

Product Number:	101863
Lot Number:	501510
Expiration:	20-Feb-2019
Platform(s):	HD-1



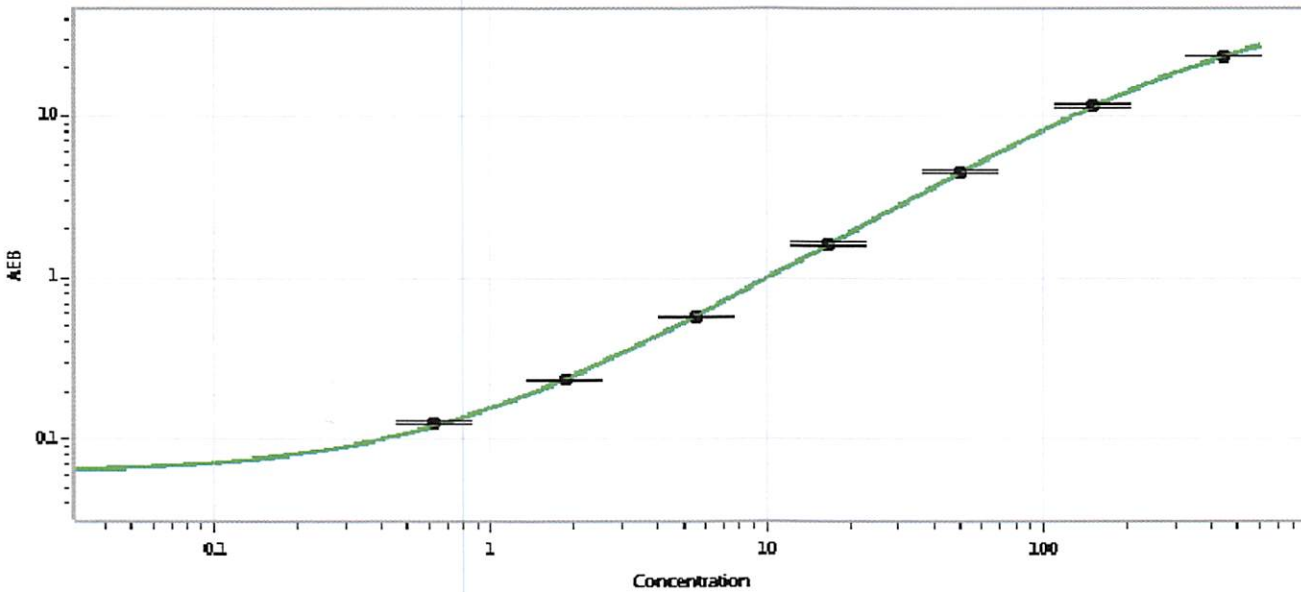
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	101904	832411
Detector Stock	101905	830508
Calibrator Concentrate	101858	MF-031518-02
Calibrator Diluent	101990	827506
Sample Diluent	101991	830408

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	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
Name

QC Supervisor
Title

[Signature] 01DEC2018
Signature/Date