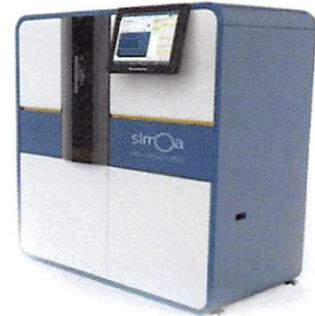


Product Number:	102570
Lot Number:	501241
Expiration:	16-Oct-2018
Platform(s):	HD-1



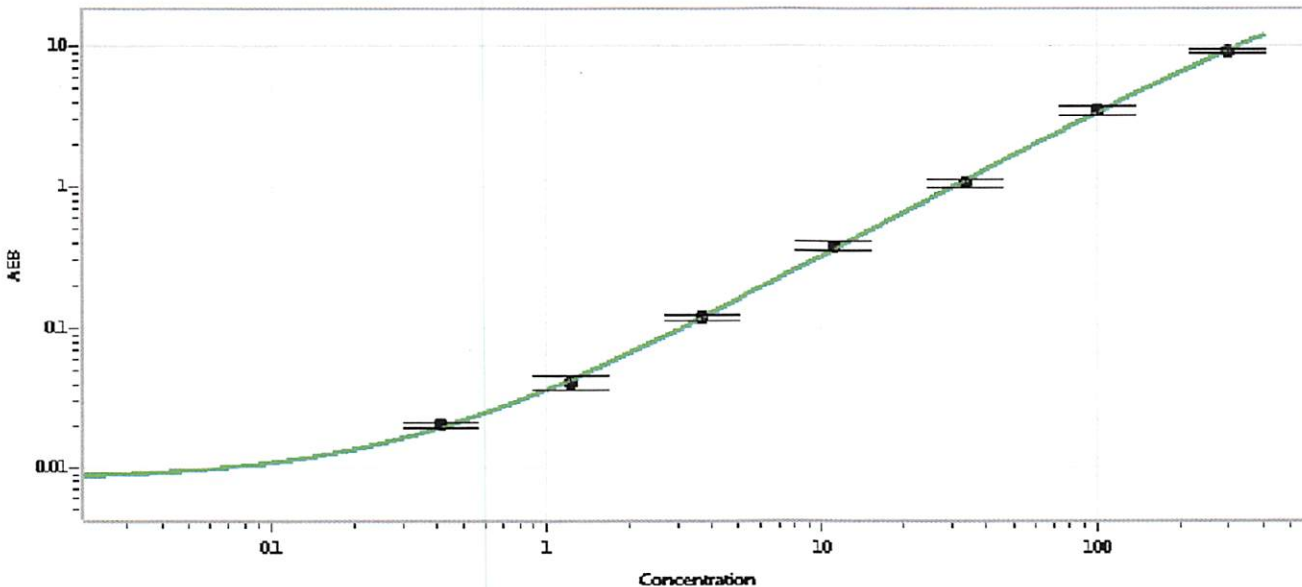
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102563	MF-050318-03
Detector Stock	102564	MF-050118-05
Detector Diluent	102590	MF-041618-03
Calibrator Concentrate	102567	MF-022217-07
Calibrator Diluent	102565	MF-041018-14
Sample Diluent	102566	MF-041018-13

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

Calibrator Concentrate Target Value	6 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
8/20/18