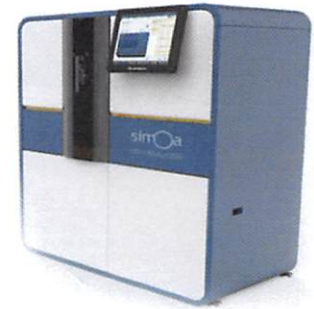


Product Number:	101664
Lot Number:	501401
Expiration:	29-Aug-2019
Platform(s):	HD-1 & SR-X



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

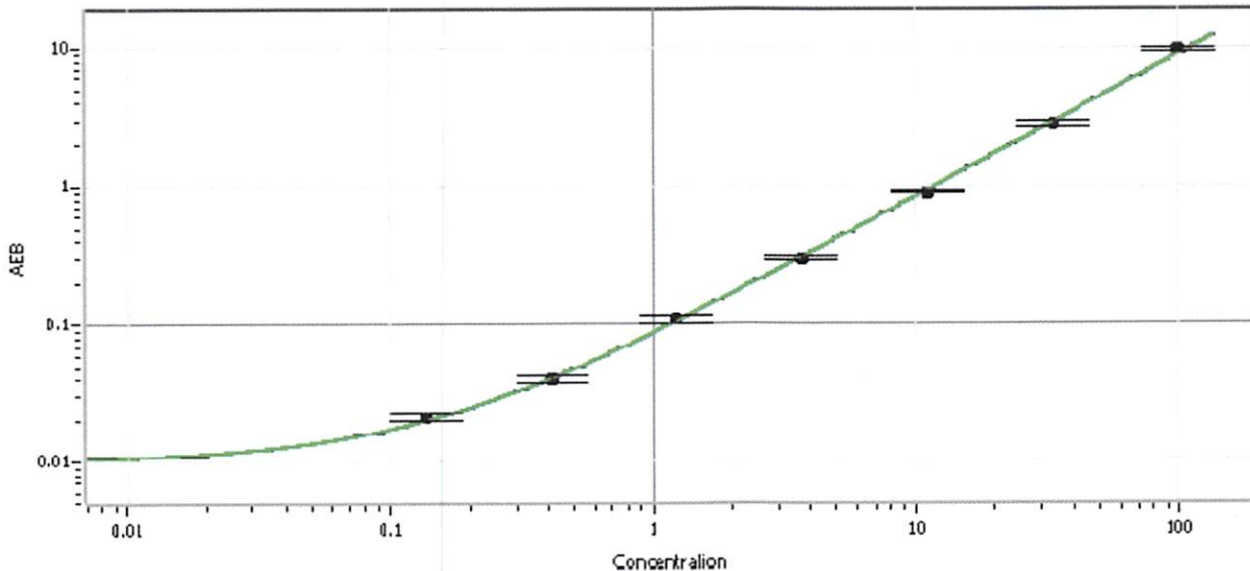
Component	Part Number	Lot Number
Bead Reagent	101666	825001
Detector Reagent	101665	824706
SBG Reagent	101667	824707
Sample Diluent	101668	824111
Calibrator Diluent	101669	824305
Calibrator Concentrate	103217	820504
RGP Reagent	103159	N/A <sup>1</sup>

<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	52520	1.11	0.740-1.33	
Control 2	52520	32.2	27.1-40.6	
Panel 1	MF-082015-01	1.18	0.942-1.53	
Panel 2	MF-081415-02	8.29	7.46-11.2	
Panel 3	MF-081415-03	10.4	9.20-15.0	
Panel 4	MF-090315-09	40.8	27.6-66.8	

Calibrator Concentrate Value Assignment (pg/mL)	2565
---	------

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 QC Supervisor  
Name Title

*[Signature]* 11 SEP 2018  
Signature/Date