

Product Number:	102100
Lot Number:	501291
Expiration:	7-Nov-2018
Platform(s):	HD-1



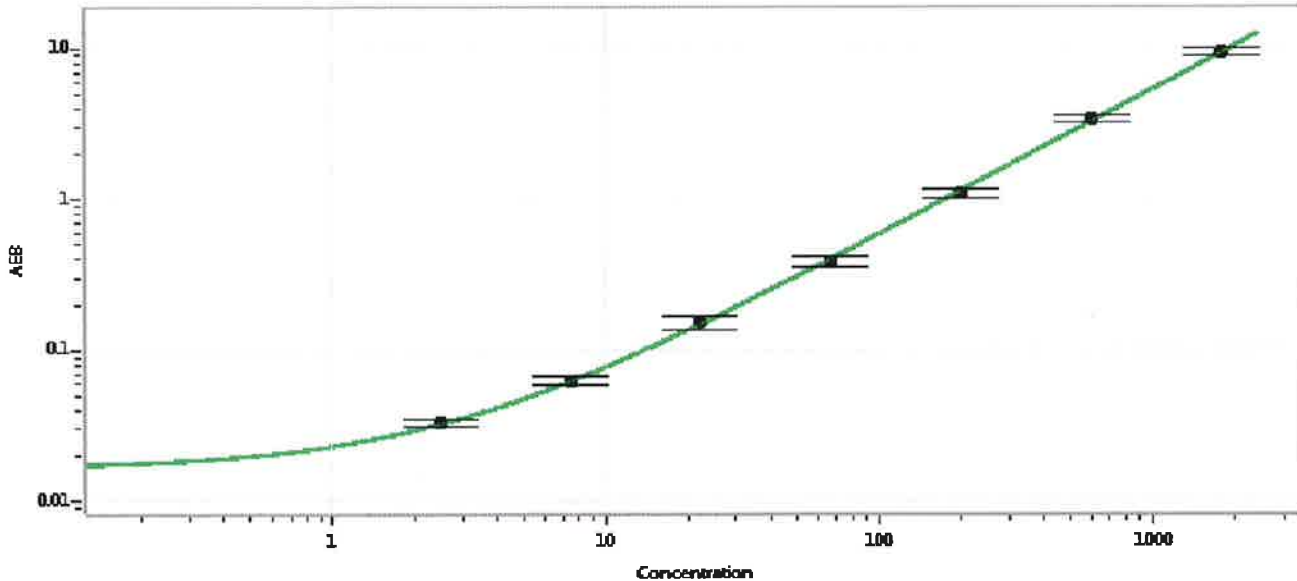
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102092	MF-052118-01
Detector Stock	102093	MF-051618-09
Calibrator Concentrate	102096	MF-060817-03
Calibrator Diluent	102094	MF-050718-03
Sample Diluent	102095	MF-050718-02

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	36 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.47
C	7.41
D	22.2
E	66.6
F	200
G	600
H	1800

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date 27 Jun 2018