

Product Number:	102635
Lot Number:	501172
Expiration:	7-Sep-2018
Platform(s):	HD-1

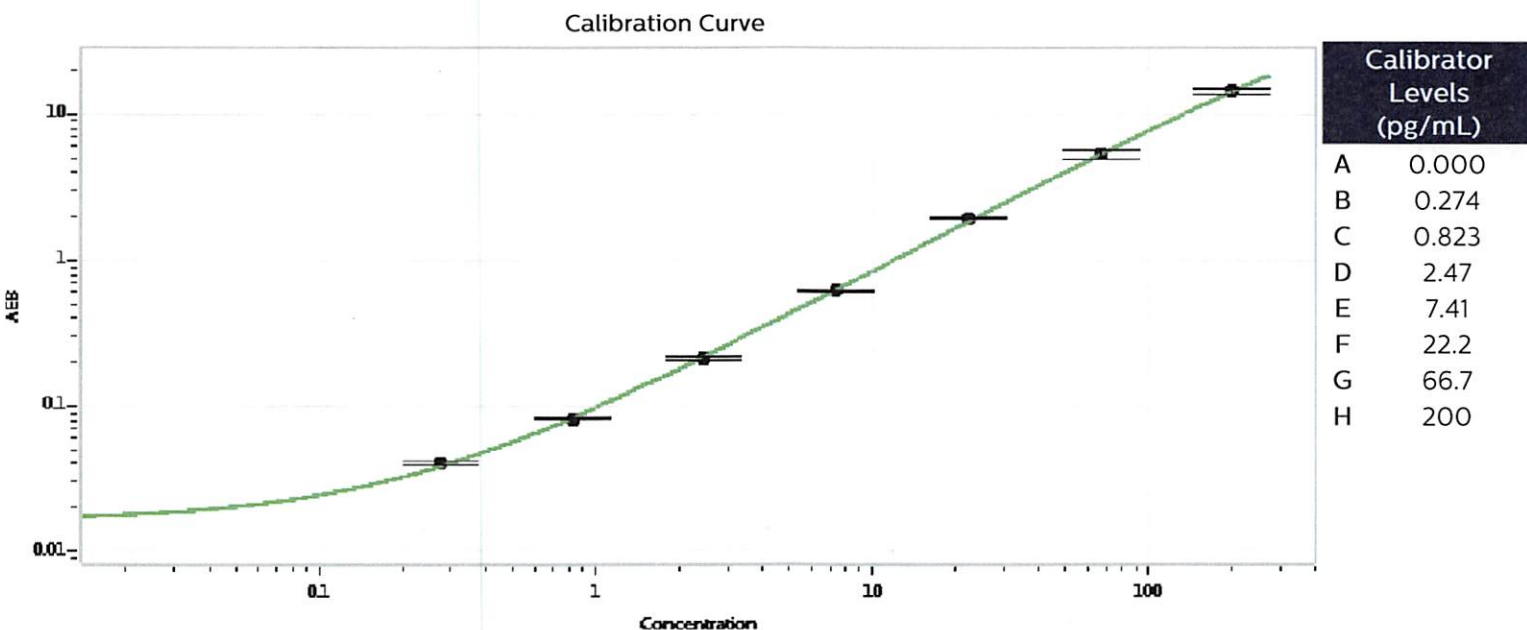
Component	Part Number	Lot Number
Bead Stock	102628	MF-031418-01
Detector Stock	102629	MF-030918-01
Calibrator Concentrate	102632	MF-032317-01
Calibrator Diluent	102630	MF-021618-09
Sample Diluent	102631	MF-030718-01



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	4 ng/mL
--	---------



Review/Approval	Sivasankari Murugan Name	QC Supervisor Title	Signature/Date
-----------------	-----------------------------	------------------------	----------------