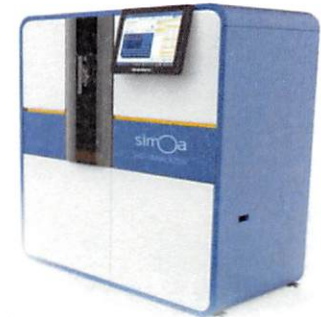


Product Number:	101863
Lot Number:	501192
Expiration:	15-Sep-2018
Platform(s):	HD-1



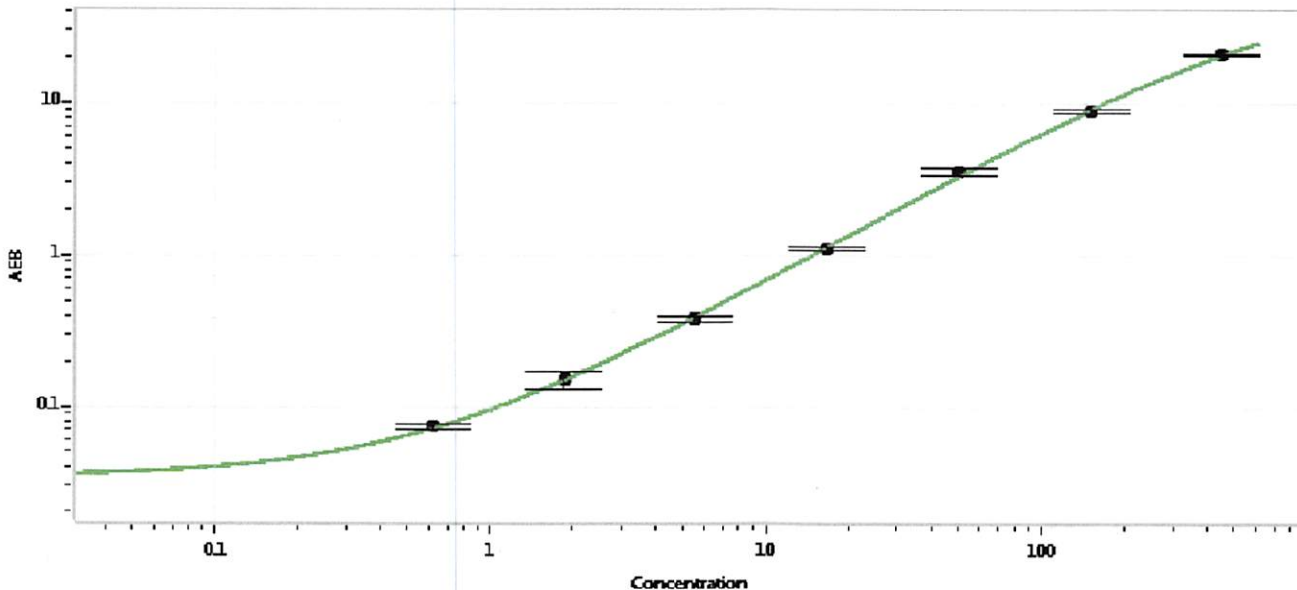
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101904	MF-032718-06
Detector Stock	101905	MF-032818-03
Calibrator Concentrate	101858	MF-031518-02
Calibrator Diluent	101990	MF-111317-02
Sample Diluent	101991	MF-031518-03

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	9 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450

Review/Approval	Sivasankari Murugan	QC Supervisor	<i>[Signature]</i> 11 APR 2018
	Name	Title	Signature/Date