

Product Number:	102599
Lot Number:	501526
Expiration:	15-Jan-2019
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102592	803204
Detector Stock	102593	807106
Calibrator Concentrate	102596	MF-011618-01
Calibrator Diluent	102594	801503
Sample Diluent	102595	801504

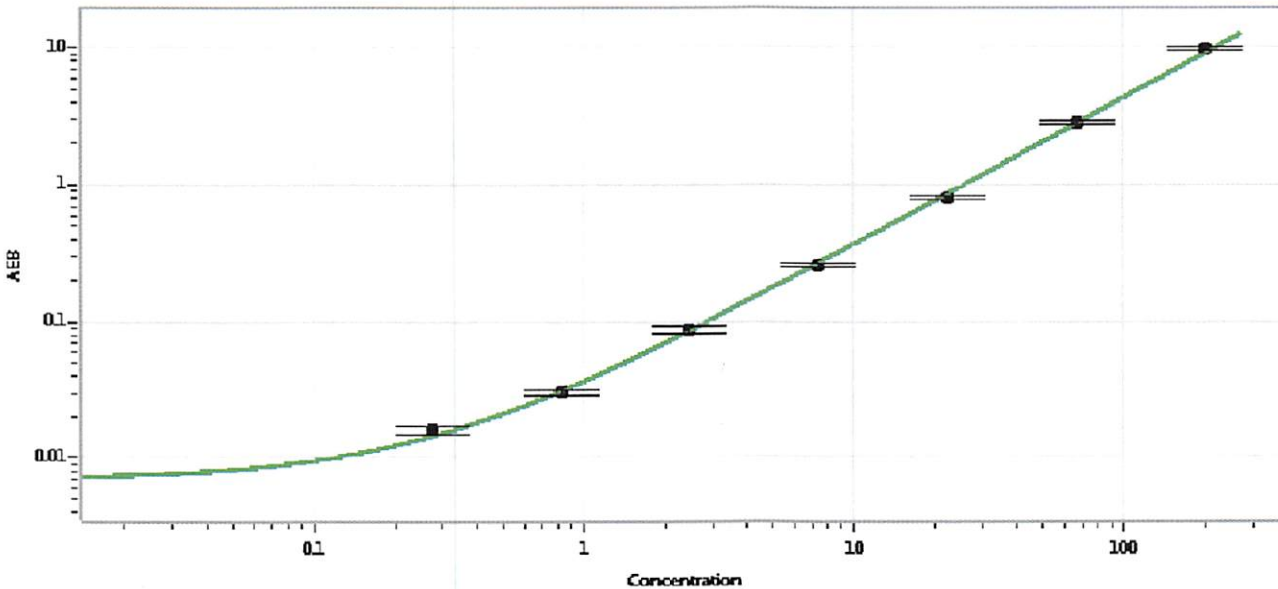


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
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

*S. Murugan* 04DEC2018  
Signature/Date