

Product Number:	102570
Lot Number:	501463
Expiration:	20-Feb-2019
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102563	828906
Detector Stock	102564	828904
Detector Diluent	102590	828507
Calibrator Concentrate	102567	829005
Calibrator Diluent	102565	826214
Sample Diluent	102566	826213

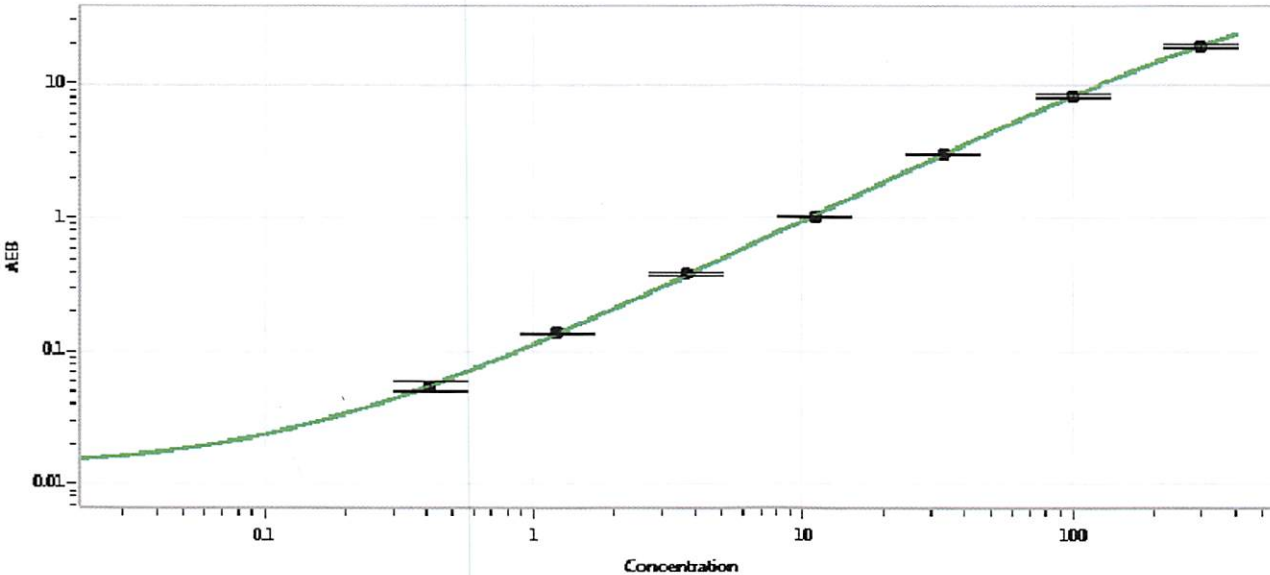


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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

Calibrator Concentrate Target Value	6 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 22 OCT 2018
Signature/Date