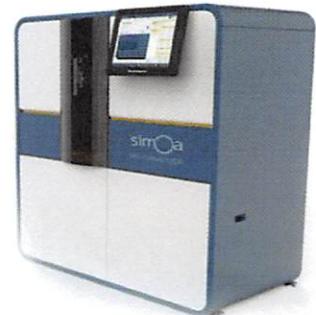


Product Number:	103093
Lot Number:	501124
Expiration:	1-May-2018
Platform(s):	HD-1 & SR-X

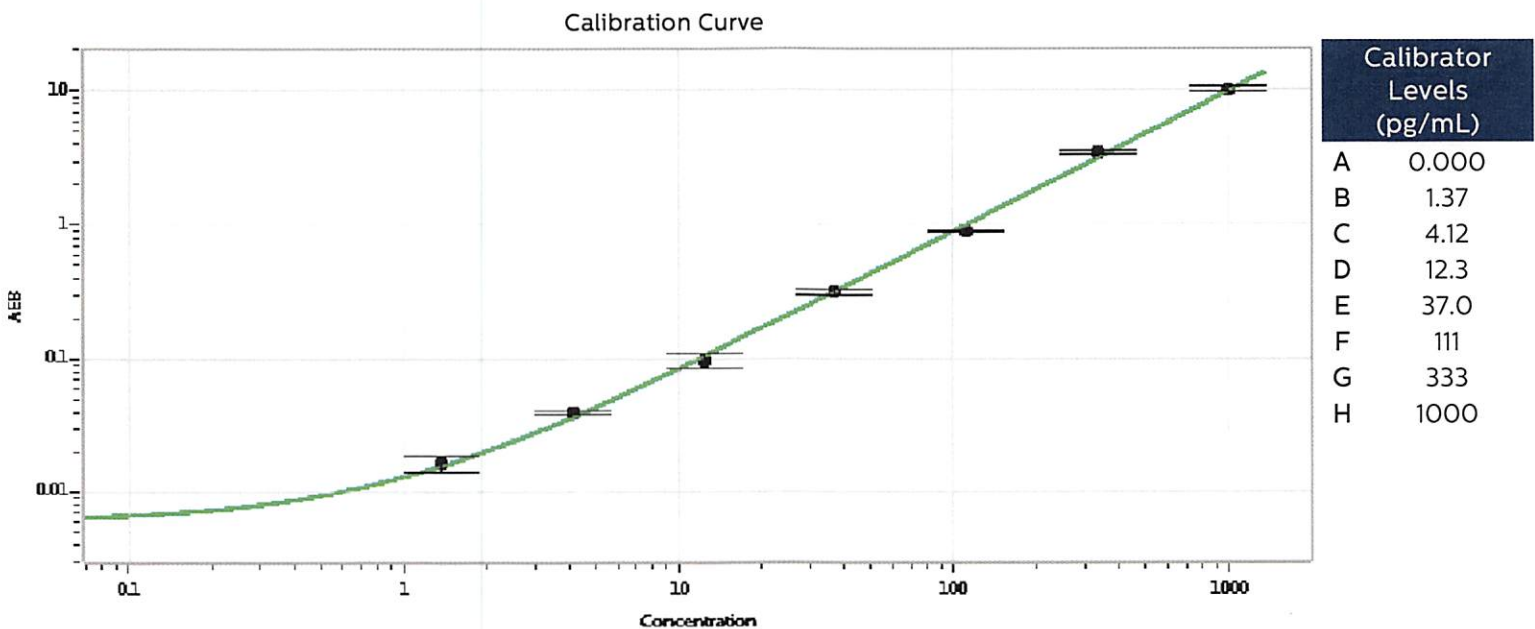


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102021	MF-122217-01
Detector Stock	103043	MF-121317-08
Calibrator Concentrate	103047	MF-010818-10
Calibrator Diluent	103044	MF-110117-06
Sample Diluent	103046	MF-103117-02

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	12 ng/mL
-------------------------------------	----------



Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 01 MAR 2018
Signature/Date