

Product Number:	102064
Lot Number:	501503
Expiration:	27-Mar-2019
Platform(s):	HD-1



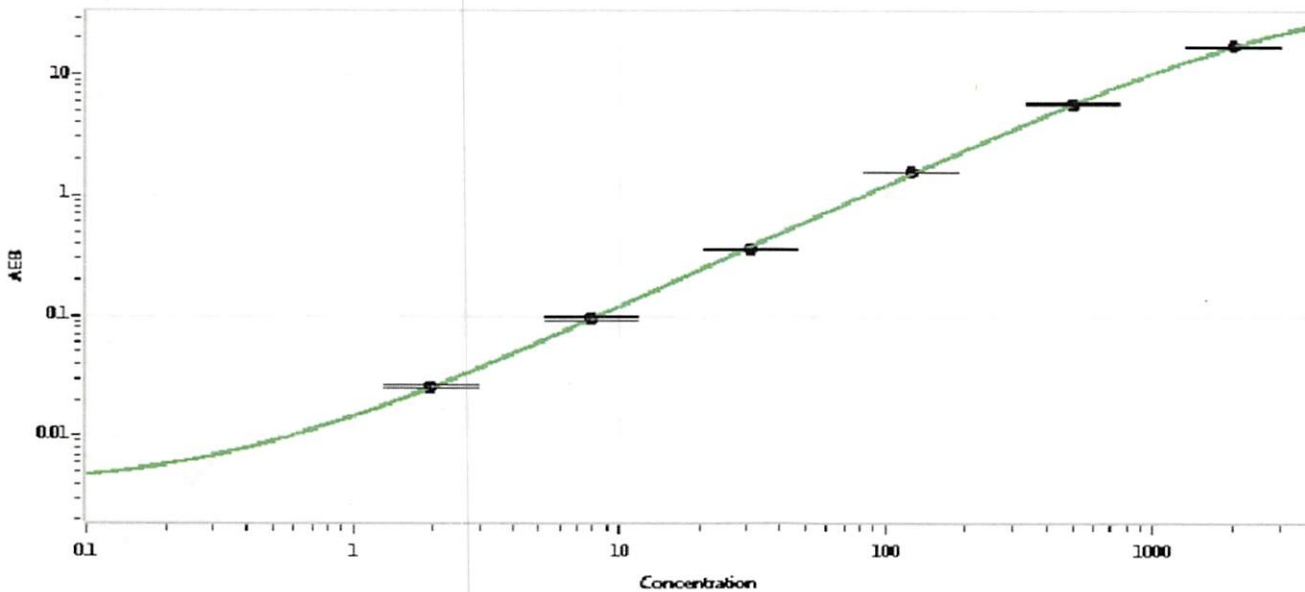
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102056	831811
Detector Stock	102057	827403
Calibrator Concentrate	102060	833709
Calibrator Diluent	102058	827006
Sample Diluent	102059	827007

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	40 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.95
C	7.81
D	31.3
E	125
F	500
G	2000

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 26DEC2018
Signature/Date