

<b>Product Number:</b>	101855
<b>Lot Number:</b>	501087
<b>Expiration:</b>	29-May-18
<b>Platform(s):</b>	HD-1

Component	Part Number	Lot Number
Bead Stock	101902	MF-011818-01
Detector Stock	101903	MF-121417-03
Calibrator Concentrate	101850	MF-121517-07
Calibrator Diluent	101988	MF-112917-12
Sample Diluent	101989	MF-120417-11

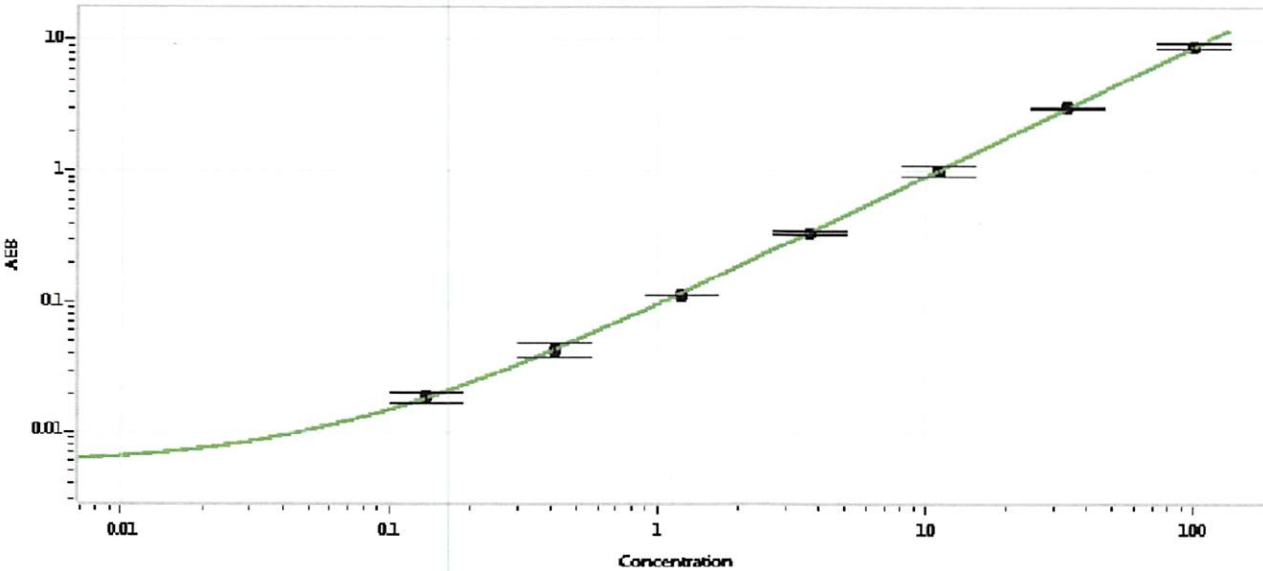


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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

<b>Calibrator Concentrate Target Value</b>	30 ng/mL
--	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.94
C	14.8
D	44.4
E	133
F	400
G	1200

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date 22 Jan 2018