

Product Number:	102713
Lot Number:	501236
Expiration:	16-Oct-2018
Platform(s):	HD-1

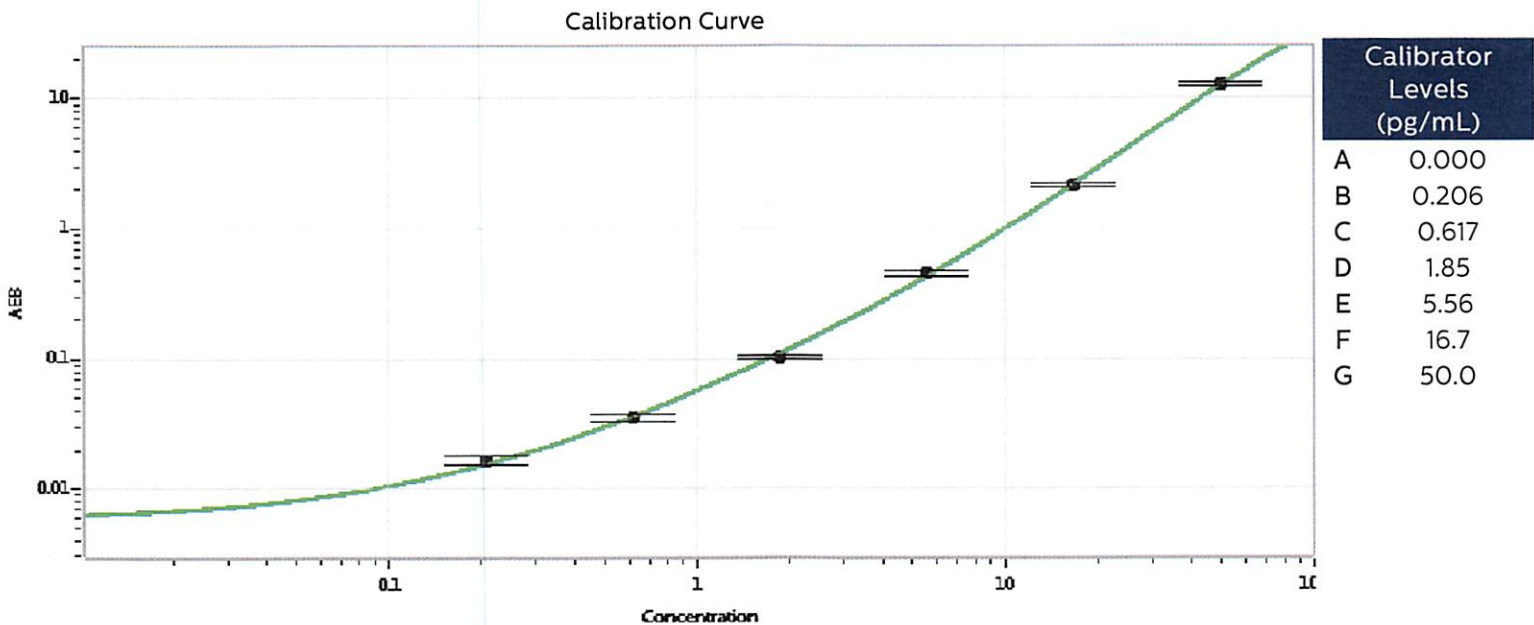


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102706	MF-043018-01
Detector Stock	102707	MF-041618-06
Calibrator Concentrate	102710	MF-041618-05
Calibrator Diluent	102708	MF-041018-15
Sample Diluent	102709	MF-041018-16

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	2 ng/mL
-------------------------------------	---------



Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date 15 May 2018