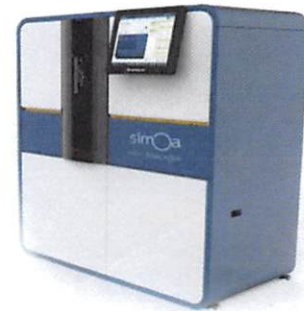


Product Number:	100988
Lot Number:	501319
Expiration:	3-Jan-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	100974	819207
Detector Reagent	100976	818401
SBG Reagent	100987	818402
Sample Diluent	100985	818002
Calibrator Diluent	100980	818404
Calibrator Concentrate	100978	73396
RGP Reagent	103159	N/A <sup>1</sup>

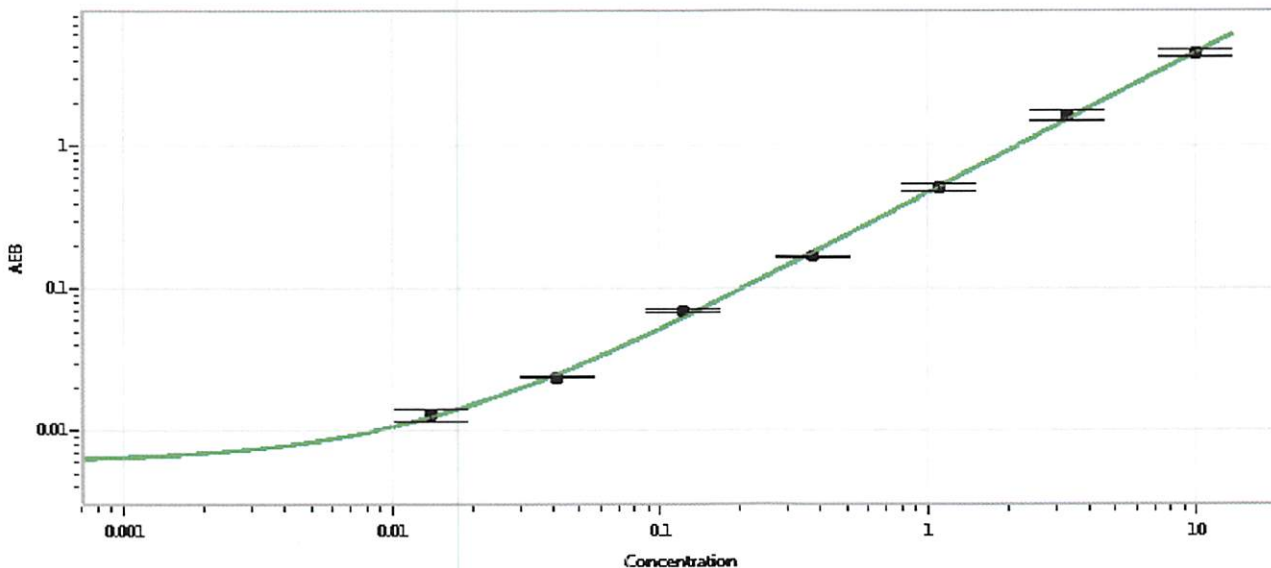
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-101816-02	0.131	0.078-0.141
Control 2	MF-101816-02	9.20	7.04-11.1
Panel 1	MF-101716-04	0.256	0.164-0.300
Panel 2	MF-101716-05	0.063	0.018-0.083
Panel 3	MF-101816-03	20.6	15.2-23.3
Panel 4	MF-050118-01	0.230	0.144-0.280

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	257
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date