

Product Number:	102475
Lot Number:	501422
Expiration:	20-Feb-2019
Platform(s):	HD-1



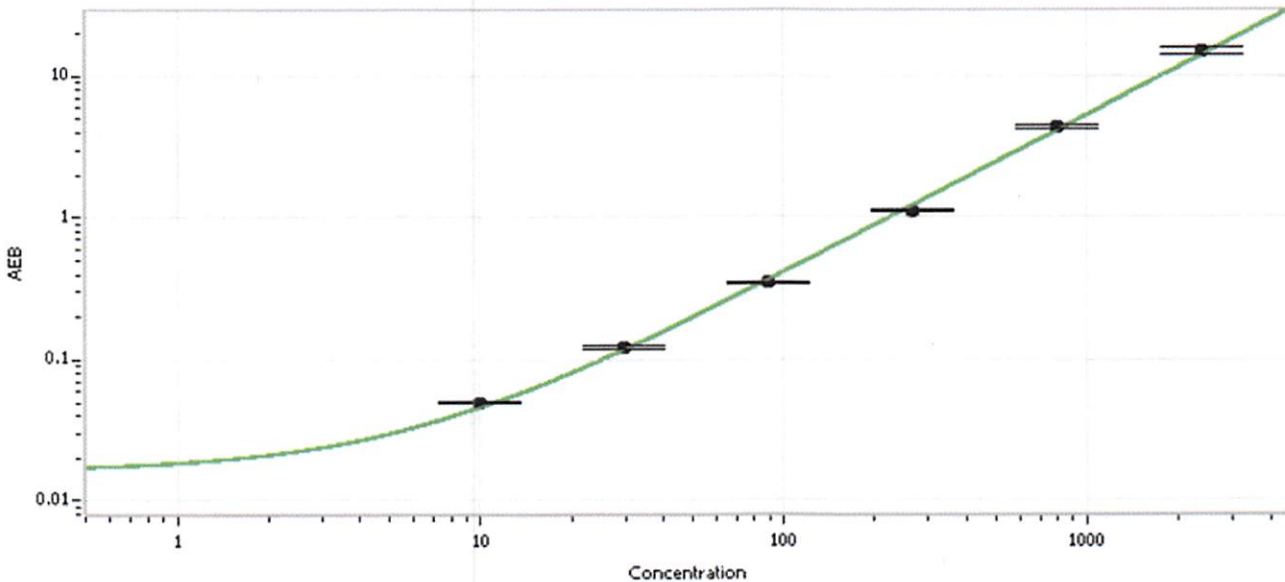
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102468	826401
Detector Stock	102469	826402
Calibrator Concentrate	102472	MF-062018-06
Calibrator Diluent	102470	823611
Sample Diluent	102471	826207

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	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.88
C	29.6
D	88.9
E	267
F	800
G	2400

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 2/5/2018
Signature/Date