

Product Number:	102504
Lot Number:	501546
Expiration:	28-Feb-2019
Platform(s):	HD-1



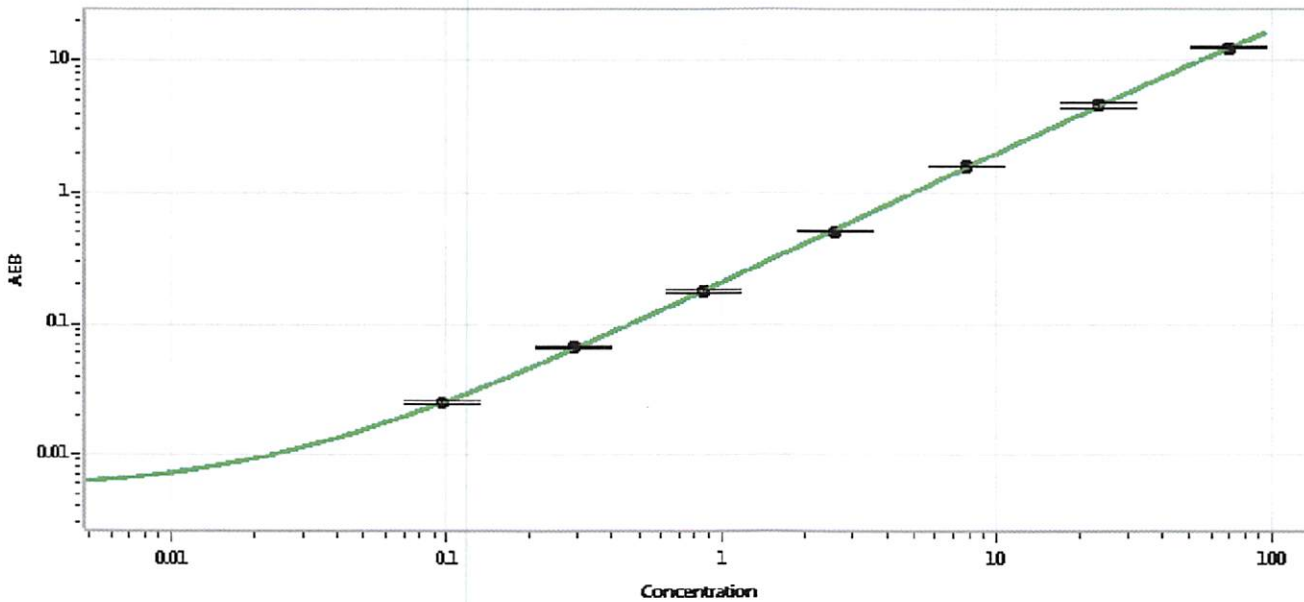
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102497	809206
Detector Stock	102498	809207
Calibrator Concentrate	102501	MF-022817-06
Calibrator Diluent	102499	808702
Sample Diluent	102500	809304

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	1.4 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.096
C	0.288
D	0.864
E	2.59
F	7.77
G	23.3
H	70.0

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 12 DEC 2018
Signature/Date