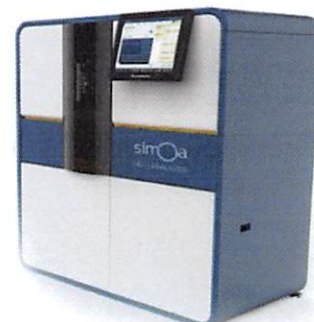


Product Number:	102394
Lot Number:	501489
Expiration:	20-Feb-2019
Platform(s):	HD-1



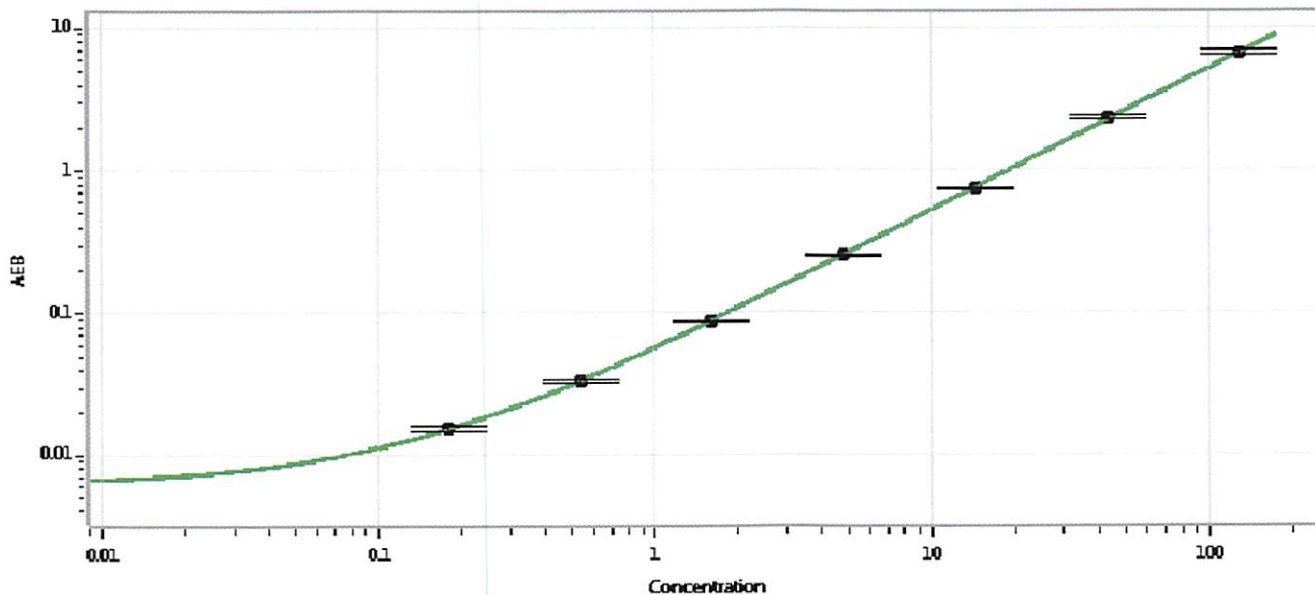
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102387	831107
Detector Stock	102388	831215
Calibrator Concentrate	102391	MF-010518-06
Calibrator Diluent	102389	829710
Sample Diluent	102390	826216

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

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

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.178
C	0.535
D	1.60
E	4.81
F	14.4
G	43.3
H	130

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
29 NOV 2018