

Product Number:	101824
Lot Number:	501135
Expiration:	24-Jul-2018
Platform(s):	HD-1



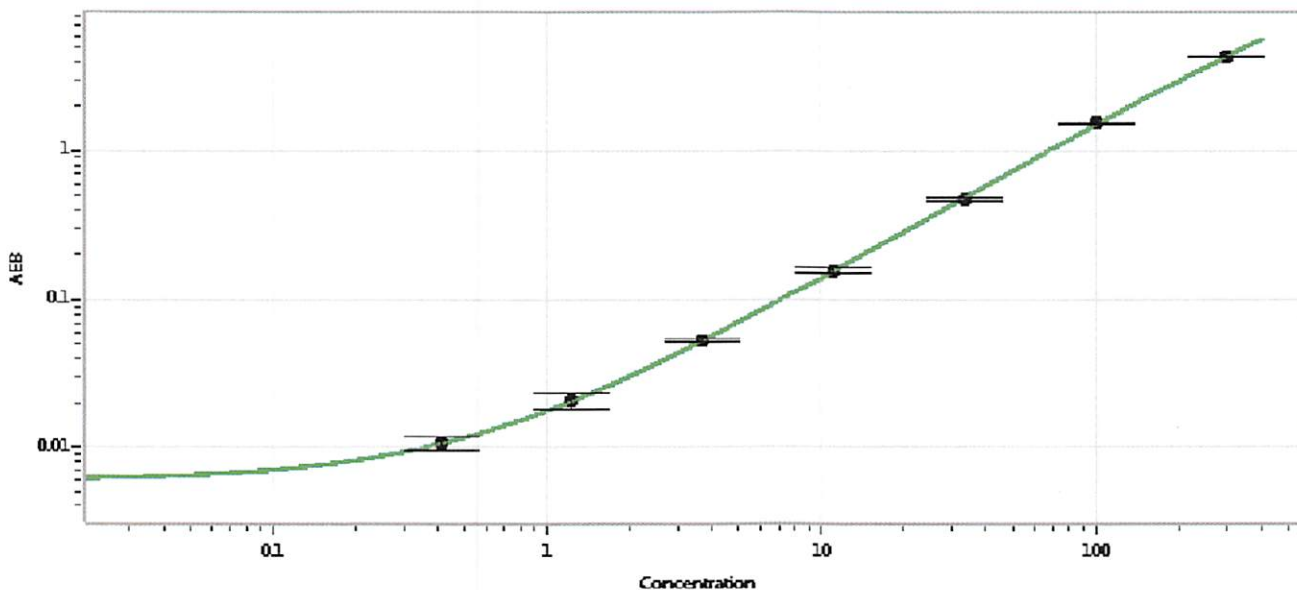
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101929	MF-021518-01
Detector Stock	101930	MF-012418-03
Calibrator Concentrate	101819	MF-030317-09
Calibrator Diluent	101943	MF-010818-12
Sample Diluent	101944	MF-091517-17

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	6000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 28/07/2018  
Signature/Date