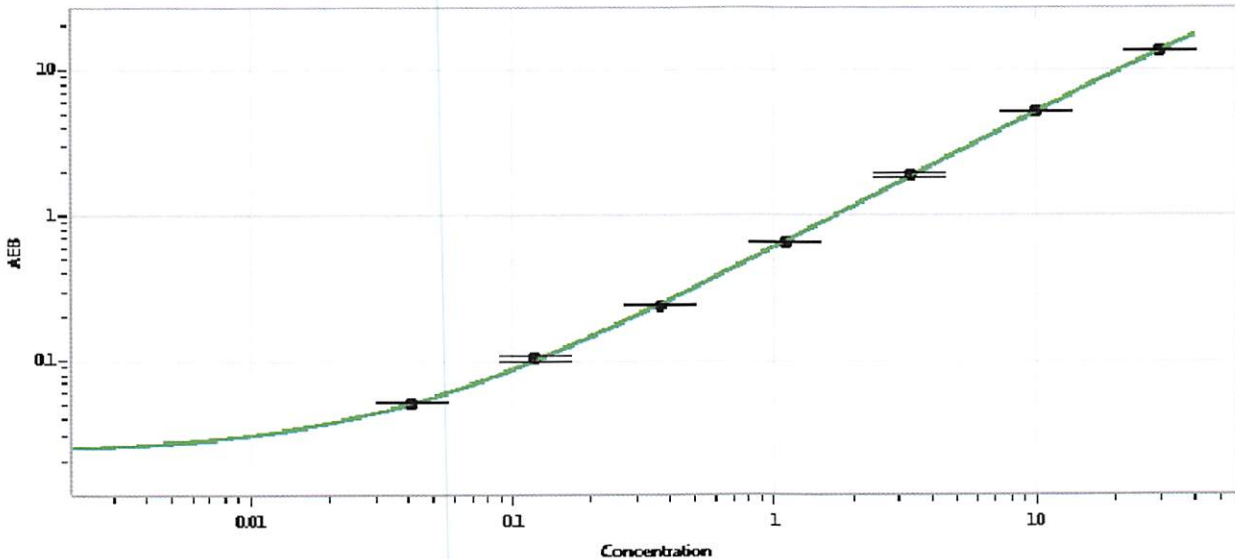
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101966	828406
Detector Stock	101967	828306
Calibrator Concentrate	101965	MF-110817-03
Calibrator Diluent	101963	823618
Sample Diluent	101964	827409

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	600 pg/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 18 OCT 2018
(Signature/Date)