

Product Number:	102599
Lot Number:	501122
Expiration:	15-Jul-2018
Platform(s):	HD-1



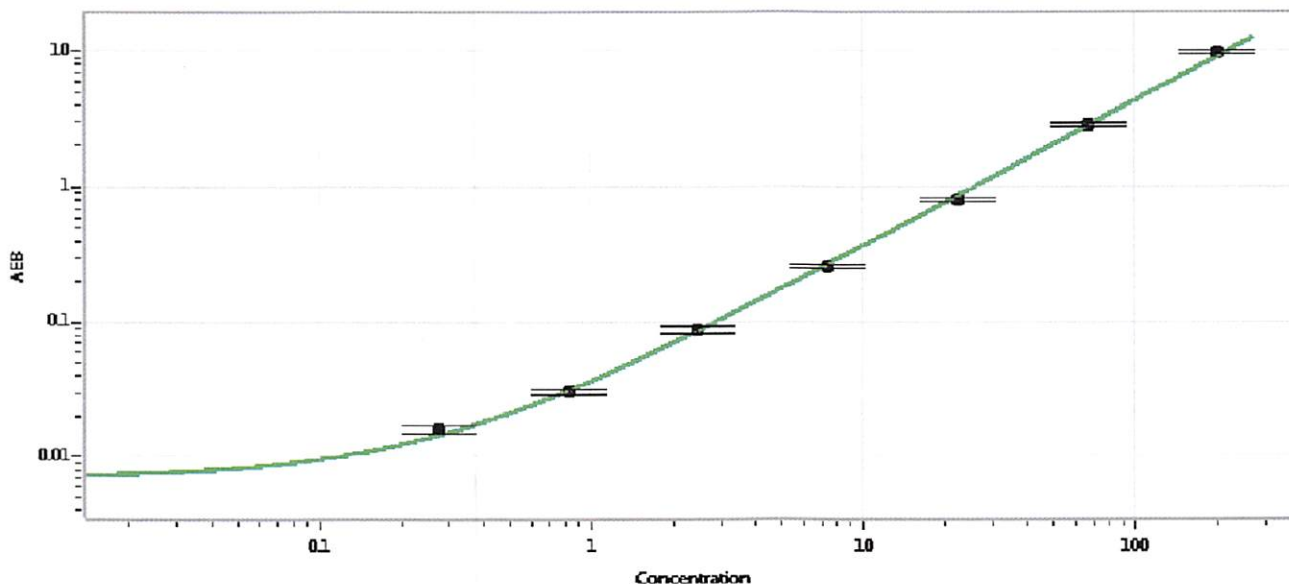
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102592	MF-020118-01
Detector Stock	102593	MF-031218-06
Calibrator Concentrate	102596	MF-011618-01
Calibrator Diluent	102594	MF-011518-10
Sample Diluent	102595	MF-011518-12

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]* 19 MAR 2018  
Signature/Date