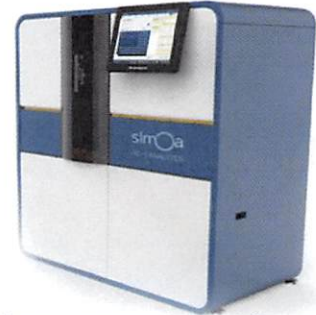


Product Number:	102343
Lot Number:	501211
Expiration:	6-Sep-2018
Platform(s):	HD-1



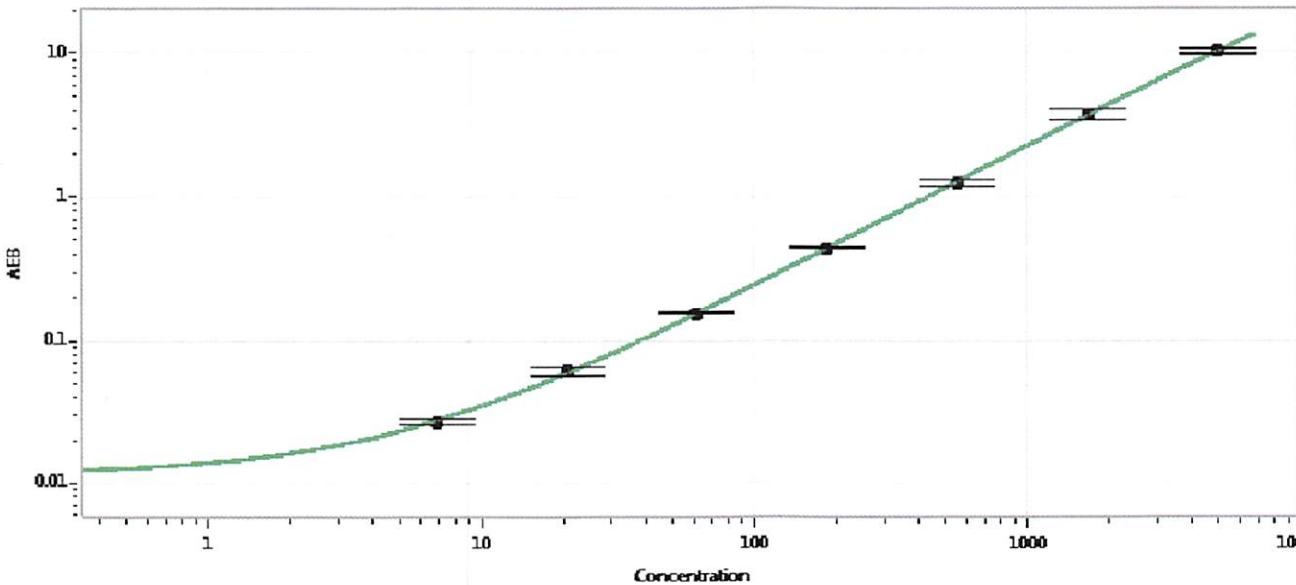
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102337	MF-040618-01
Detector Stock	102338	MF-041018-01
Calibrator Concentrate	102342	MF-102016-02
Calibrator Diluent	102339	MF-030618-09
Sample Diluent	102340	MF-040518-03

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	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
Name

QC Supervisor
Title

[Signature] 17 APR 2018
Signature/Date