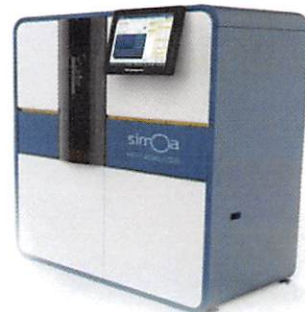


Product Number:	102656
Lot Number:	501164
Expiration:	12-Jul-2018
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102652	MF-011218-03
Detector Stock	102653	MF-030218-05
Detector Diluent	102843	MF-010518-13
Sample Diluent	102655	MF-010618-03
Reference Calibrators	102845	80600
Controls	102847 & 102848	80601



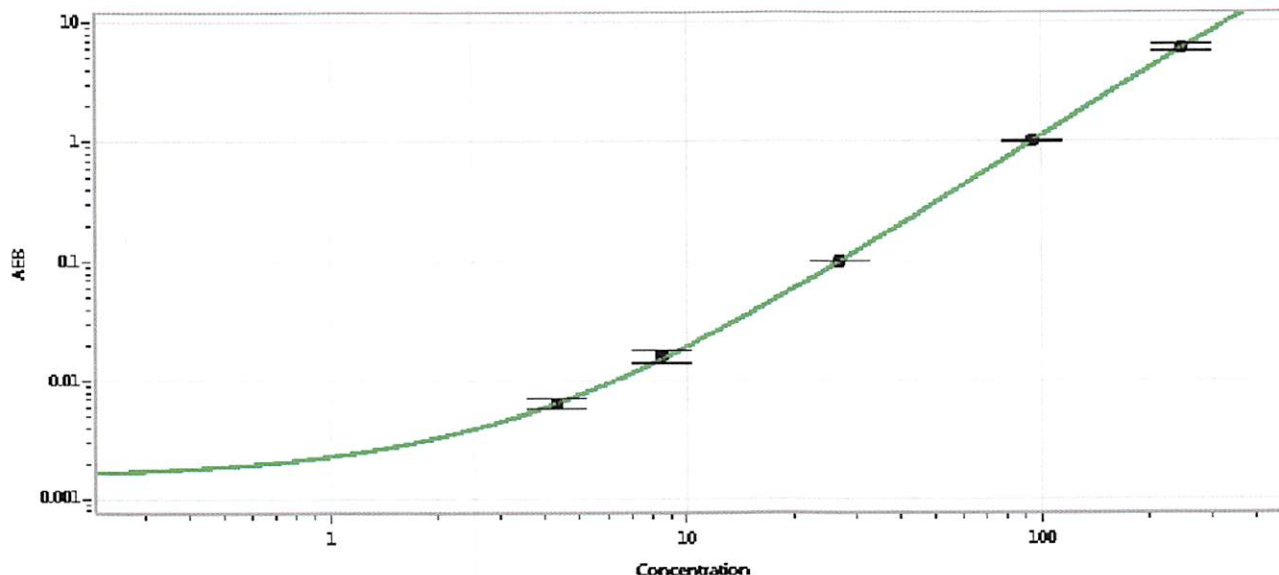
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Concentrate	101957
RGP Reagent	103159

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80601	4.67	4.34-7.08
Control 2	80601	155	129-193

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.33
C	8.53
D	26.9
E	94.6
F	248

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 16 MAR 2018
Signature/Date