

Product Number:	101855
Lot Number:	501440
Expiration:	27-Mar-19
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	101902	827103
Detector Stock	101903	827011
Calibrator Concentrate	101850	MF-121517-07
Calibrator Diluent	101988	823616
Sample Diluent	101989	827408

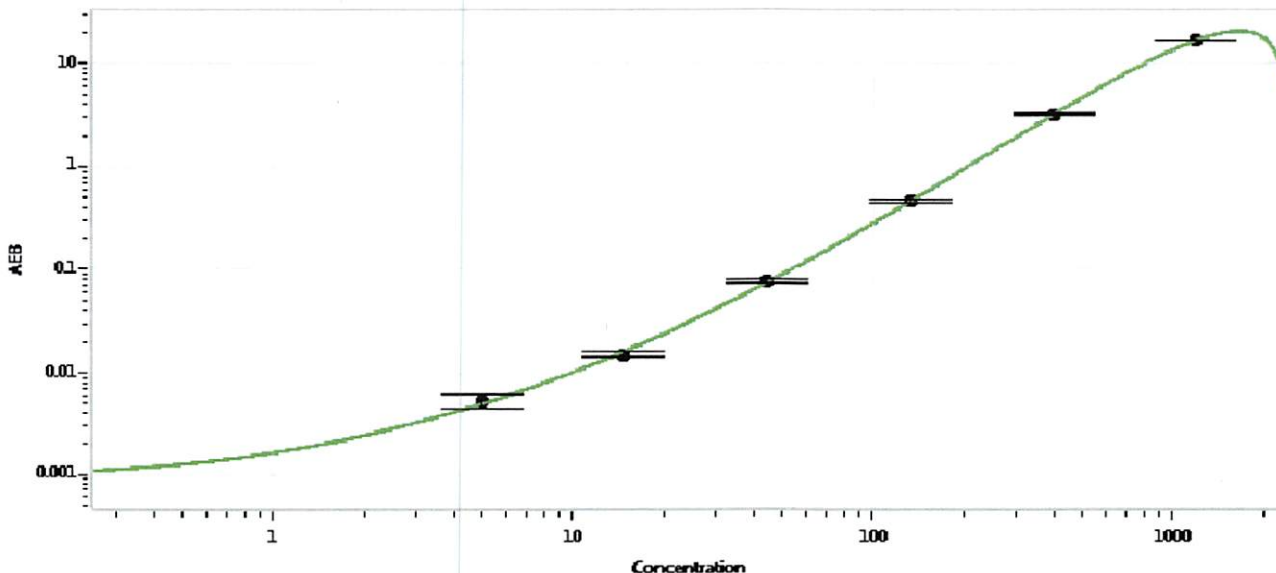


Data below represents results generated on the Simoa™ HD-1

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	30 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.94
C	14.8
D	44.4
E	133
F	400
G	1200

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 15 OCT 2018
Signature/Date