

Product Number:	101588
Lot Number:	501384
Expiration:	10-Feb-2019
Platform(s):	HD-1 & SR-X



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

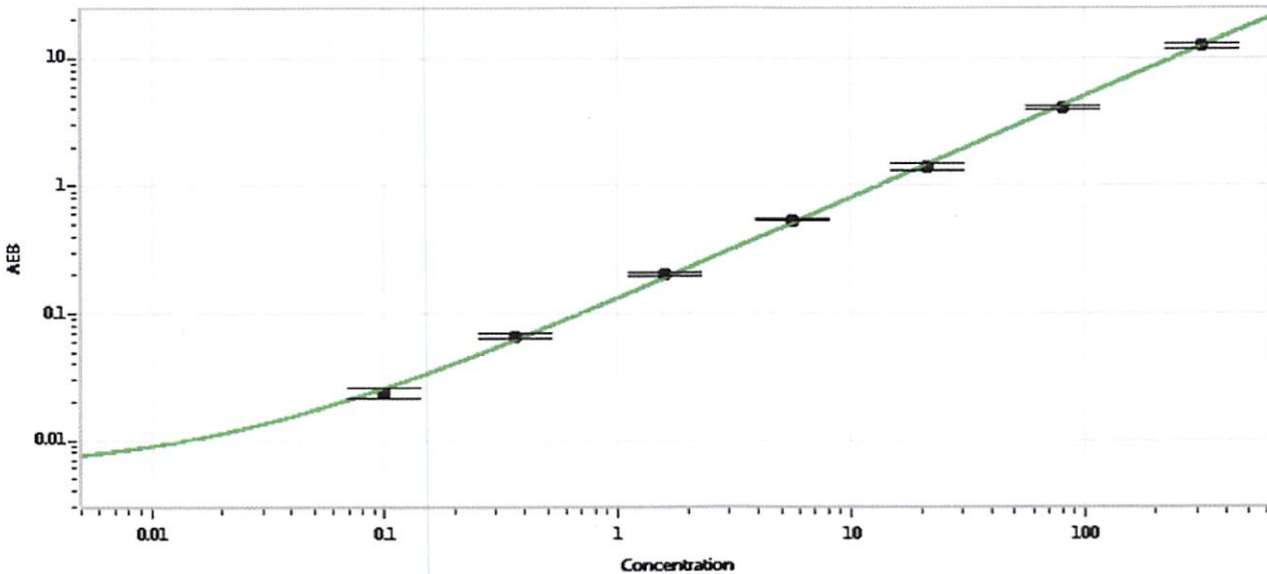
Component	Part Number	Lot Number
Bead Reagent	101590	822606
Detector Reagent	101589	822603
SBG Reagent	101591	823607
Sample Diluent	101592	822209
Calibrators	101110	80330
Controls	100551-100552	823402
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ 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	2.04	1.75-2.96	
Control 2	823402	128	110-164	
Panel 1	MF-031918-01	1.08	0.675-1.22	
Panel 2	MF-031918-02	6.60	4.84-7.27	
Panel 3	MF-031918-03	124	105-157	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.099
C	0.366
D	1.60
E	5.61
F	21.3
G	80.5
H	317

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date