

Product Number:	103093
Lot Number:	501521
Expiration:	1-Feb-2019
Platform(s):	HD-1 & SR-X



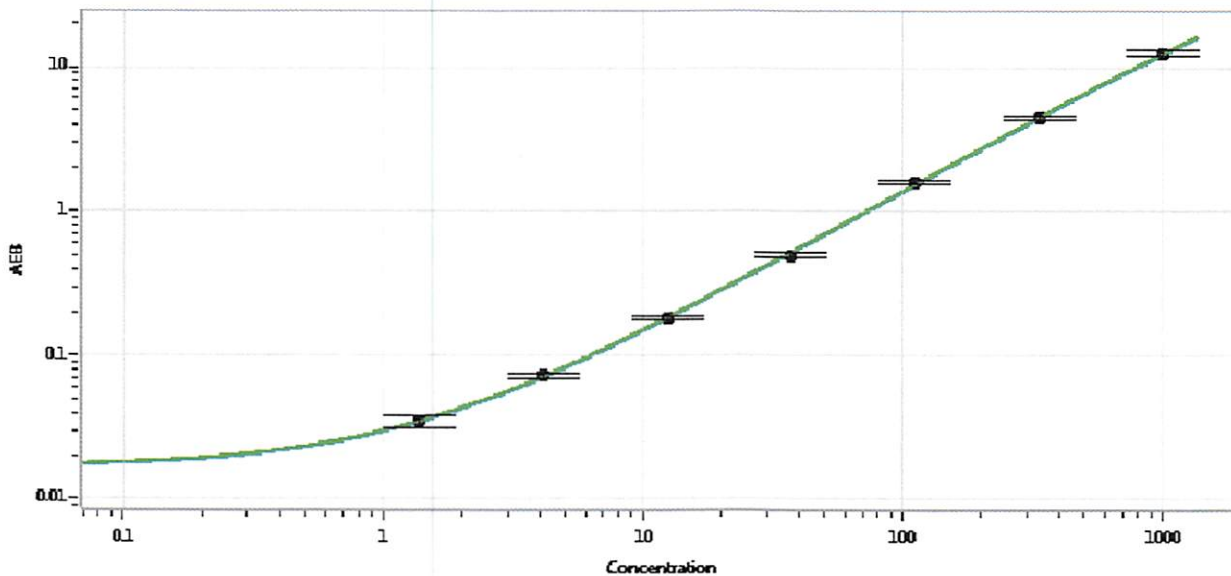
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102021	832011
Detector Stock	103043	821310
Calibrator Concentrate	103047	MF-010818-10
Calibrator Diluent	103044	832008
Sample Diluent	103046	831805

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	12 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 05DEC2018
Signature/Date