

Product Number:	101588
Lot Number:	502009
Expiration:	26-Mar-2020
Platform(s):	HD-1/HD-x & SR-X



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

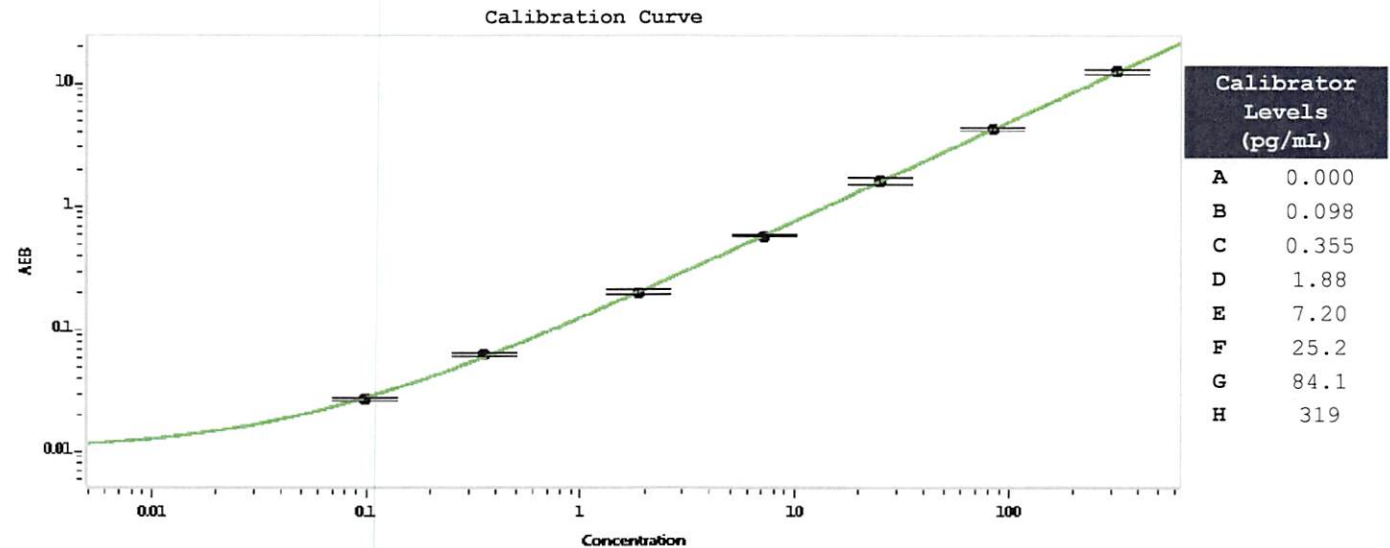
Component	Part Number	Lot Number
Bead Reagent	101590	910813
Detector Reagent	101589	909802
SBG Reagent	101591	908515
Sample Diluent	101592	908516
Calibrators	101110	922101, 907901
Controls	100551-100552	907902
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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	3.28	2.78-4.17
Control 2	907902	164	134-202
Panel 1	MF-031918-01	0.955	0.642-1.22
Panel 2	MF-031918-02	5.59	4.78-7.17
Panel 3	MF-031918-03	111	91.2-160

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



Review/Approval: Patrizia Stadler Assoc. Director, QC p.stadler 09 Sep 2019
 Name Title Signature/Date