

Product Number:	102491
Lot Number:	501402
Expiration:	28-Feb-2019
Platform(s):	HD-1



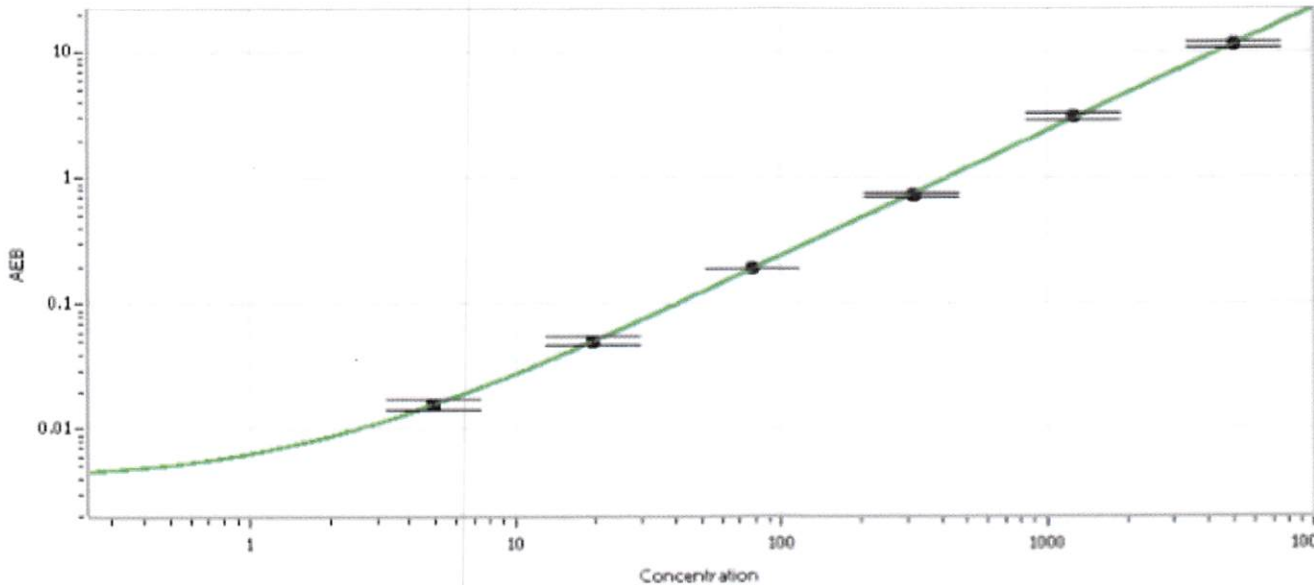
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102484	825306
Detector Stock	102485	824902
Detector Diluent	103084	824314
Calibrator Concentrate	102488	MF-100217-02
Calibrator Diluent	102486	824304
Sample Diluent	102487	824205

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	100 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.88
C	19.5
D	78.1
E	313
F	1250
G	5000

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 12 SEP 2018  
Signature/Date