

Product Number:	102635
Lot Number:	501366
Expiration:	26-Jan-2019
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102628	821106
Detector Stock	102629	820706
Calibrator Concentrate	102632	822101
Calibrator Diluent	102630	819812
Sample Diluent	102631	820803

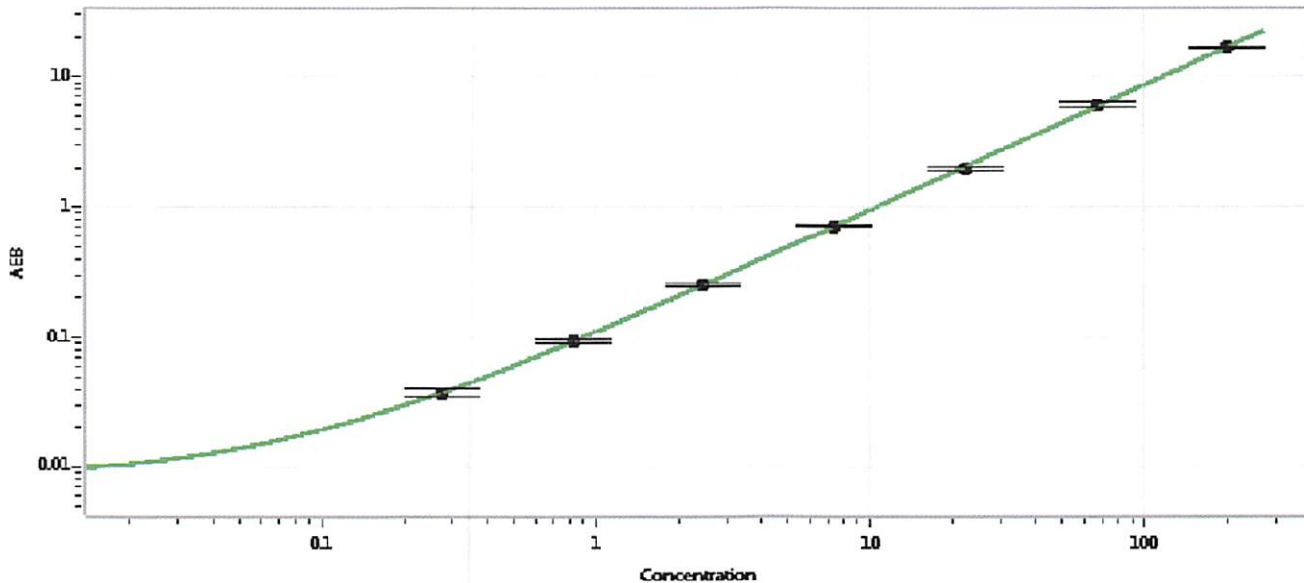


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Concentrate	101957
RGP Reagent	103159

Calibrator Concentrate Target Value	4 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
13 Aug 2018