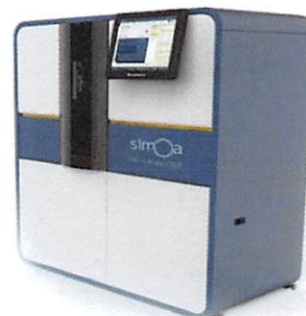


Product Number:	102109
Lot Number:	501232
Expiration:	2-Oct-2018
Platform(s):	HD-1

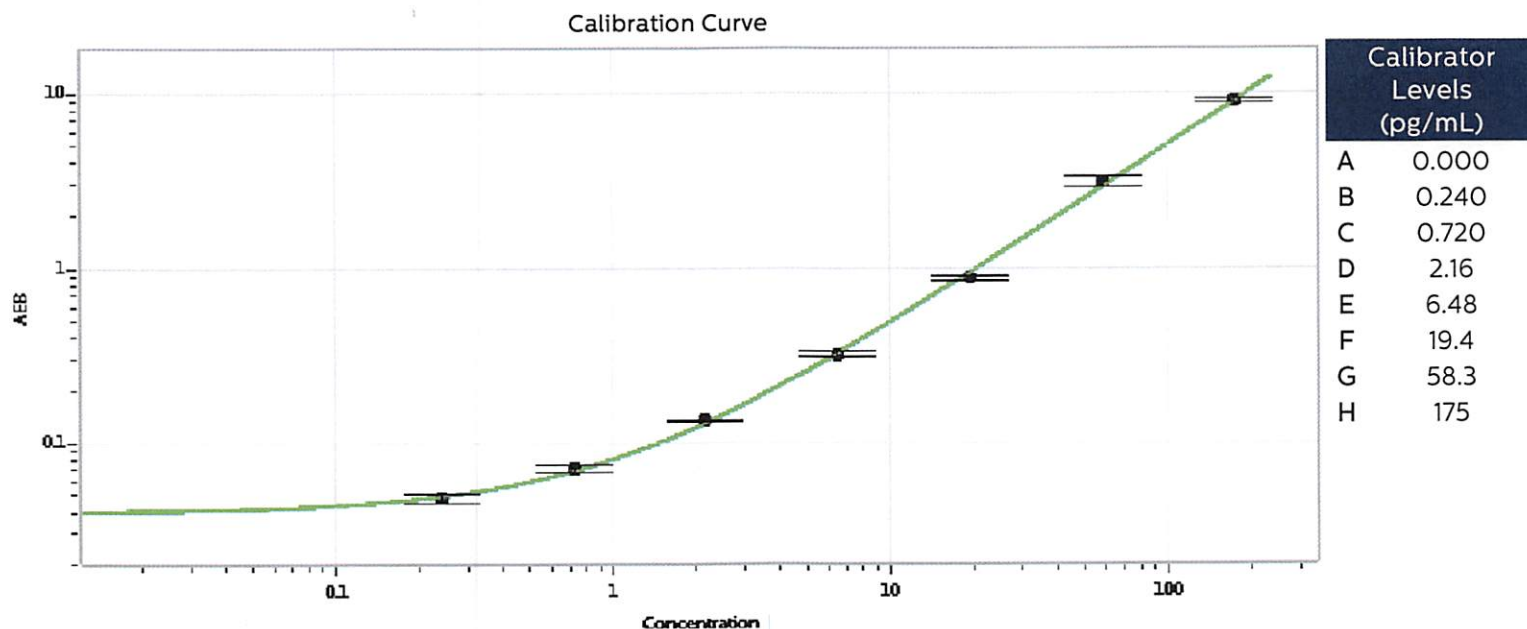
Component	Part Number	Lot Number
Bead Stock	102101	MF-052418-05
Detector Stock	102102	MF-040218-01
Calibrator Concentrate	102105	MF-040218-07
Calibrator Diluent	102103	MF-032018-06
Sample Diluent	102104	MF-012018-03



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

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	3500 pg/mL
-------------------------------------	------------



Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date 14 Jun 2018