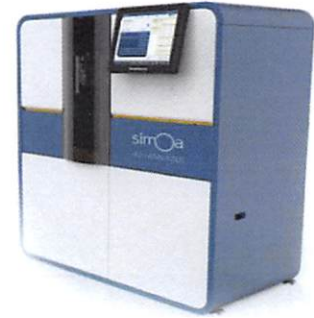


Product Number:	100906
Lot Number:	501434
Expiration:	24-Mar-2019
Platform(s):	HD-1



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

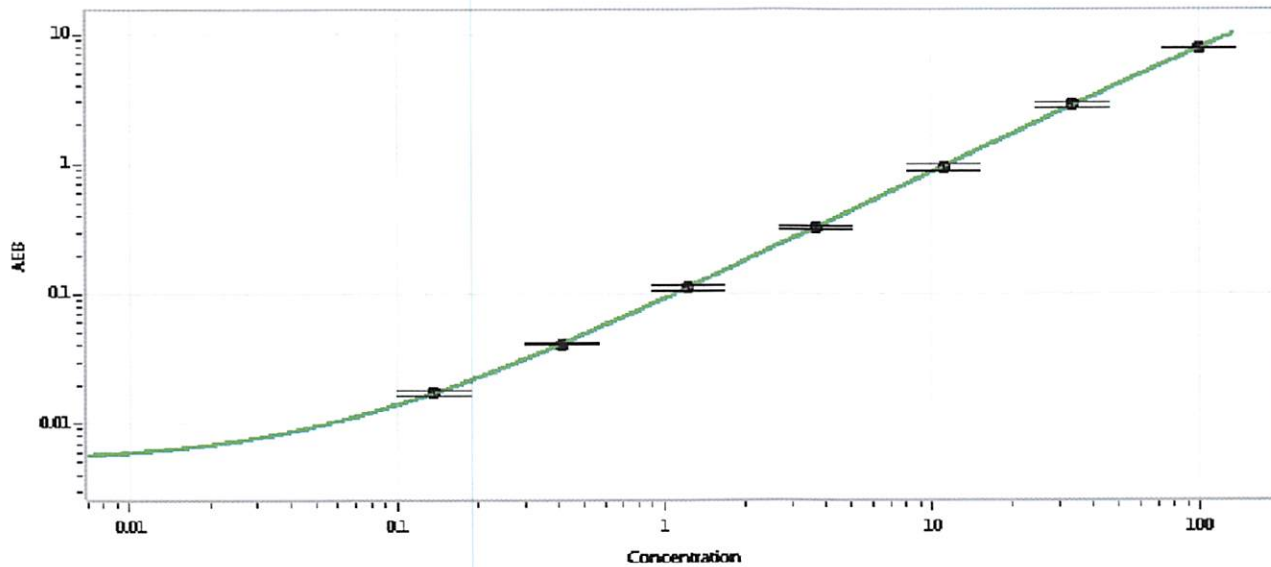
Component	Part Number	Lot Number
Bead Reagent	100911	827104
Detector Reagent	100908	826807
SBG Reagent	100914	826704
Sample Diluent	103321	826210
Calibrator Diluent	100918	823612
Calibrator Concentrate	100919	80200
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	MF-051717-02	1.98	1.77-2.65	
Control 2	MF-051717-02	50.3	38.7-58.1	
Panel 1	MF-032217-12	65.0	49.4-74.1	
Panel 2	MF-051817-02	42.9	32.9-53.6	
Panel 3	MF-051817-03	175	130-195	

Calibrator Concentrate Value Assignment (pg/mL)	2053
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 11/02/2018
Signature/Date