

Product Number:	102343
Lot Number:	501500
Expiration:	20-Feb-2019
Platform(s):	HD-1



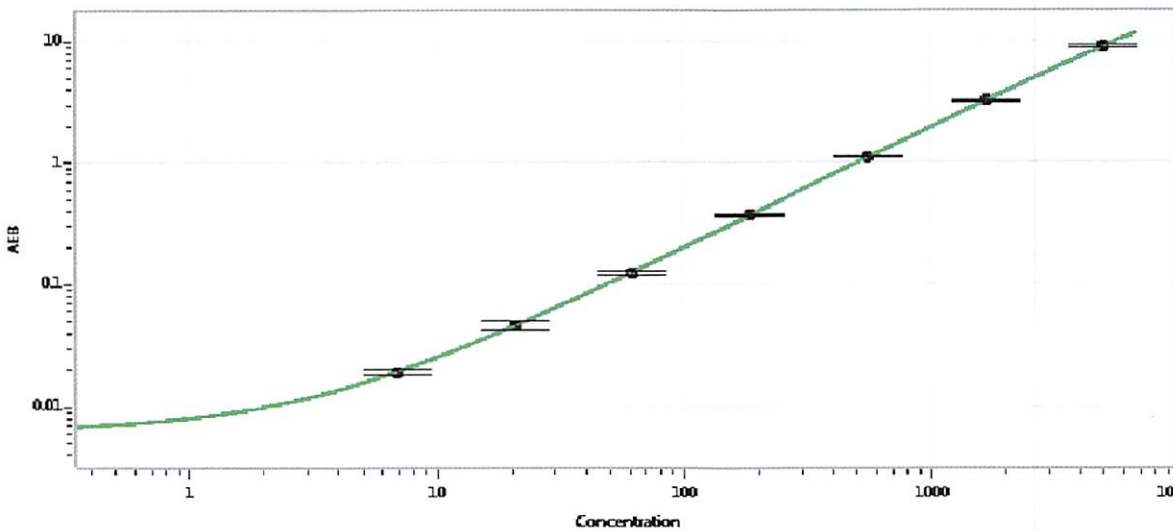
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102337	831904
Detector Stock	102338	831206
Calibrator Concentrate	102342	831815
Calibrator Diluent	102339	827412
Sample Diluent	102340	830901

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	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

[Signature] 16/NOV/2018
Signature/Date