

Product Number:	101588
Lot Number:	502541
Expiration:	23-May-2021
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	026604
Detector Reagent	101589	014508
SBG Reagent	101591	014403
Sample Diluent	101592	014402
Calibrators	101110	922101, 922802
Control 1	100551	922102
Control 2	100552	922102
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

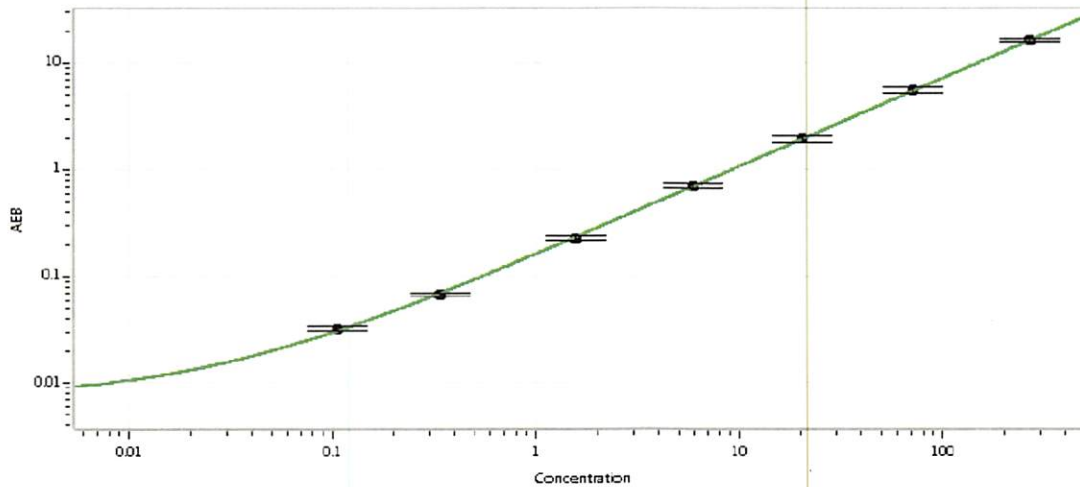
¹ RGP is not Kit Lot Specific

† 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	2.72	2.15-4.40
Control 2	922102	129	101-181
Panel 1	025501	1.35	1.04-2.56
Panel 2	025501	6.20	4.80-10.8
Panel 3	027401	161	137-274

* 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	0.106
C	0.339
D	1.58
E	5.88
F	20.3
G	70.6
H	268

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 21Oct2020
Signature/Date