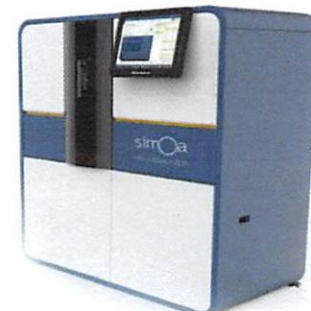


Product Number:	102669
Lot Number:	501336
Expiration:	4-Oct-2018
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102662	MF-062618-04
Detector Stock	102663	819908
Calibrator Concentrate	102666	MF-020718-03
Calibrator Diluent	102664	MF-040418-03
Sample Diluent	102665	MF-062618-03

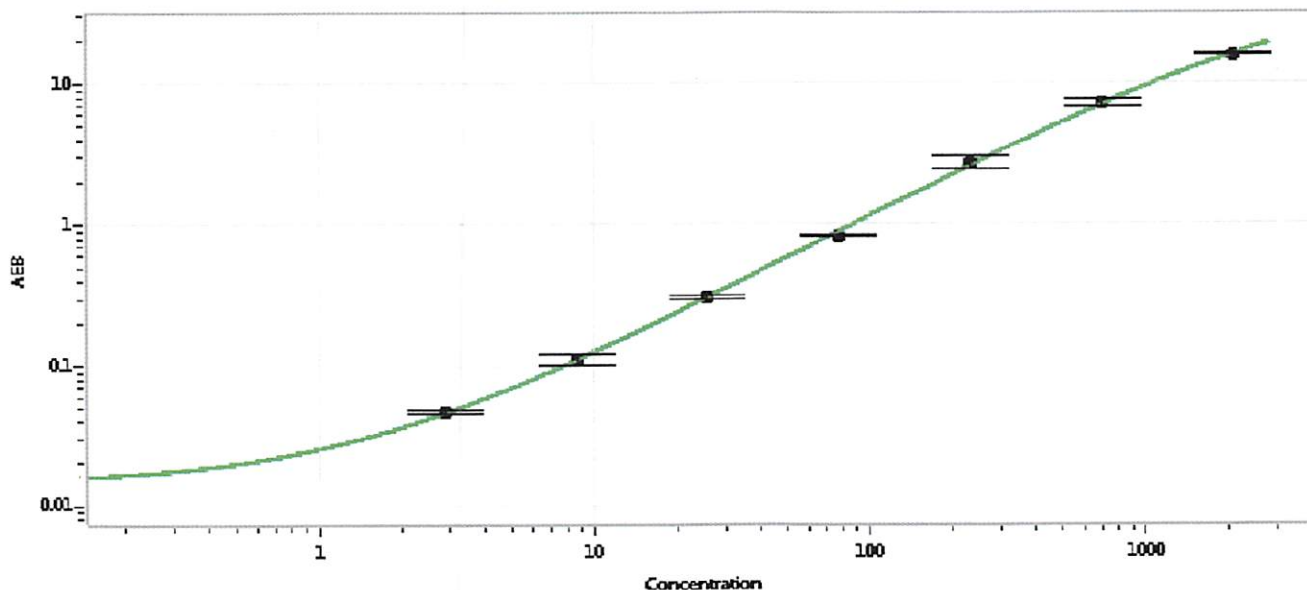


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Concentrate	101957
RGP Reagent	103159

Calibrator Concentrate Target Value	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.88
C	8.64
D	25.9
E	77.8
F	233
G	700
H	2100

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 22 Jul 2018
Signature/Date