

Product Number:	102794
Lot Number:	501133
Expiration:	10-Aug-2018
Platform(s):	HD-1



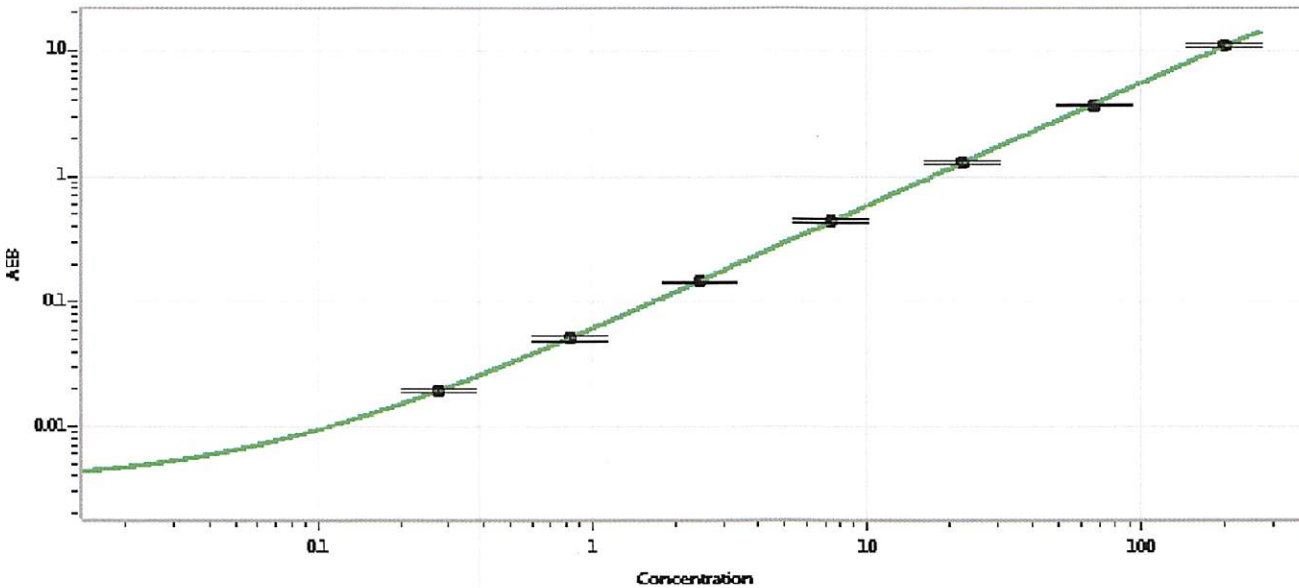
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102787	MF-021418-01
Detector Stock	102788	MF-021018-07
Calibrator Concentrate	102791	MF-062817-04
Calibrator Diluent	102789	MF-021018-05
Sample Diluent	102790	MF-021018-06

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	4 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 23 FEB 2018
Signature/Date