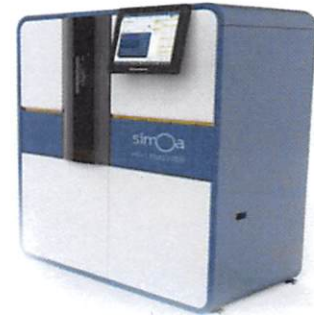


Product Number:	100200
Lot Number:	501358
Expiration:	21-Dec-2018
Platform(s):	HD-1 & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	101089	820601
Detector Reagent	101091	821308
SBG Reagent	101101	819104
Sample Diluent	101099	818005
Calibrator Diluent	101095	816308
Calibrator Concentrate	101093	63560
RGP Reagent	103159	N/A <sup>1</sup>

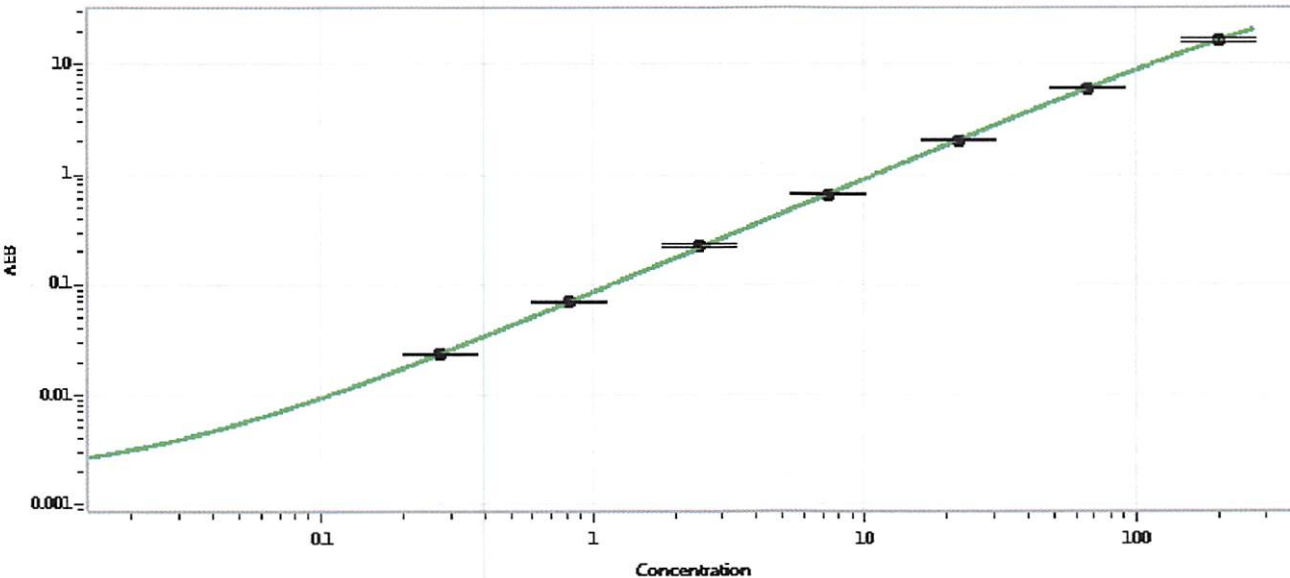
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092215-03	10.1	6.99-11.4
Control 2	MF-092215-03	111	67.3-112
Panel 1	MF-012116-06	2.43	1.17-2.99
Panel 2	MF-093015-02	4.57	1.45-5.96
Panel 3	MF-093015-03	43.9	24.1-59.5

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

Calibrator Concentrate Value Assignment (pg/mL)	1,751
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 04/11/2018  
Signature/Date