

Product Number:	101588
Lot Number:	501897
Expiration:	16-Jan-2020
Platform(s):	HD-1/HD-x & 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

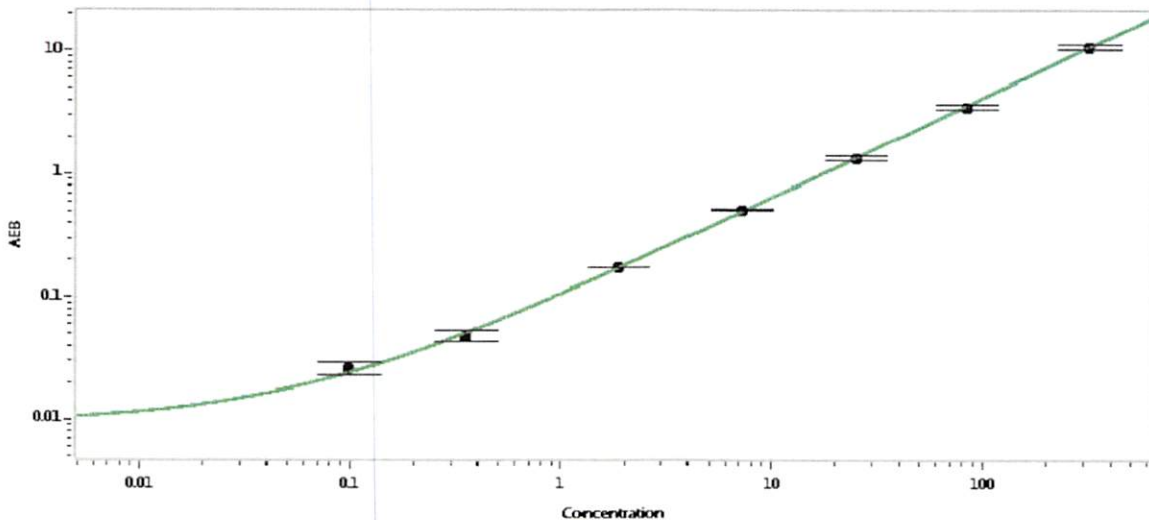
¹ RGP is not Kit Lot Specific

† 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

P. Stadler 17 Jul 2019

Name

Title

Signature/Date