

Product Number:	102039
Lot Number:	501530
Expiration:	14-May-2019
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Stock	102028	832309
Detector Stock	102029	833106
Calibrator Concentrate	102032	MF-042618-03
Calibrator Diluent	102030	829712
Sample Diluent	102031	831813

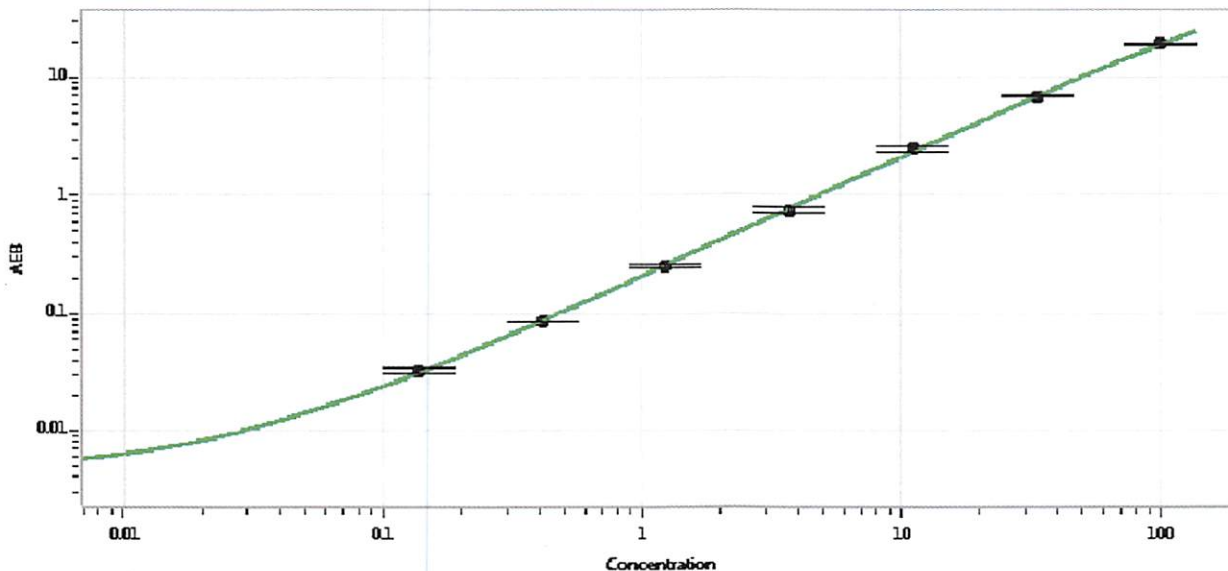


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	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] 06 DEC 2018
Signature/Date