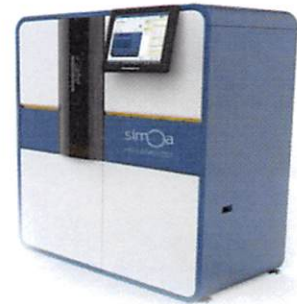


Product Number:	101588
Lot Number:	501456
Expiration:	10-Feb-2019
Platform(s):	HD-1 & SR-X



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

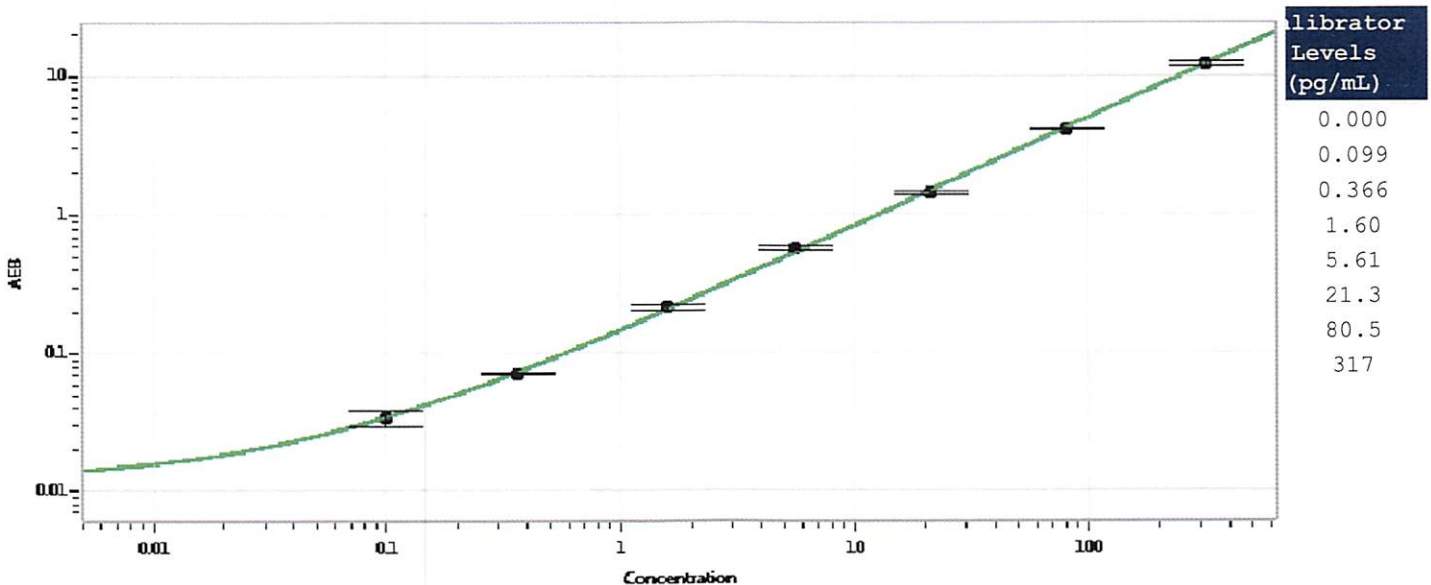
Component	Part Number	Lot Number
Bead Reagent	101590	822606
Detector Reagent	101589	827607
SBG Reagent	101591	823607
Sample Diluent	101592	822209
Calibrators	101110	80330
Controls	100551-100552	823402
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Troponin-I Advantage SR-X Kit Instructions (KI-0098).

	Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	823402	1.98	1.76-2.96	
Control 2	823402	127	110-164	
Panel 1	MF-031918-01	0.968	0.675-1.22	
Panel 2	MF-031918-02	5.53	4.84-7.27	
Panel 3	MF-031918-03	136	105-157	

Calibration Curve



Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 10/2/18  
Signature/Date