

Product Number:	103093
Lot Number:	501452
Expiration:	20-Feb-2019
Platform(s):	HD-1 & SR-X



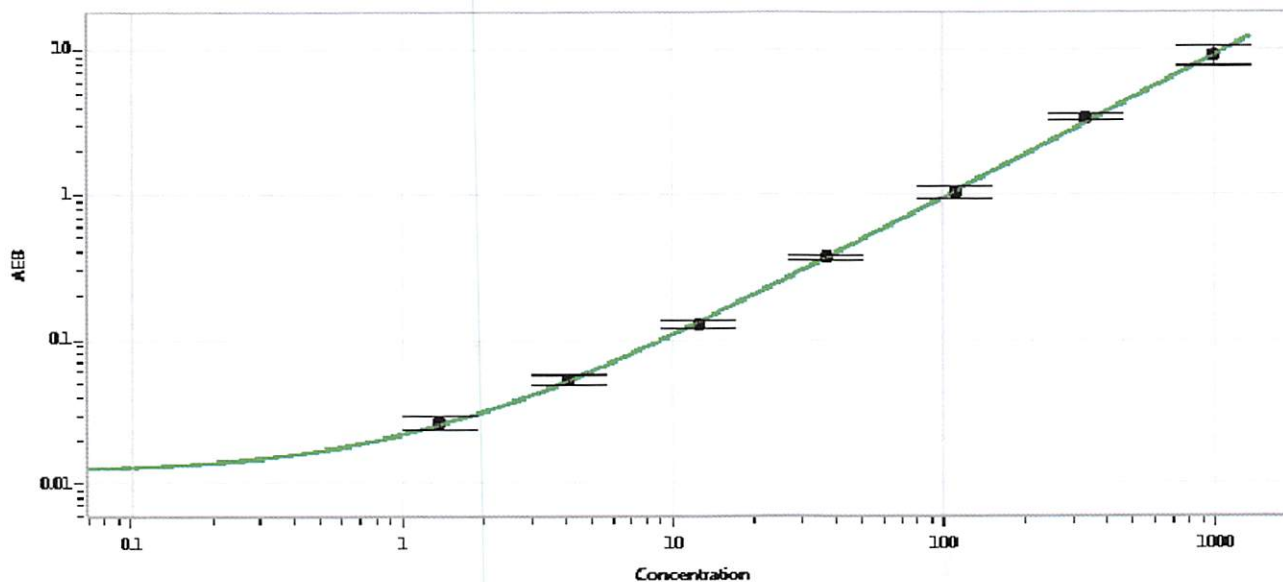
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102021	828505
Detector Stock	103043	828401
Calibrator Concentrate	103047	MF-010818-10
Calibrator Diluent	103044	828402
Sample Diluent	103046	828301

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	12 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]* 18 OCT 2018  
Signature/Date