

Product Number:	102318
Lot Number:	501344
Expiration:	11-Jan-2019
Platform(s):	HD-1



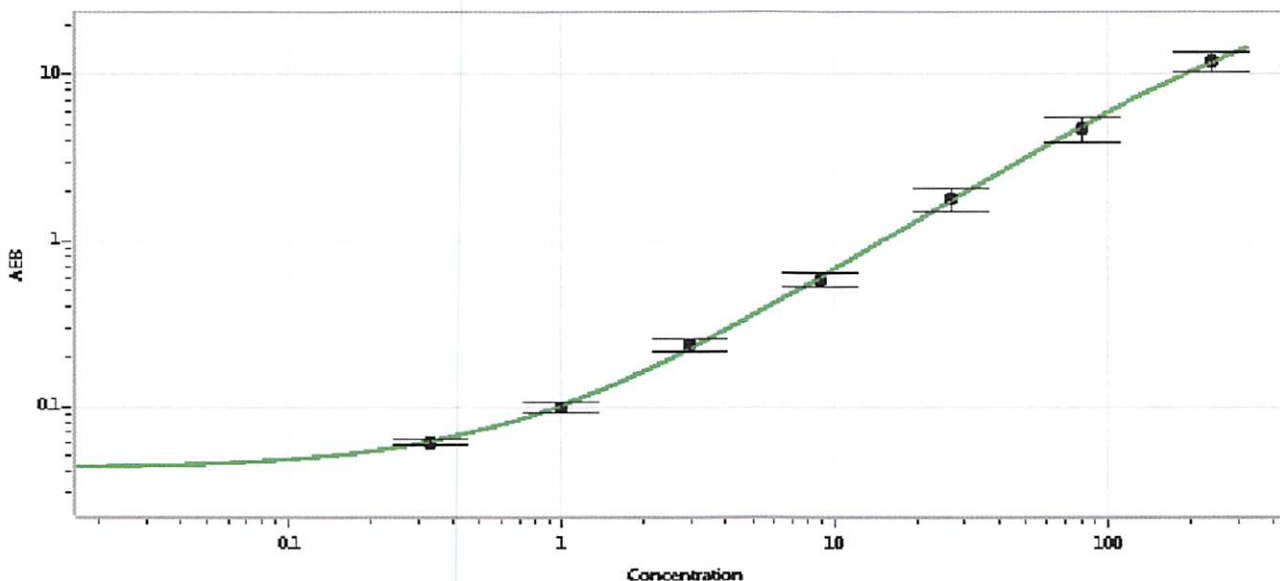
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	819806
Detector Stock	102310	820502
Calibrator Concentrate	102314	818414
Calibrator Diluent	102311	819201
Sample Diluent	102312	819202

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	4,800 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
2/6/2018
Signature/Date