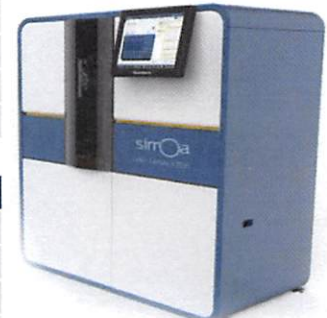


Product Number:	102700
Lot Number:	501414
Expiration:	30-Nov-2018
Platform(s):	HD-1



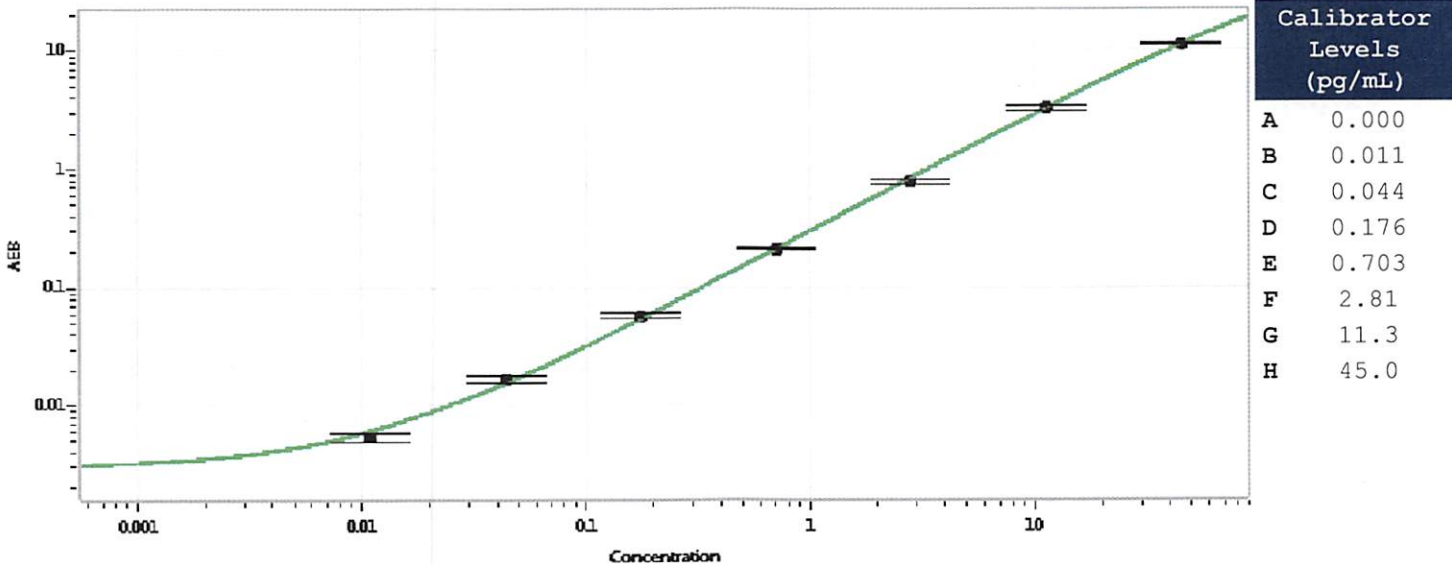
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102693	MF-053018-06
Detector Stock	102694	MF-060118-08
Calibrator Concentrate	102697	MF-060418-01
Calibrator Diluent	102695	MF-041018-20
Sample Diluent	102696	MF-041018-19

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	1 ng/mL
-------------------------------------	---------

Calibration Curve



Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

Sivasankari Murugan
Signature/Date
18SEP2018