

Product Number:	101588
Lot Number:	501979
Expiration:	13-Jun-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	920310
Detector Reagent	101589	920713
SBG Reagent	101591	920311
Sample Diluent	101592	920313
Calibrators	101110	922101, 922802
Controls	100551-100552	922102
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

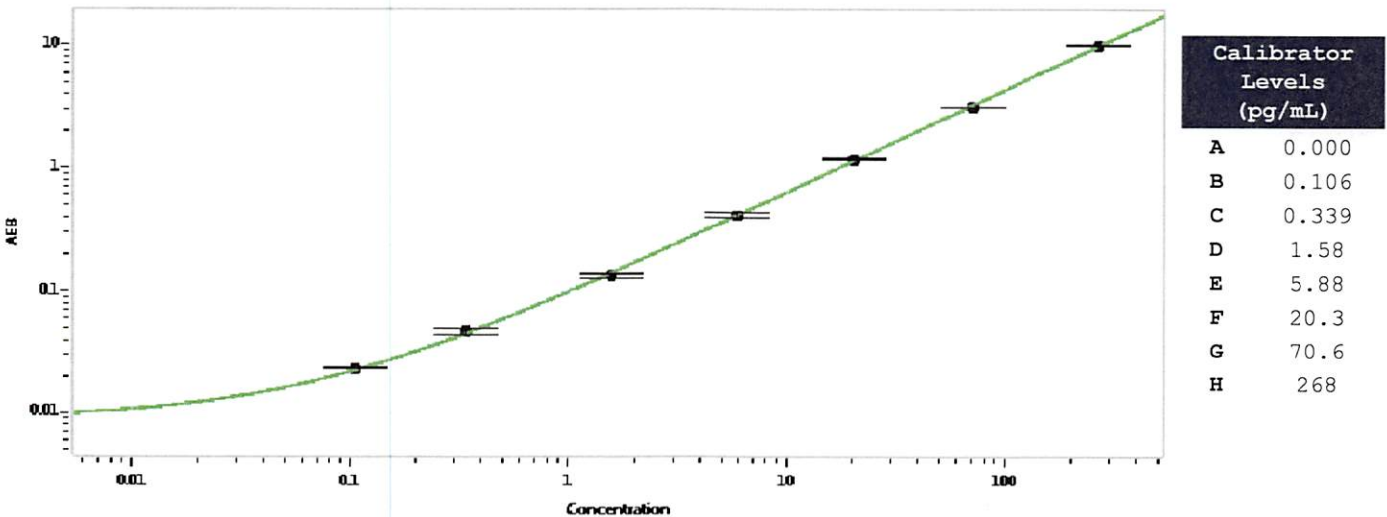
¹ RGP is not Kit Lot Specific

† 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	922102	3.47	2.52-4.37
Control 2	922102	138	111-181
Panel 1	MF-031918-01	0.895	0.642-1.22
Panel 2	MF-031918-02	5.31	4.78-7.17
Panel 3	MF-031918-03	98.1	91.2-160

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler
Signature/Date

29 Aug 2019