

Product Number:	102517
Lot Number:	501412
Expiration:	3-Jan-2019
Platform(s):	HD-1 & SR-X



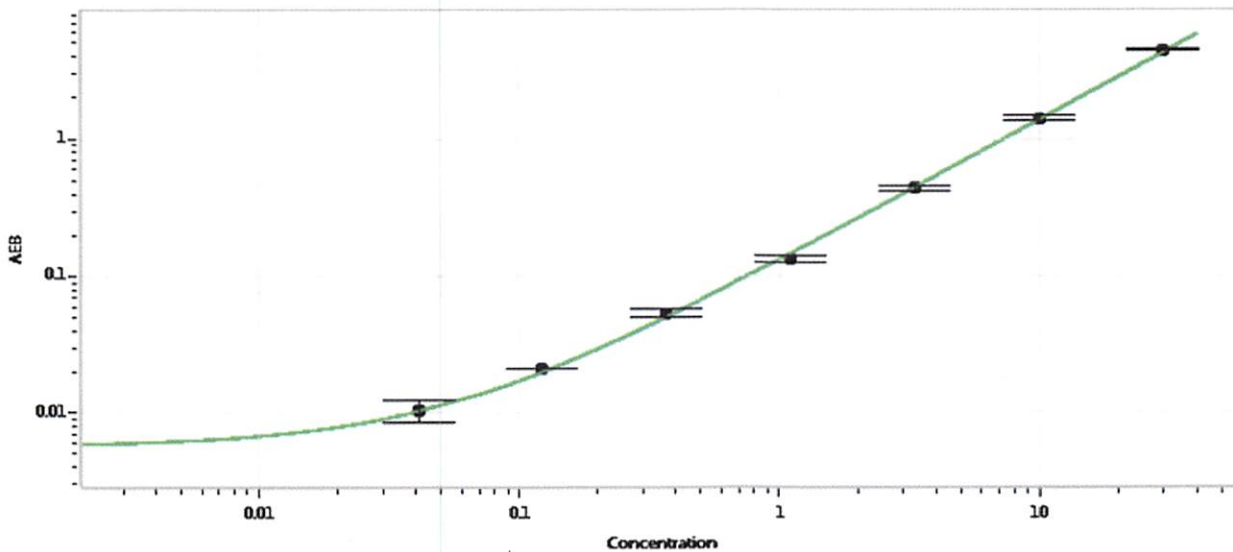
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102510	822607
Detector Stock	102511	822601
Calibrator Concentrate	102514	MF-022818-09
Calibrator Diluent	102512	818418
Sample Diluent	102513	818417

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]*  
Signature/Date  
17 SEP 2018