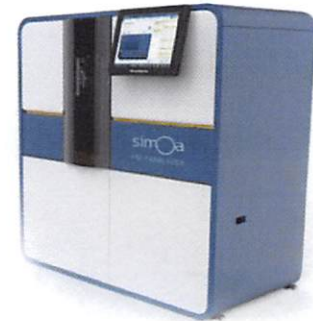


Product Number:	100196
Lot Number:	501410
Expiration:	12-Mar-2019
Platform(s):	HD-1



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

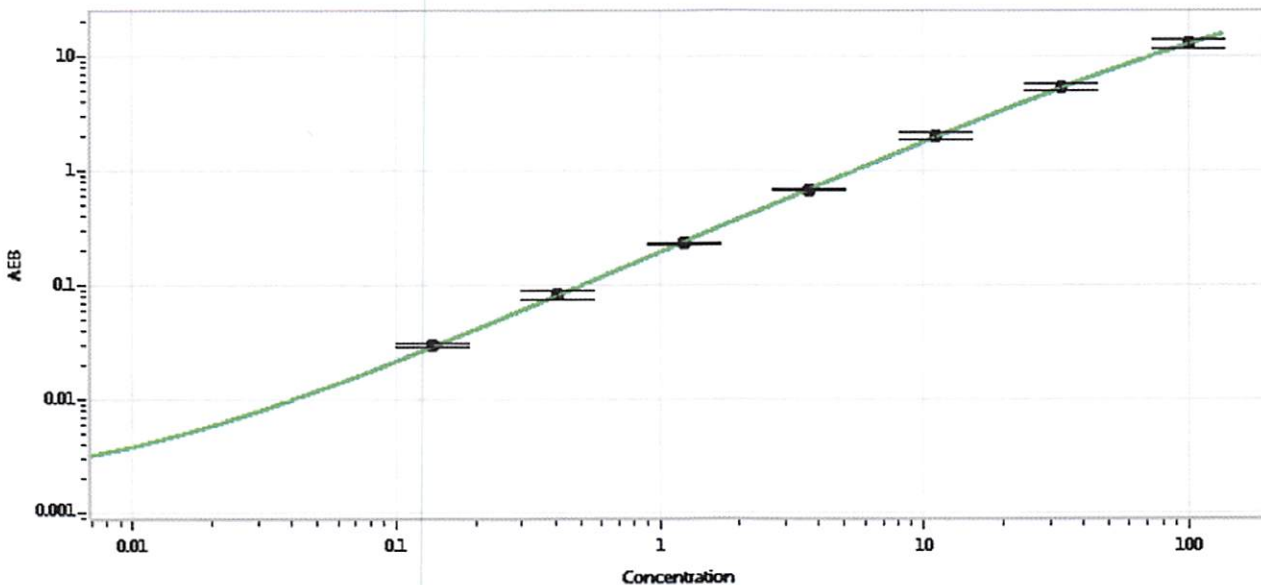
Component	Part Number	Lot Number
Bead Reagent	100391	825607
Detector Reagent	100315	825510
SBG Reagent	100316	825502
Sample Diluent	100317	818011
Calibrator Diluent	100665	825507
Calibrator Concentrate	100664	72701
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	52521	97.4	75.6-118	
Control 2	52521	9.07	7.72-11.6	
Panel 1	MF-012116-05	1.02	0.871-1.73	
Panel 2	MF-090815-18	6.73	4.45-9.13	
Panel 3	MF-090815-19	83.6	53.9-94.9	

Calibrator Concentrate Value Assignment (pg/mL)	2154
---	------

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

*S. Murugan* 20SEP2018  
Signature/Date