

Product Number:	101419
Lot Number:	501075
Expiration:	28-Apr-18
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	101899	MF-011518-04
Detector Stock	101900	MF-011518-03
Calibrator Concentrate	101420	MF-042816-02
Calibrator Diluent	101418	MF-112917-15
Sample Diluent	101416	MF-122117-07

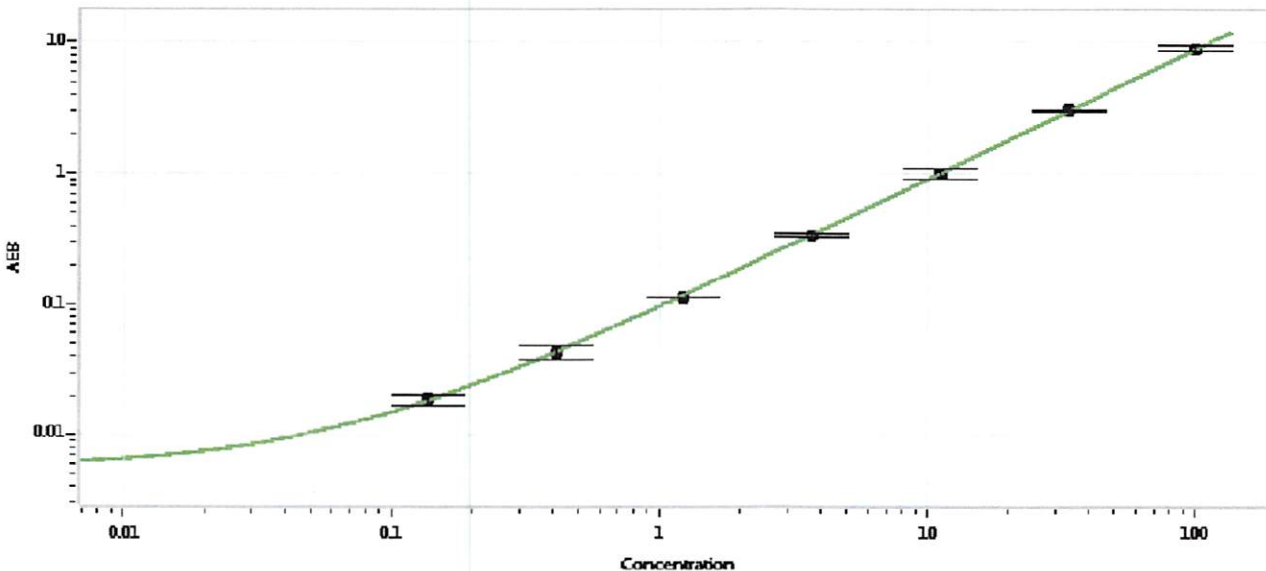


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	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] 22 Jan 2018
Signature/Date