

Product Number:	102635
Lot Number:	501426
Expiration:	26-Jan-2019
Platform(s):	HD-1



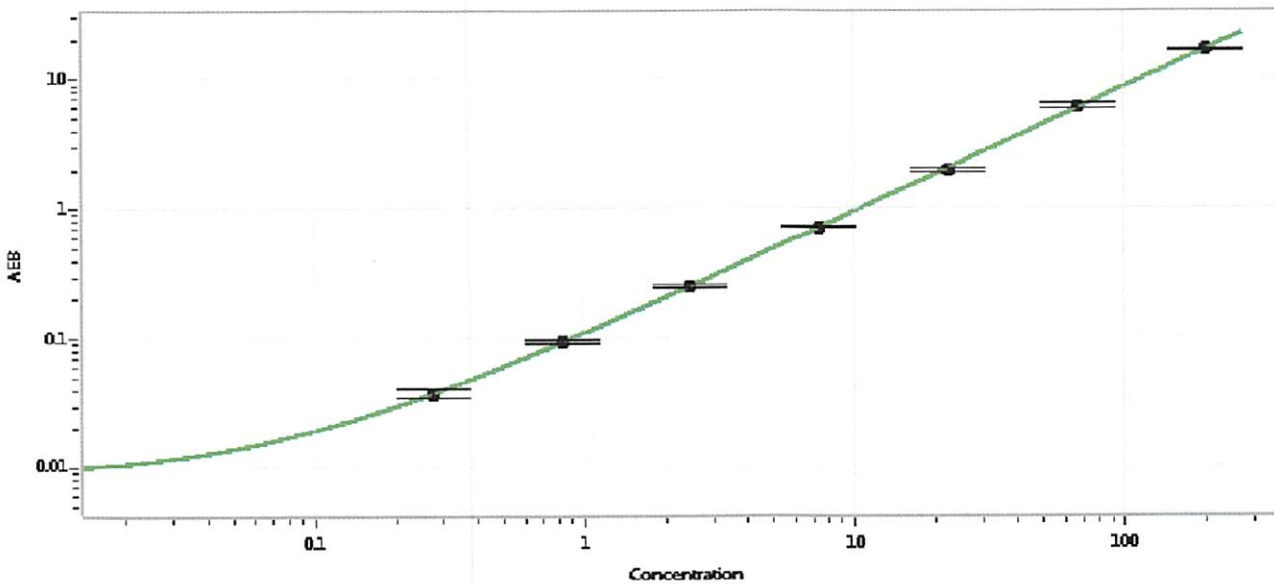
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102628	821106
Detector Stock	102629	820706
Calibrator Concentrate	102632	822101
Calibrator Diluent	102630	819812
Sample Diluent	102631	820803

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	4 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 26 SEP 2018  
Signature/Date