

Product Number:	101871
Lot Number:	501465
Expiration:	20-Feb-2019
Platform(s):	HD-1



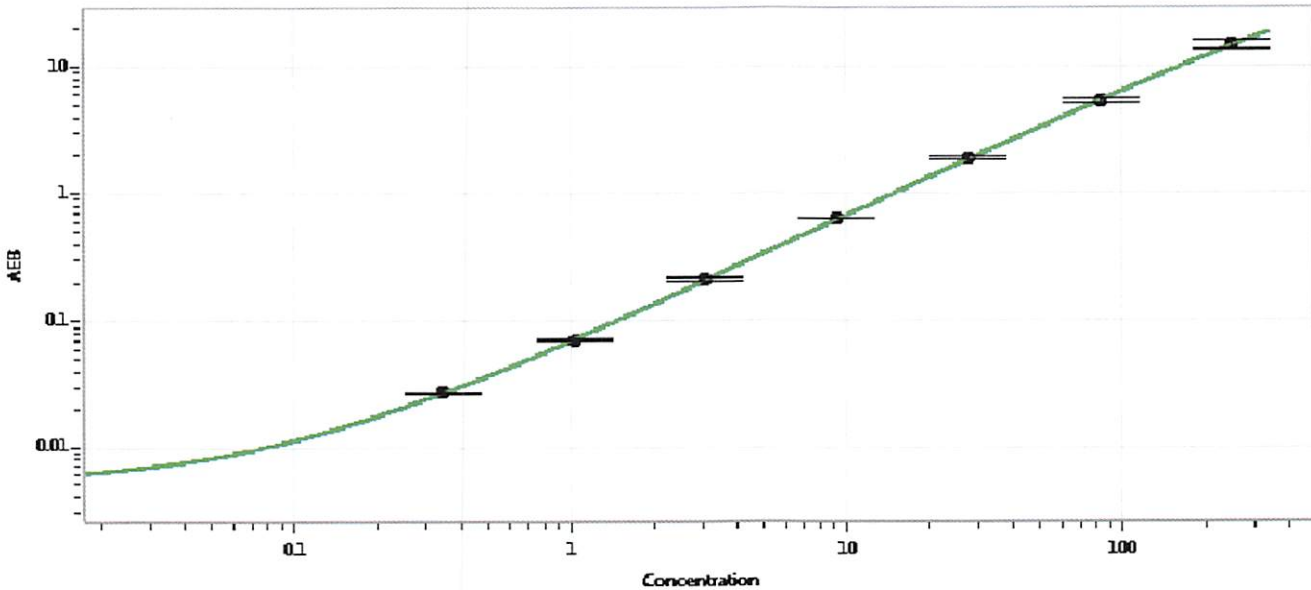
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101893	828905
Detector Stock	101894	828603
Calibrator Concentrate	101866	829602
Calibrator Diluent	101959	827611
Sample Diluent	101958	827410

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	5 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.343
C	1.03
D	3.09
E	9.26
F	27.8
G	83.3
H	250

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 25 OCT 2018
Signature/Date