

Product Number:	102343
Lot Number:	501408
Expiration:	20-Oct-2018
Platform(s):	HD-1



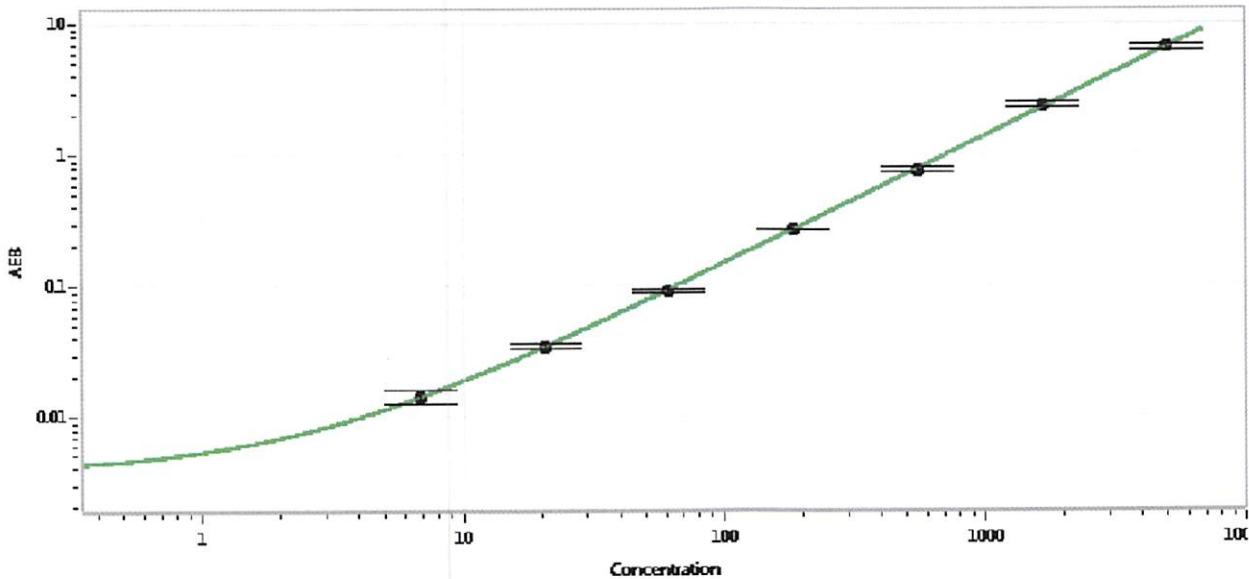
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102337	822504
Detector Stock	102338	819013
Calibrator Concentrate	102342	MF-102016-02
Calibrator Diluent	102339	819009
Sample Diluent	102340	819010

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	100 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	6.86
C	20.6
D	61.7
E	185
F	556
G	1667
H	5000

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

S. Murugan 12SEP2018
Signature/Date