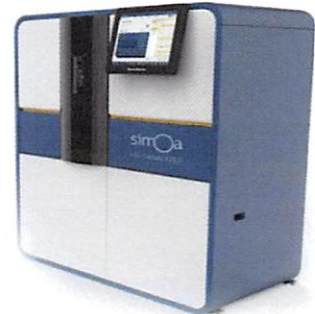


Product Number:	101832
Lot Number:	501196
Expiration:	28-Sep-2018
Platform(s):	HD-1



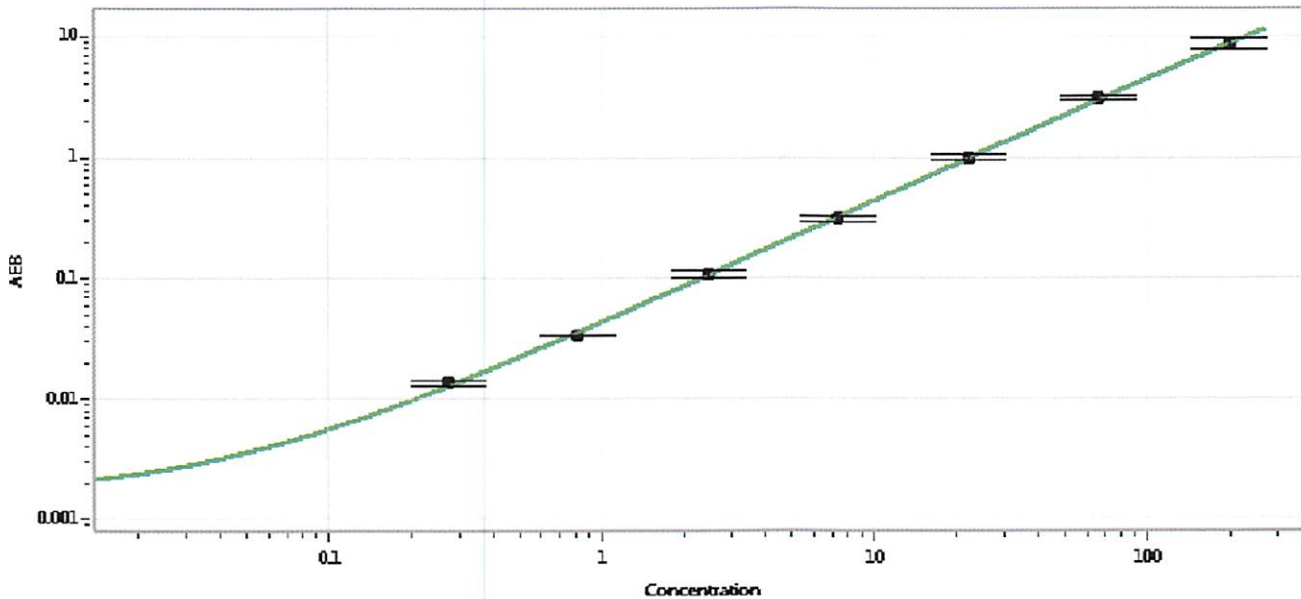
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101927	MF-040218-08
Detector Stock	101928	MF-032818-07
Calibrator Concentrate	101827	MF-041618-09
Calibrator Diluent	101945	MF-032018-05
Sample Diluent	101946	MF-120117-06

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

Calibrator Concentrate Target Value	4000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sivasankari Murugan QC Supervisor  
Name Title

*[Signature]* 20 APR 2018  
Signature/Date