

Product Number:	102082
Lot Number:	501288
Expiration:	18-Oct-18
Platform(s):	HD-1



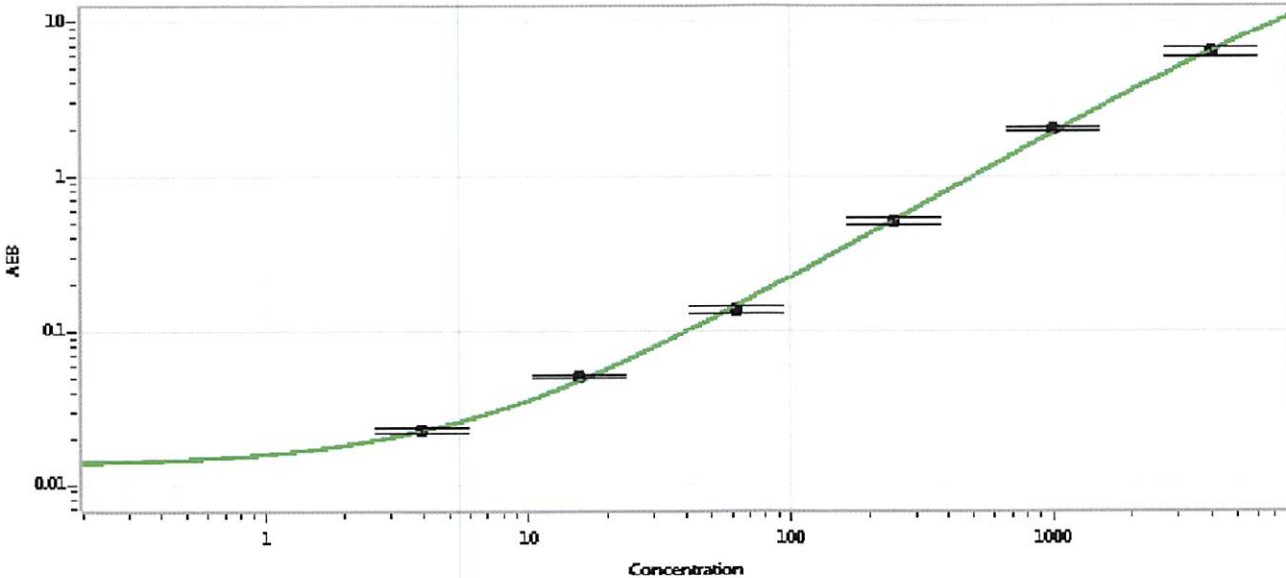
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102074	MF-061218-03
Detector Stock	102075	MF-061518-02
Detector Diluent	102306	MF-061318-04
Calibrator Concentrate	102078	MF-101816-11
Calibrator Diluent	102076	MF-061218-02
Sample Diluent	102077	MF-061118-08

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Concentrate	101957
RGP Reagent	103159

Calibrator Concentrate Target Value	40,000 pg/mL
-------------------------------------	--------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	3.91
C	15.6
D	62.5
E	250
F	1000
G	4000

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date 22 Jun 2018