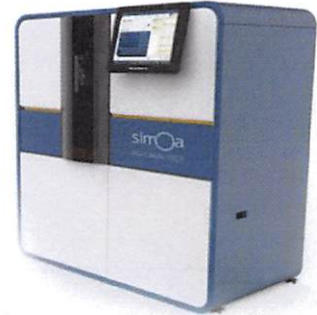


Product Number:	102475
Lot Number:	501345
Expiration:	19-Dec-2018
Platform(s):	HD-1



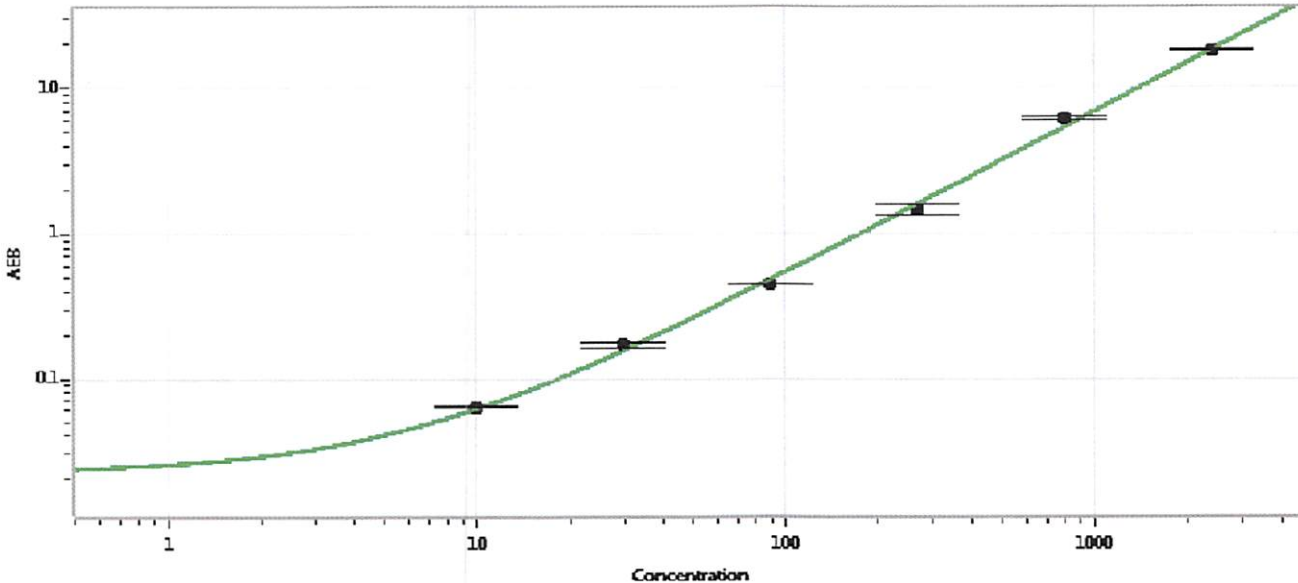
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102468	MF-061918-08
Detector Stock	102469	820602
Calibrator Concentrate	102472	MF-062018-06
Calibrator Diluent	102470	MF-050318-04
Sample Diluent	102471	MF-041018-21

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	9.88
C	29.6
D	88.9
E	267
F	800
G	2400

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 26/12/2018  
Signature/Date