

Product Number:	103071
Lot Number:	501315
Expiration:	3-Jan-2019
Platform(s):	HD-1 & SR-X



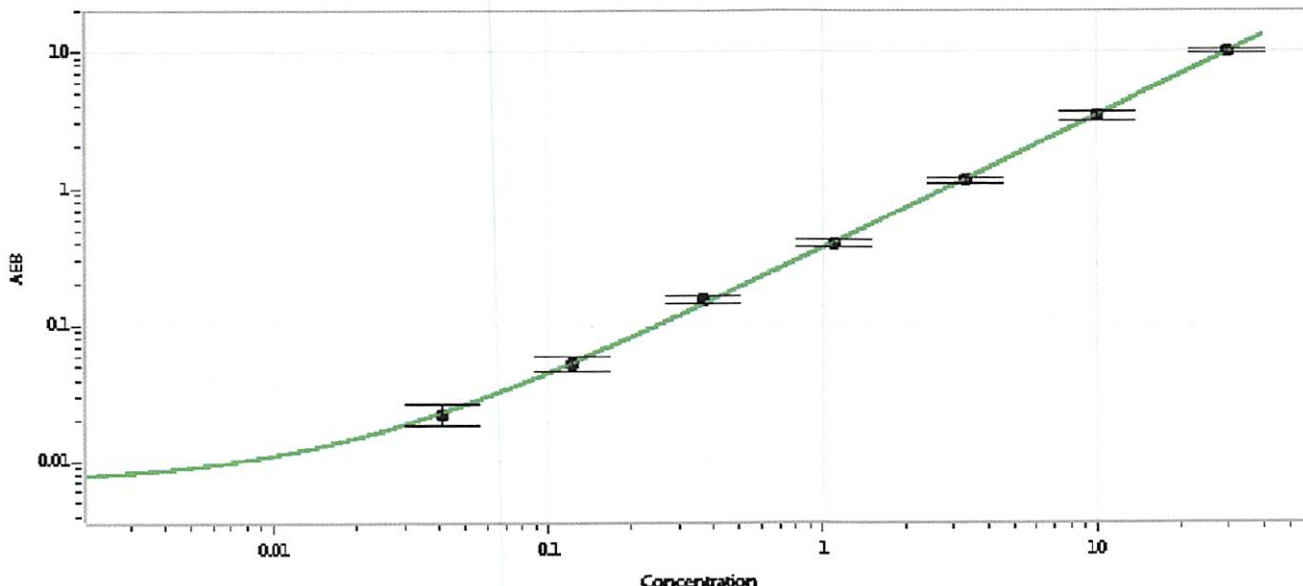
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	103064	818406
Detector Stock	103065	818409
Calibrator Concentrate	103068	MF-020418-01
Calibrator Diluent	103066	816305
Sample Diluent	103067	818003

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] 12 Jul 2018
Signature/Date