

Product Number:	102039
Lot Number:	501394
Expiration:	7-Feb-2019
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Stock	102028	821904
Detector Stock	102029	821902
Calibrator Concentrate	102032	MF-042618-03
Calibrator Diluent	102030	819817
Sample Diluent	102031	821903

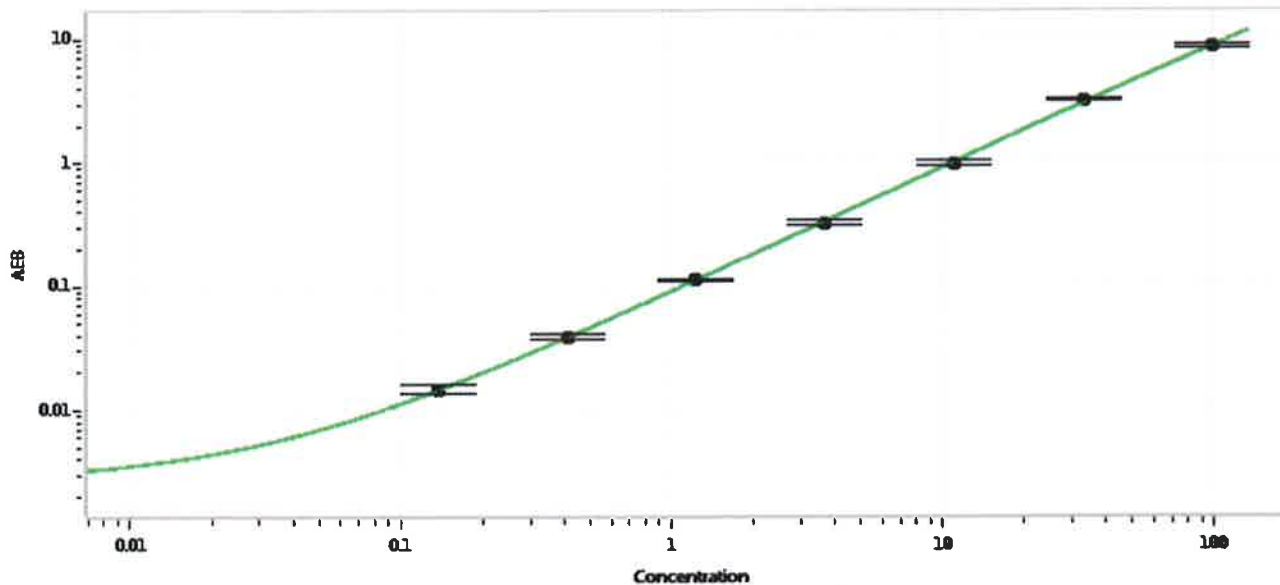


Data below represents results generated on the Simoa™ HD-1

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	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
10 SEP 2018