

Product Number:	101588
Lot Number:	501073
Expiration:	8-Jul-2018
Platform(s):	HD-1



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

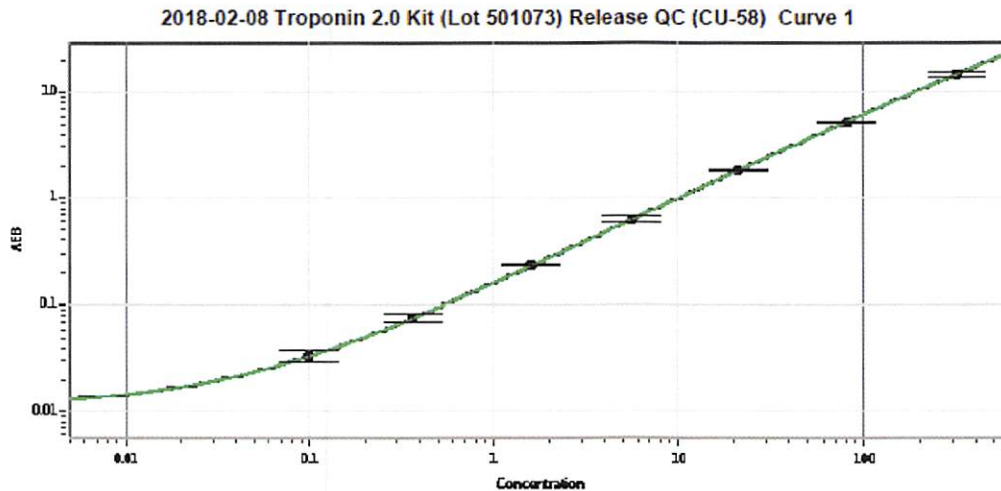
Component	Part Number	Lot Number
Bead Reagent	101590	80090
Detector Reagent	101589	80082
SBG Reagent	101591	80083
Sample Diluent	101592	80084
Calibrators	101110	80330
Controls	100551-100552	80331
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80331	0.219	0.111-0.300
Control 2	80331	142	121-202
Panel 1	MF-021716-05	2.15	1.66-2.50
Panel 2	MF-021716-06	5.16	4.18-6.27
Panel 3	MF-021716-07	124	95.7-144

* 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	0.099
C	0.366
D	1.60
E	5.61
F	21.3
G	80.5
H	317

Review/Approval

Sivasankari Murugan QC Supervisor
Name Title

[Signature] 16 FEB 2018
Signature/Date