

Product Number:	102343
Lot Number:	501372
Expiration:	20-Oct-2018
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102337	822504
Detector Stock	102338	819013
Calibrator Concentrate	102342	MF-102016-02
Calibrator Diluent	102339	819009
Sample Diluent	102340	819010

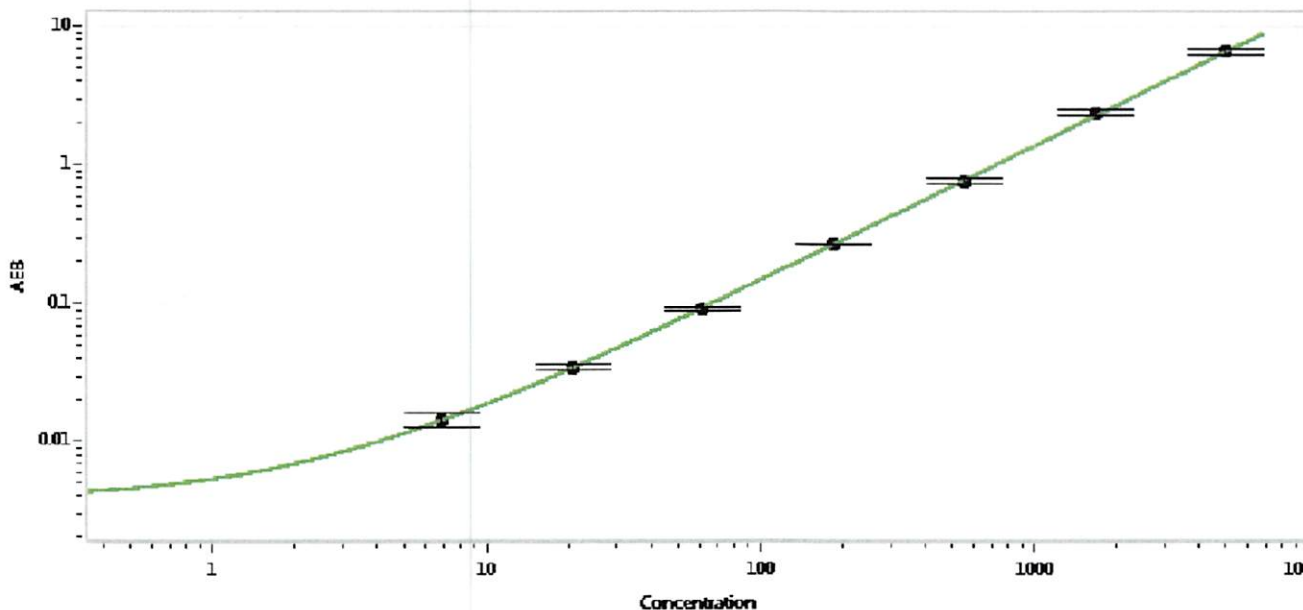


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Lot Number
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Concentrate	101957
RGP Reagent	103159

Calibrator Concentrate Target Value	100 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	6.86
C	20.6
D	61.7
E	185
F	556
G	1667
H	5000

Review/Approval

Sivasankari Murugan QC Supervisor  
Name Title

*[Signature]* 9/18/2018  
Signature/Date