

Product Number:	102100
Lot Number:	501520
Expiration:	7-May-2019
Platform(s):	HD-1



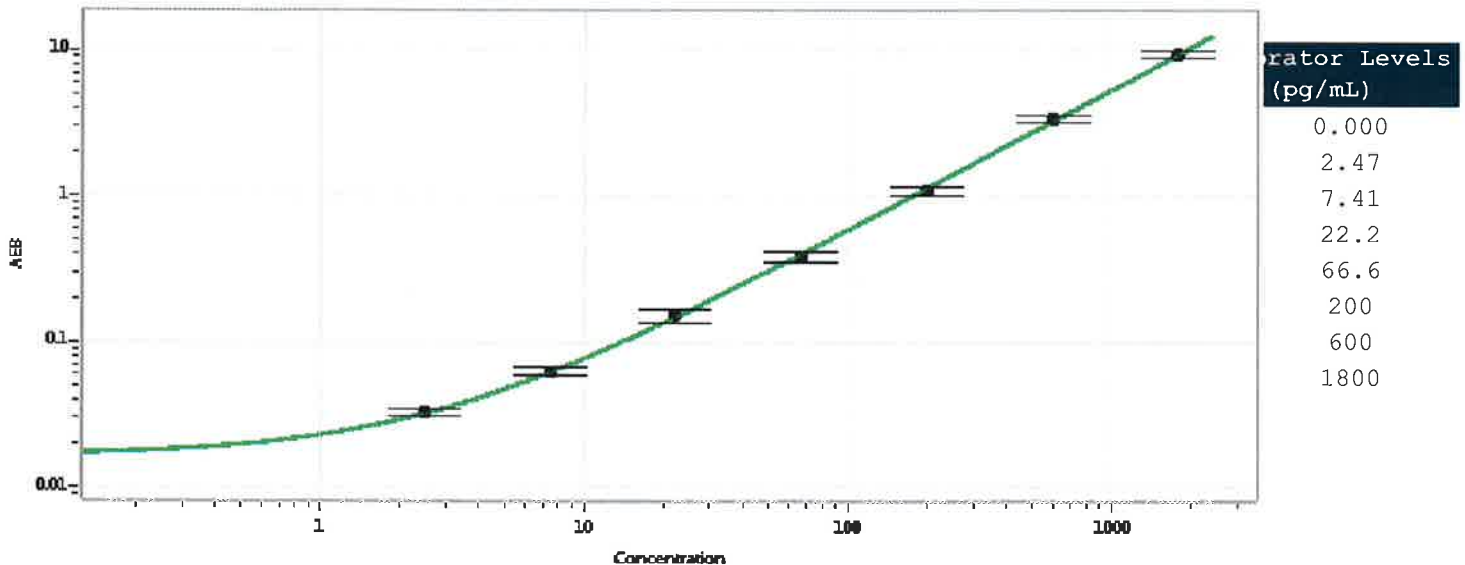
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102092	814103
Detector Stock	102093	813601
Calibrator Concentrate	102096	MF-060817-03
Calibrator Diluent	102094	812705
Sample Diluent	102095	812704

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	36 ng/mL
-------------------------------------	----------

Calibration Curve



Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 050602018
Signature/Date