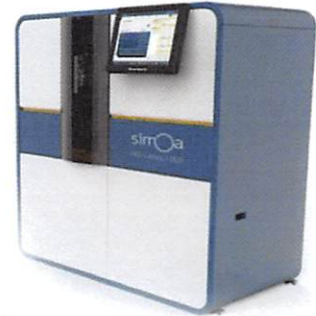


Product Number:	102336
Lot Number:	501146
Expiration:	19-Jul-2018
Platform(s):	HD-1

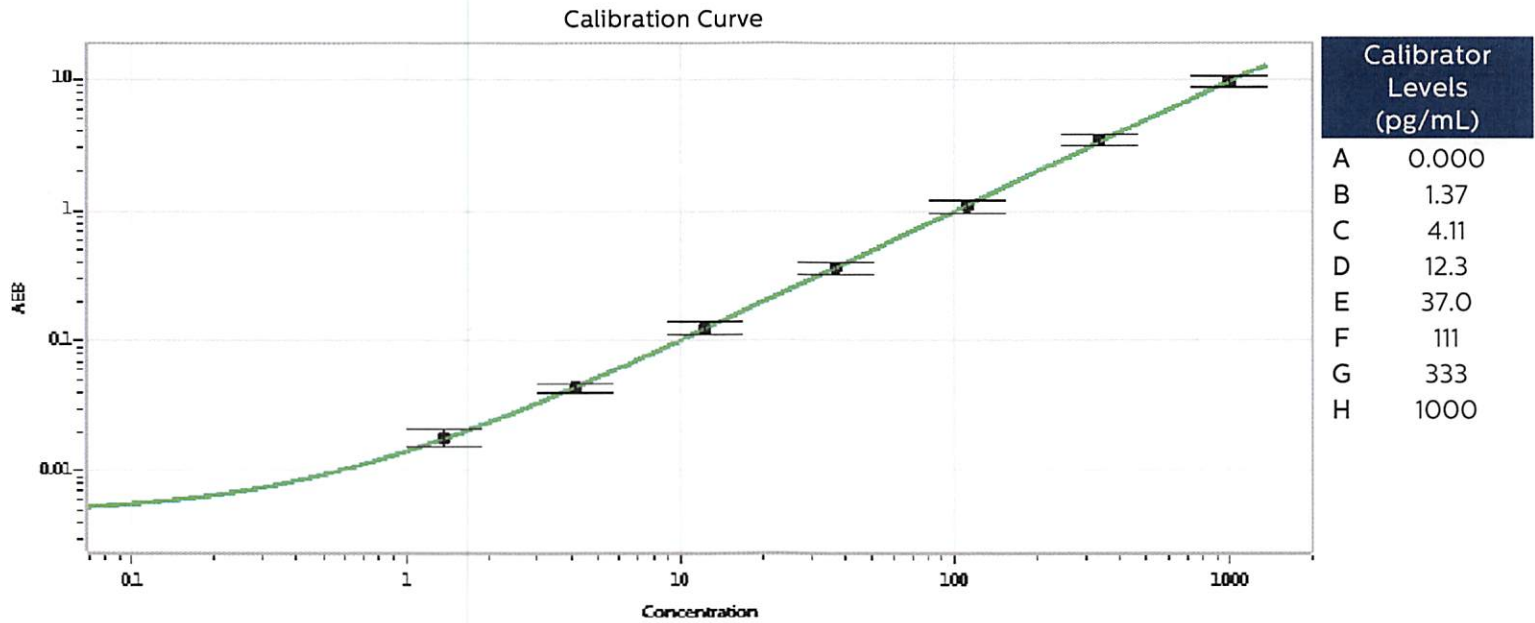


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	MF-022018-04
Detector Stock	102331	MF-022018-02
Calibrator Concentrate	102335	MF-092917-03
Calibrator Diluent	102332	MF-011918-10
Sample Diluent	102333	MF-022218-01

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	20 ng/mL
-------------------------------------	----------



Review/Approval: Sivasankari Murugan Name QC Supervisor Title [Signature] Signature/Date 01 MAR 2018