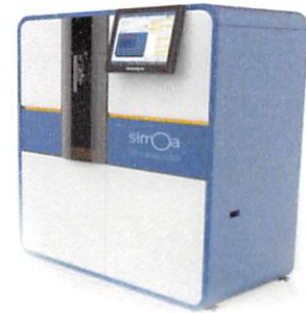


Product Number:	101212
Lot Number:	501117
Expiration:	29-Jul-2018
Platform(s):	HD-1



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

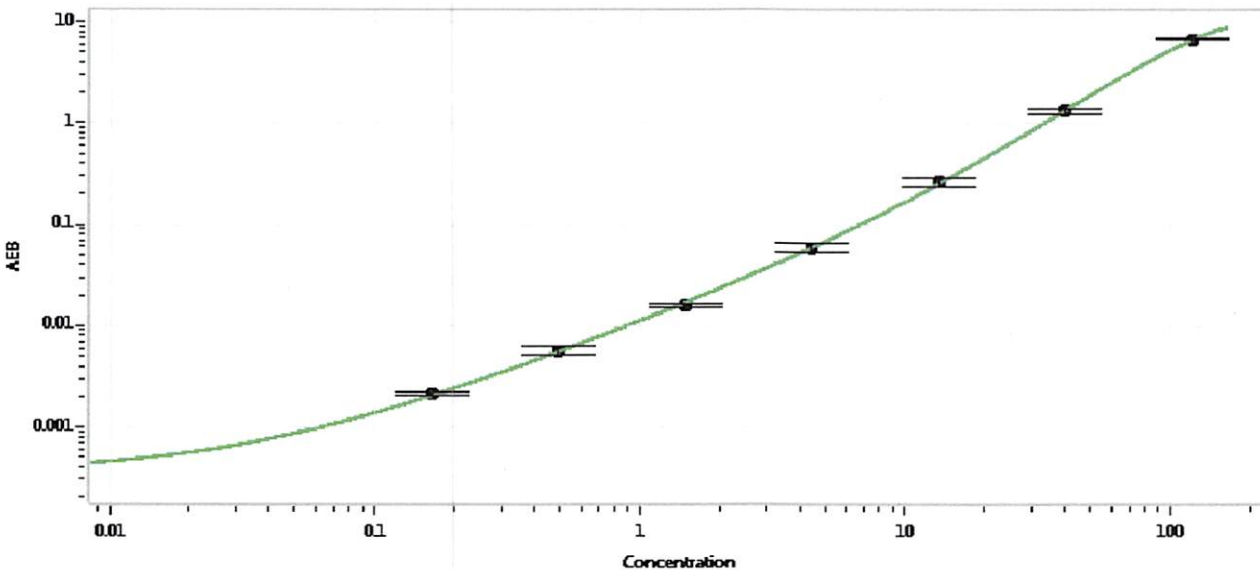
Component	Part Number	Lot Number
Bead Reagent	101198	80332
Detector Reagent	101200	80293
SBG Reagent	101211	80292
Sample Diluent	101209	80294
Calibrator Diluent	101204	80295
Calibrator Concentrate	101202	72780
RGP Reagent	103159	N/A ¹

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 2	MF-111617-02	69.9	55.6-83.4	
Control 3	72780-C2	244	206-339	
Panel 1	2005	84.7	64.3-143	
Panel 3	1984	44.4	33.1-70.9	

Calibrator Concentrate Value Assignment (pg/mL)	2032
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
15 FEB 2018