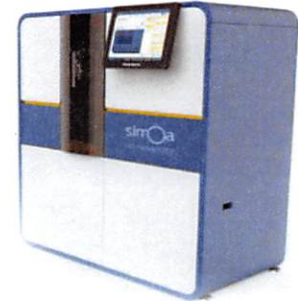


Product Number:	101588
Lot Number:	501763
Expiration:	16-Jan-2020
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101590	901608
Detector Reagent	101589	908801
SBG Reagent	101591	901910
Sample Diluent	101592	901607
Calibrators	101110	906404, 907901
Controls	100551-100552	907902
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

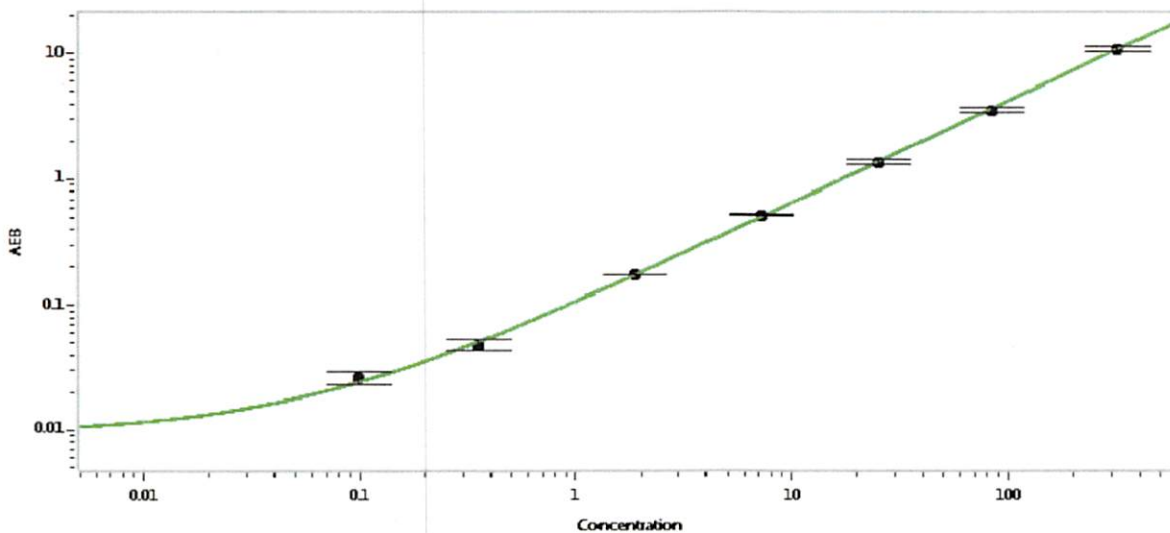
¹ Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	907902	3.51	2.79-4.18
Control 2	907902	159	134-204
Panel 1	MF-031918-01	0.820	0.634-1.23
Panel 2	MF-031918-02	5.42	4.78-7.17
Panel 3	MF-031918-03	99.5	92.0-162

* 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	<i>P. Stadler</i>	09 Apr 2019
	Name	Title	Signature/Date	