

Product Number:	102336
Lot Number:	501418
Expiration:	20-Feb-2019
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102330	826103
Detector Stock	102331	824901
Calibrator Concentrate	102335	826303
Calibrator Diluent	102332	819018
Sample Diluent	102333	824805

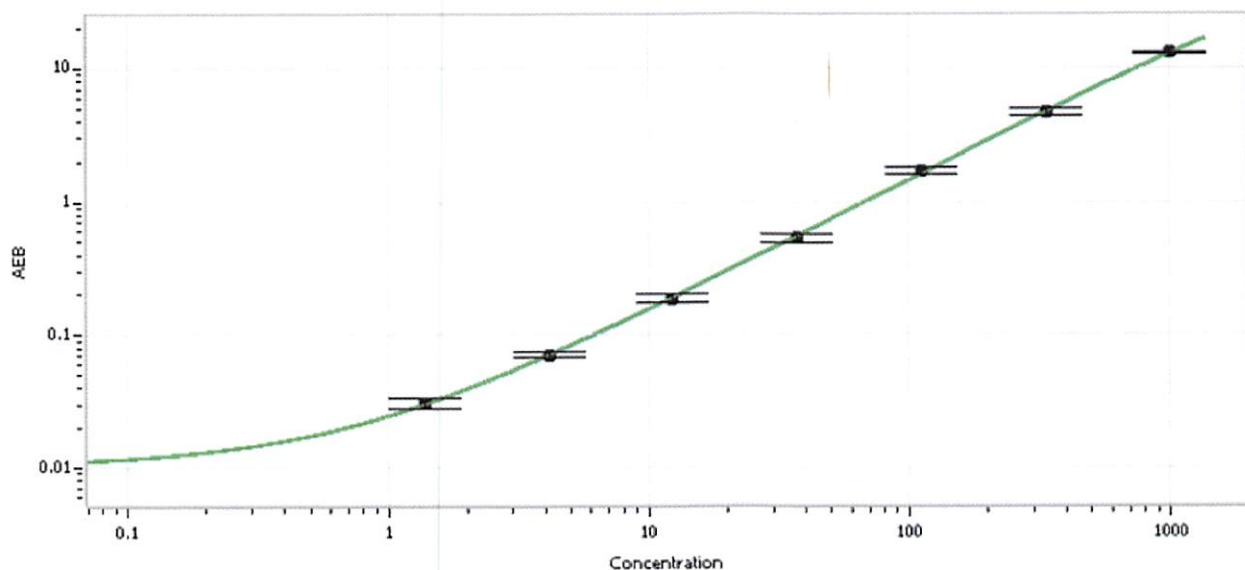


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159

Calibrator Concentrate Target Value	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date 05/NOV/2018