

<b>Product Number:</b>	102082
<b>Lot Number:</b>	501046
<b>Expiration:</b>	8-Jul-18
<b>Platform(s):</b>	HD-1



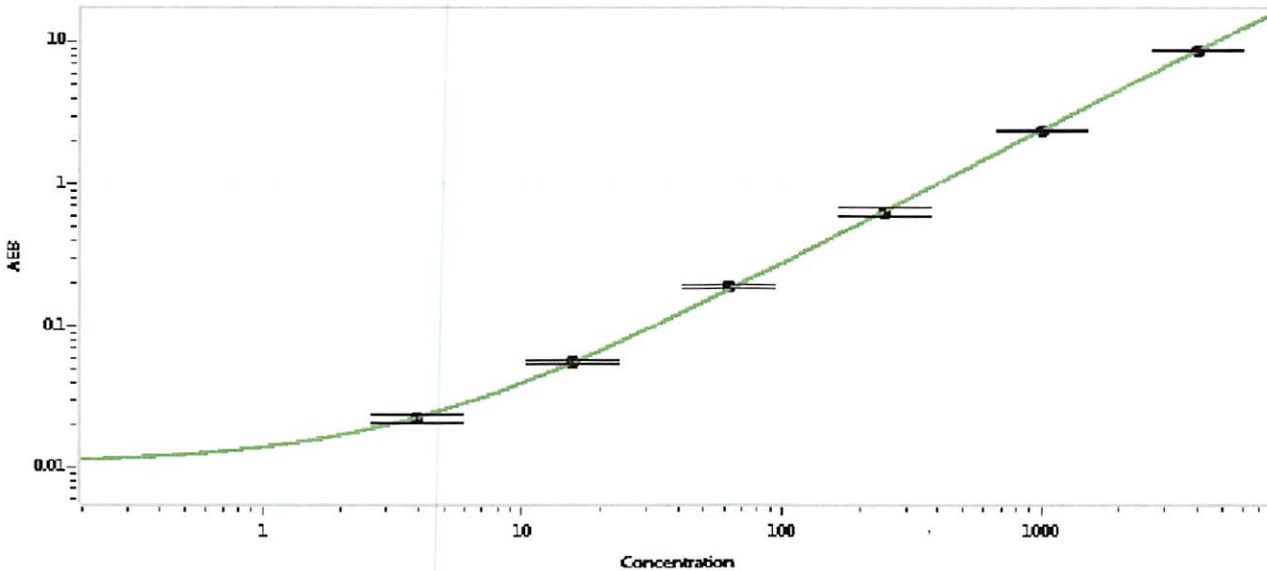
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102074	MF-011218-05
Detector Stock	102075	MF-010918-04
Detector Diluent	102306	MF-010818-11
Calibrator Concentrate	102078	MF-101816-11
Calibrator Diluent	102076	MF-010818-13
Sample Diluent	102077	MF-010918-01

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

<b>Calibrator Concentrate Target Value</b>	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]* 22 Jan 2018  
Signature/Date