

Product Number:	102517
Lot Number:	501163
Expiration:	26-Aug-2018
Platform(s):	HD-1



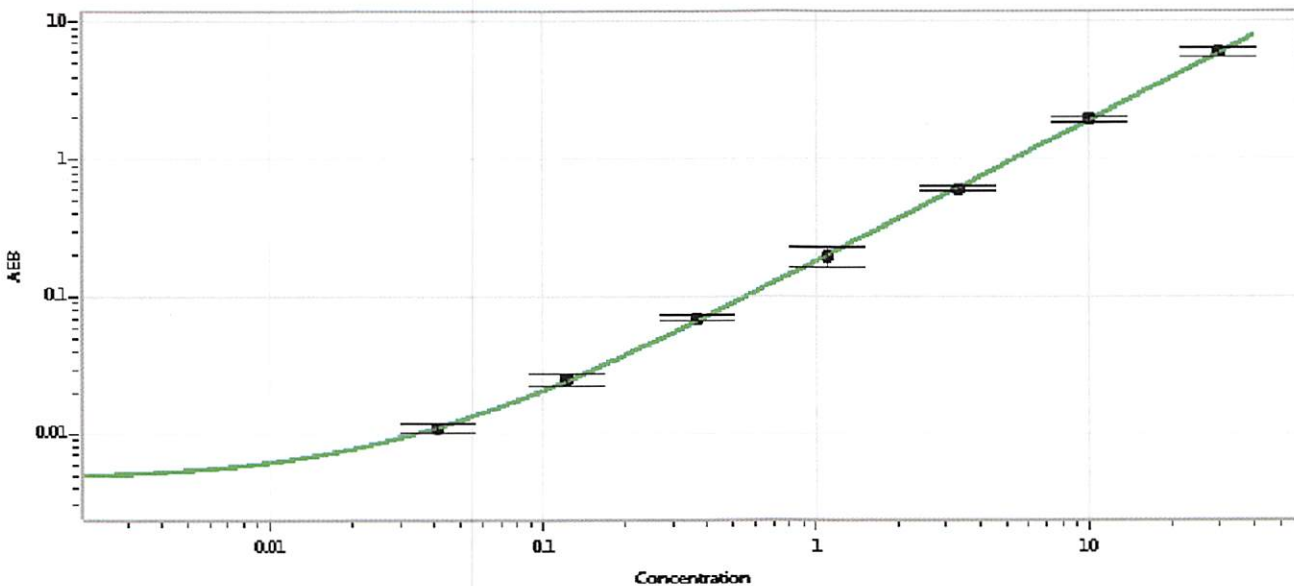
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102510	MF-030218-06
Detector Stock	102511	MF-030218-01
Calibrator Concentrate	102514	MF-022818-09
Calibrator Diluent	102512	MF-022618-02
Sample Diluent	102513	MF-022618-01

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	600 pg/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 07 MAR 2018
Signature/Date