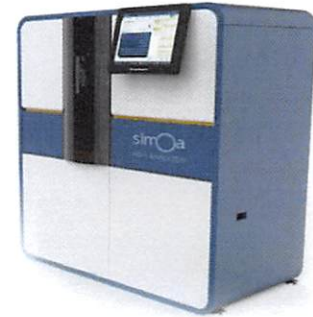


Product Number:	101580
Lot Number:	501307
Expiration:	10-Apr-2018
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101582	818407
Detector Reagent	101581	81786
SBG Reagent	101583	81782
Sample Diluent	101584	81006
Calibrator Diluent	101585	81781
Calibrator Concentrate	100655	71510
RGP Reagent	103159	N/A ¹

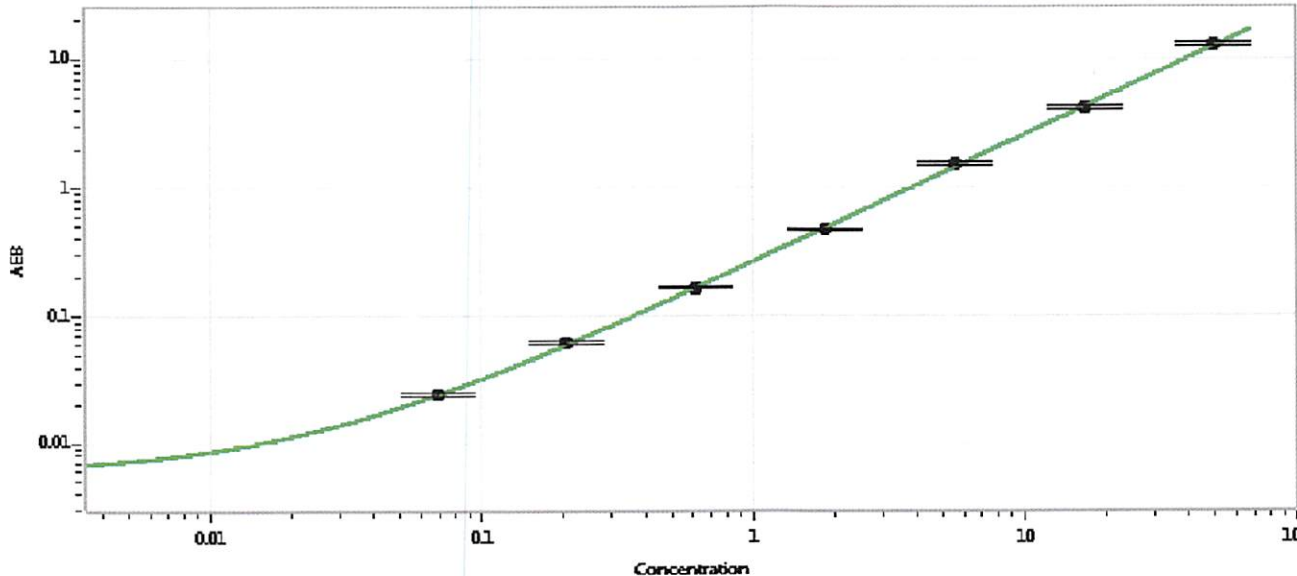
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	2.21	2.05-3.08
Control 2	MF-051617-02	44.0	35.8-53.7
Panel 1	MF-051017-03	25.3	20.9-32.4
Panel 2	MF-032217-11	1.31	1.17-2.07
Panel 3	MF-050117-05	1.98	1.19-2.11

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

Calibrator Concentrate Value Assignment (pg/mL)	1205
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 12/14/2018
Signature/Date