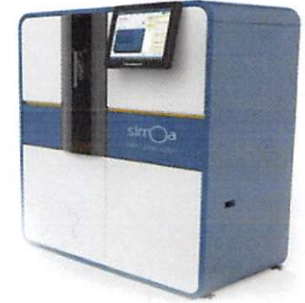


Product Number:	101588
Lot Number:	501787
Expiration:	26-Mar-2020
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101590	910813
Detector Reagent	101589	909802
SBG Reagent	101591	908515
Sample Diluent	101592	908516
Calibrators	101110	906404, 907901
Controls	100551-100552	907902
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

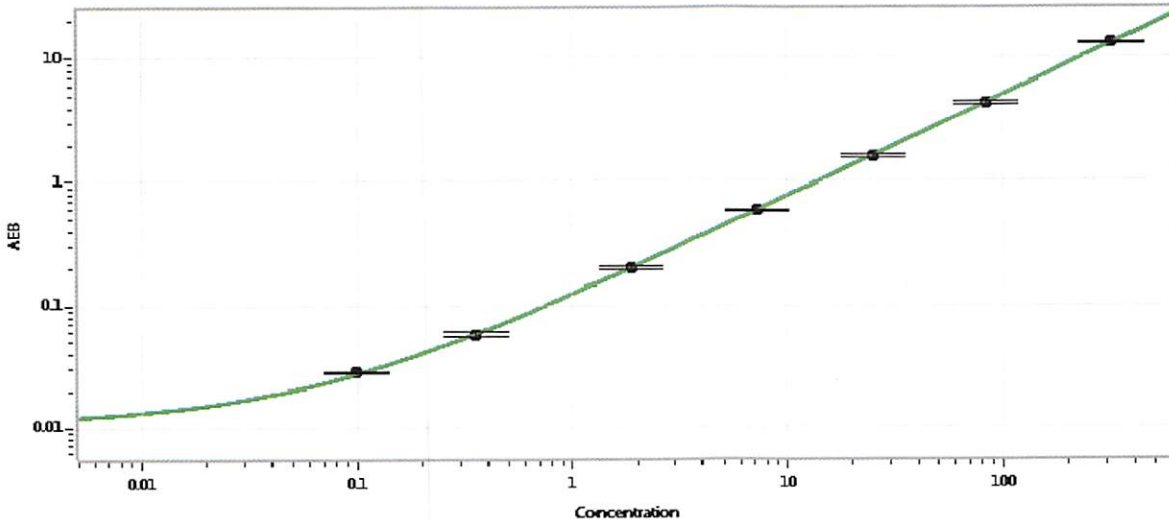
<sup>1</sup> Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	907902	3.60	2.80-4.20
Control 2	907902	166	135-202
Panel 1	MF-031918-01	0.934	0.642-1.22
Panel 2	MF-031918-02	5.95	4.78-7.17
Panel 3	MF-031918-03	115	91.2-160

\* 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	0.098
C	0.355
D	1.88
E	7.20
F	25.2
G	84.1
H	319

Review/Approval: Patrizia Stadler Assoc. Director, QC P. Stodler 25 Apr 2019

Name Title Signature/Date