

Product Number:	102583
Lot Number:	501362
Expiration:	16-Jan-2019
Platform(s):	HD-1



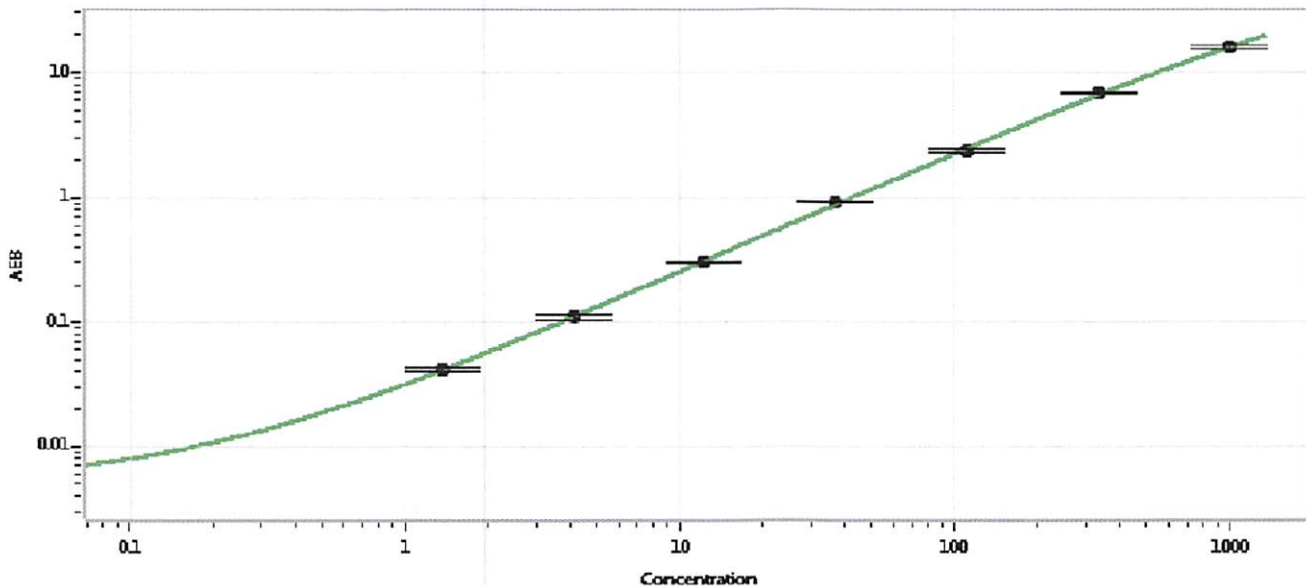
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102576	820802
Detector Stock	102577	819707
Calibrator Concentrate	102580	MF-051617-03
Calibrator Diluent	102578	816309
Sample Diluent	102579	819705

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	10 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval: Sivasankari Murugan Name, QC Supervisor Title, [Signature] Signature/Date, 13 AUG 2018 Date