

Product Number:	101808
Lot Number:	501490
Expiration:	20-Feb-2019
Platform(s):	HD-1



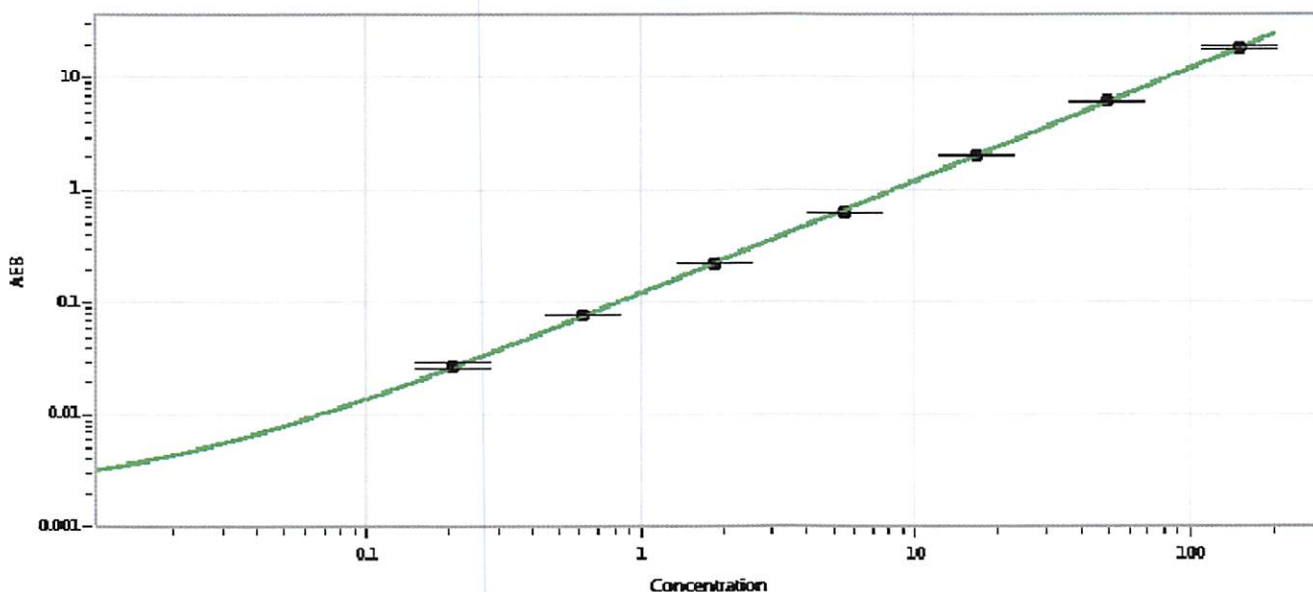
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	101922	831606
Detector Stock	101923	827603
Calibrator Concentrate	101803	831707
Calibrator Diluent	101925	827502
Sample Diluent	101926	827604

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	3 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0
H	150

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 29 Nov 2018
Signature/Date