

Product Number:	102100
Lot Number:	501405
Expiration:	7-Nov-2018
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102092	MF-052118-01
Detector Stock	102093	MF-051618-09
Calibrator Concentrate	102096	MF-060817-03
Calibrator Diluent	102094	MF-050718-03
Sample Diluent	102095	MF-050718-02

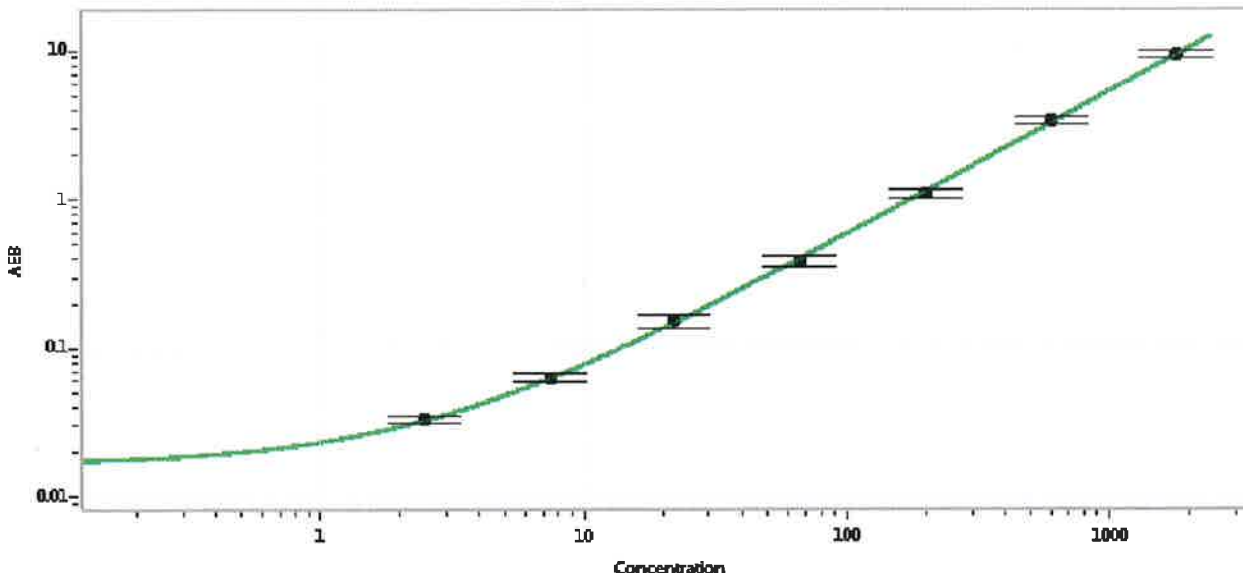


Data below represents results generated on the

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

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
11/SEP/2018
Signature/Date