

Product Number:	102336
Lot Number:	501418
Expiration:	5-Mar-2019
Platform(s):	HD-1



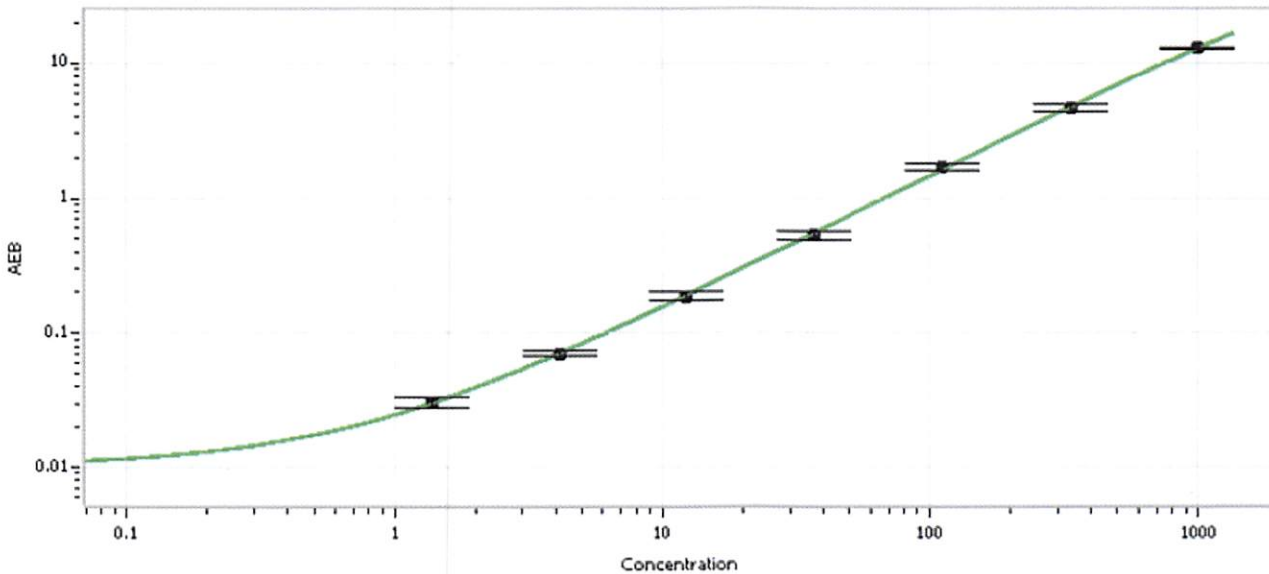
Data below represents results generated on the Simoa™ HD-1 Analyzer.

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

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	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