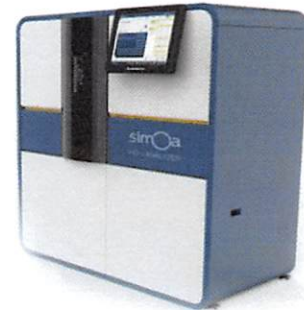


Product Number:	100198
Lot Number:	501254
Expiration:	23-Oct-2018
Platform(s):	HD-1



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

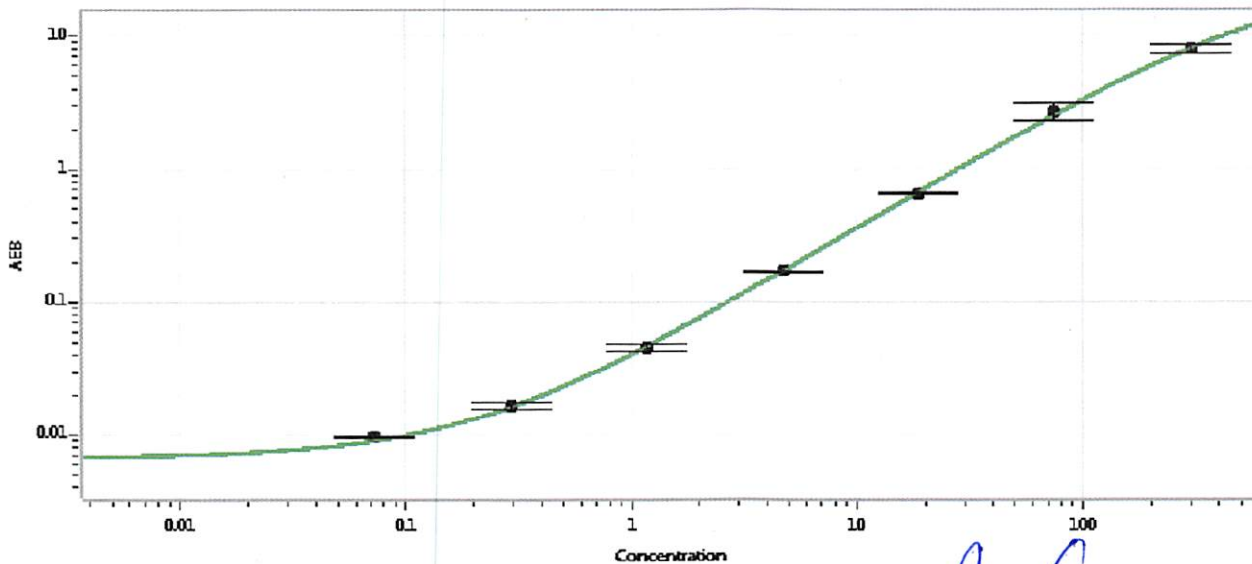
Component	Part Number	Lot Number
Bead Reagent	100815	81151
Detector Reagent	100816	81152
SBG Reagent	100817	81353
Sample Diluent	100818	81133
Calibrator Diluent	100829	81134
Calibrator Concentrate	100828	73380
RGP Reagent	103159	N/A ¹

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-091217-02	10.1	7.76-11.6	
Control 2	MF-091217-02	172	147-220	
Panel 1	MF-061617-01	5.94	3.46-6.45	
Panel 2	MF-061617-02	52.1	42.0-62.9	
Panel 3	MF-061617-03	339	279-459	

Calibrator Concentrate Value Assignment (pg/mL)	15,328
---	--------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.073
C	0.293
D	1.17
E	4.69
F	18.8
G	75.0
H	300

James Drummey
Senior Director, Manufacturing Operations
On behalf of:

Review/Approval

Sivasankari Murugan QC Supervisor
Name Title

[Signature]
May 18, 2018
Signature/Date